



NightSHADE evo LB 985N
***In vivo* Imaging System with indiGO Software**

Operating Manual
73037BA2

Rev. No.: 00, 09/2022

Not for use in in-vitro diagnostic (IVD) procedures.

The information in this guide is subject to change without notice.

DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, BERTHOLD TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

This instrument is not designed or intended for use with installations or equipment in hazardous environments. Servicing of the instrument must only be performed by Berthold Technologies Field Service Engineers or service staff authorized by Berthold Technologies.

Please contact our Service Center at service@berthold.com if you have any operational issues.

Berthold Technologies GmbH & Co. KG

Calmbacher Str. 22
75323 Bad Wildbad, Germany
www.berthold.com

Telephone +49 7081 177-0
Fax +49 7081 177-100
bio@berthold.com

Contents

1	PREFATORY COMMENTS	7
1.1	History of Changes	7
1.2	Typographical conventions	7
1.3	Warning signs at the instrument	7
2	SAFETY INSTRUCTIONS	9
2.1	Consignes de Sécurité	12
2.2	Sicherheitshinweise	15
3	WARRANTY AND CUSTOMER SERVICE	18
3.1	Warranty Statement	18
3.2	Customer Service	19
4	QUICK START	20
4.1	Installation	21
4.2	First Acquisitions	40
4.3	Image Processing and Evaluation Functions	59
4.4	Creating a Report	69
4.5	Using the Scheduler	73
5	SYSTEM DESCRIPTION	77
5.1	Intended Use	78
5.2	Design of the Acquisition Device	79
5.3	Acquisition Chamber	81
6	INDIGO SOFTWARE DOCUMENTATION	87
6.1	Graphical User Interface	88
6.2	Application Toolbar	88
6.3	Identity Context	89
6.4	Identities and Measurements	91
6.5	Contrast Slider	91
6.6	Data Slider	94
6.7	Toolbar	97
6.8	Status Bar	101
6.9	IndiGO Menu	102
6.10	Tools Menu	104
6.11	Service Menu	106
6.12	View Menu	112
6.13	Batch Image Export	112
7	TROUBLE SHOOTING	113
7.1	Hardware	113
7.2	Software	113
8	CLEANING AND MAINTENANCE	114
8.1	Camera and Lens	115
8.2	Darkbox	116
9	TECHNICAL DATA	119
10	APPENDIX	120
10.1	Accessories	120

10.2 Index

128

1 Prefatory Comments

1.1 History of Changes

Revision	Date	Chapter / Page	Change	Author
Rev. 00	09/2022	---	Initial release	R. Kembügler

1.2 Typographical conventions

The following typographical conventions will be used in this operating manual:



Caution! Danger!

☒ **Important information**

Menus and **options** in the **software** are printed in bold type **Measurement, View**

The **selection of a menu item** is written as a sequence of commands, separated by a vertical line

The names of **dialog boxes** and **parameters** are also printed in bold type

Buttons are printed in bold type inside angular brackets

Enumerations are emphasized by

Action **steps** are identified by

Example

Device | Configuration

Filter Setting

<Start>, <Stop>

☒

☒

1.3 Warning signs at the instrument

	<p>Wear eye protection When using the LED panel an eye protection (DIN EN 166) is required to prevent from high intensity light caused by the LED lights.</p>
	<p>Caution Danger This sign indicates a general danger (e.g. switched power socket inside the instrument).</p>
	<p>Dangerous electrical voltage The switched power socket inside the instrument carries high voltage. Be careful when handling this device.</p>
	<p>Hot surface The excitation light source can become hot. Be careful when handling these device.</p>

2 Safety Instructions



Caution! This sign alerts you to important operating procedures with a potential danger of damaging the equipment and endangering your safety on disobeying. Refer to the user and instrument manuals for precautionary instructions.



This operating manual includes information and warnings that have to be observed by the user in order to ensure safe operation of the instruments.



Please do always act according to the following safety instructions, before as well as during operation of the system! Before set up and operation of the instrument it is necessary to read the instructions below as neither safe operation of the instrument nor safety of the user are guaranteed otherwise. Failure to follow the instructions may invalidate the warranty.

The instrument has been manufactured in accordance with the safety requirements for electronic and medical measuring systems. If the law lays down regulations on the installation and/or operation of sample measuring system, then it is the operator's responsibility to adhere to them.

The instrument is designed in a way that no harmful noise levels are emitted by the instrument itself. If the user detects unusual noise from the instrument the service must be contacted.

The manufacturer has done everything possible to guarantee that the equipment functions safely, both electrically and mechanically. The user has to make sure that the instrument will be set up and installed properly to guarantee safe operation.

The instruments are tested by the manufacturer and supplied in a condition that allows safe and reliable operation. Improper use of the instrument can disable safety features and could lead to an unsafe operation of the instrument



- This equipment must be installed and used in accordance with the manufacturer's recommendations. Installation must be performed by properly trained and authorized personnel.
- The over-voltage category is II.
- The instrument may only be operated by personnel who have been trained on the use of the system. It is strongly recommended that all users read this manual prior to use.



- Never reach into the instrument while the unit is in operation, neither with parts of your body nor with other instruments.**
- Remove the transportation lock before switching on the instrument.**
- Use the instrument only for the designated application.
- The instrument is designed for indoor use only (dry environment).
- The instrument may not be operated at altitudes above 2000 m above sea level.
- The instrument is designed to be operated within a temperature range of 10 to 28 °C. Operating the instrument at higher temperatures will lead to problems with the camera cooling (safety shutdown of the camera cooler)
- The instrument is designed to be operated at a maximum relative humidity of 80 % for temperatures up to 28 °C (non condensing).

- ❑ BERTHOLD TECHNOLOGIES assumes no liability for any damages, including those to third parties, caused by improper use or handling of the instrument.
- ❑ The user is responsible for connecting the instrument in accordance with the valid regulations for electrical instruments.
- ❑ Set the instrument up to ensure easy access to the mains switch.
- ❑ The mains supply voltage fluctuations must not exceed +/- 10 %. Operating voltage is AC 100- 240V +/- 10 %.
- ❑ The instrument is provided with a 3-pole grounded plug. If your wall outlet does not allow connection of a 3-pole plug, have a suitable wall outlet installed by qualified personnel or use an adapter for safe grounding. Please observe the safety specifications of the grounded plug. Only use the provided power cable or a cable that fits the specifications of the power cable provided with the instrument.



- ❑ The instruments are designed according to the IEC 61010-1 or EN 610 10-1 regulations for electrical measuring systems.
- ❑ To disconnect the instrument from power the appliance coupler has to be removed from the mains. Ensure that the distance to the wall is high enough to allow an easy removal of the mains cable.
- ❑ Do not open any instrument doors as long as the instrument is in operation.
- ❑ Service and repair work must be carried out by certified personnel only. After any service work the secure operation of the instrument must be tested.
- ❑ The operator may only perform the maintenance work described in this user guide.
- ❑ There are no exchangeable electrical components in the instrument. In case of malfunction call authorized service personnel.
- ❑ Use only parts described in this manual for servicing.
- ❑ Disconnect power supply before opening the instrument.
- ❑ Pull the power cord to disconnect instrument from power supply.
- ❑ Turn instrument off before pulling the power cord.
- ❑ Upon removal of the front and top parts of the housing no safety measures are in effect. Be aware of any moving parts. The interior of the instrument may reach temperatures that can cause burns. Some parts of the instrument may remain hot without visual indication for some time after the power has been turned off.



- ❑ The internal battery can explode if not properly inserted. The internal battery must be replaced with the original spare part (ID 17391) by certified service personnel. Used batteries must be recycled based on the manufacturer's directions.
- ❑ If you can see that the instrument has become unsafe to use, switch it off and disconnect it from power supply.



- ❑ If liquid gets inside the instruments, pull the power cord. Clean the unit or have it cleaned by an authorized service center.
- ❑ Protect yourself from electrostatic charge, as discharge could damage sensitive instrument parts, especially sensitive parts of the computer and electronics boards.

- ❑ When the lid is opened (e.g. filter change) ESD can no longer be guaranteed. To avoid any damages to the electronic parts it is recommended to take precautions (touching the metal case of a safety grounded object, wearing a grounding strap, etc.).
- ❑ The system always has to be primed with solutions recommended by the kit manufacturer.
- ❑ Use only reagents recommended by the kit manufacturer.
- ❑ Use reagents only in accordance with the kit manufacturer's instructions.
- ❑ Do not use any flammable or explosive solutions or liquids whose mixture is flammable or explosive.
- ❑ For instrument cleaning, please refer to the respective sections in this manual.
- ❑ Reliable instrument function can be guaranteed only when original spare parts are used.
- ❑ Pollution degree is 2.
- ❑ The instrument may only be used in rooms with a max. pollution degree of 2.
- ❑ This instrument must not be used in areas with potentially explosive atmosphere.

The tests and service work recommended by the manufacturer has to be performed to make sure that the operator remains safe and that the instrument continues to work correctly. Any service and maintenance work not described in this user guide has to be performed by authorized service personnel. Use the instruments only for the designated application.

2.1 Consignes de Sécurité

Attention! Ce symbole d'alarme, vous avertit de prêter attention aux consignes opératoires. En effet si vous ne suivez pas ces instructions, il peut y avoir un risque d'endommagement du matériel et également vous faire courir des risques pour votre propre sécurité. Il est impératif de respecter les instructions du mode d'emploi et de les respecter.



Ce mode d'emploi contient des informations et avertissements qui doivent être suivis par l'utilisateur afin de garantir un fonctionnement sûr des instruments.



Il est impératif de respecter les consignes de sécurité suivantes non seulement avant la mise en service mais aussi pendant le fonctionnement de l'appareil! Avant l'installation et mise en service de l'instrument tous les utilisateurs des appareils sont tenus de lire d'abord ces instructions de service, autrement ni le fonctionnement correct de l'appareil ni la sécurité de l'utilisateur peuvent être garantis. Ne pas suivre ces instructions de service peut invalider la garantie.

Le appareil a été fabriqué conformément aux prescriptions de sécurité en vigueur pour les appareils de mesure électroniques et médicaux. Si l'installation et/ou l'utilisation des appareils de mesure de prélèvements-échantillons sont/est soumise(s) à des réglementations prescrites par la loi, il appartient à l'utilisateur de les respecter.

Le constructeur a fait tout le nécessaire pour assurer le fonctionnement sûr des appareils (du point de vue électrique, électronique et mécanique). L'utilisateur est tenu de veiller à ce que les appareils soient installés correctement afin d'éviter toute altération de leur utilisation sûre.

Les appareils sont contrôlés à l'usine et livrés dans un état assurant la sécurité de fonctionnement.



- Les appareils ne peuvent être mis en service que par des personnes autorisées et ne peuvent être utilisés que par un personnel formé. Tous les utilisateurs qui travaillent avec les appareils doivent d'abord lire ce mode d'emploi.
- La catégorie de surtension est II.
- Les appareils ne doivent être utilisés que par du personnel formé à cet effet. Il est recommandé à tous les utilisateurs de lire ce mode d'emploi avant d'utiliser l'appareil.
- Retirer les sécurités de transport avant la mise en marche.
- Les appareils ne doivent être utilisés que pour l'usage prévu.
- Berthold Technologies n'assume aucune garantie, y compris pour les dommages causés à des tiers, en cas de manipulation non conforme des appareils.
- Les appareils ne doivent être utilisés qu'à l'intérieur de locaux fermés (environnement sec).
- Les appareils ne doivent pas être utilisés à des altitudes supérieures à 2000 m au-dessus du niveau de la mer.

- ❑ Les appareils sont conçus pour fonctionner dans une plage de température de 10 à 28 °C. Des températures ambiantes plus élevées affectent le refroidissement de la caméra (arrêt de sécurité du refroidissement de la caméra).
- ❑ Les appareils sont conçus pour fonctionner avec une humidité relative maximale de 80 % (jusqu'à 28 °C) (sans condensation).
- ❑ L'alimentation électrique ne doit pas dépasser ± 10 % de la valeur nominale. La plage de tension est de 100 - 240 V AC ± 10 %.
- ❑ Il est de la responsabilité de l'utilisateur de s'assurer que les appareils sont installés conformément aux réglementations électriques locales.



- ❑ Les appareils sont équipés d'un câble d'alimentation à 3 pôles. Il s'agit d'un équipement de sécurité. Si la prise de courant ne prend pas en charge une connexion à 3 pôles, un électricien spécialisé doit installer une prise de courant à 3 pôles appropriée ou fournir un adaptateur approprié pour la mise à la terre de la connexion. Ne détruisez jamais les dispositifs de sécurité du raccordement à la terre. Utilisez uniquement le câble d'alimentation fourni ou un câble de même spécification et de même certification.

- ❑ Les appareils sont conformes aux prescriptions des normes CEI 61010-1 et EN 61010-1 relatives aux appareils de mesure électriques.



- ❑ Ne pas ouvrir l'appareil lorsqu'il est en service.
- ❑ Les travaux de service et de réparation ne doivent être effectués que par des spécialistes. Après les travaux, le fonctionnement sûr doit être contrôlé.
- ❑ Seuls les travaux de maintenance décrits dans le manuel peuvent être effectués par l'utilisateur.
- ❑ Lors des travaux d'entretien, seules les pièces indiquées doivent être utilisées.
- ❑ Avant d'ouvrir l'appareil, l'alimentation électrique doit être coupée.
- ❑ Pour déconnecter complètement l'appareil du secteur, il est possible de débrancher le câble d'alimentation. Il faut veiller à ce que la distance par rapport au mur soit suffisante pour que le câble d'alimentation puisse être retiré sans problème.



- ❑ Éteindre l'appareil avant de le débrancher.
- ❑ Lorsque l'appareil est ouvert, les mesures de sécurité ne fonctionnent plus. Attention aux composants mobiles ! L'intérieur des appareils peut atteindre des températures qui peuvent provoquer des brûlures. Certaines pièces peuvent rester chaudes sans signe visible, même après l'arrêt de l'appareil.
- ❑ Tous les appareils et accessoires livrés doivent être raccordés au réseau avec mise à la terre. Utiliser une fiche de sécurité !



- ❑ Attention : la batterie interne peut exploser si elle est mal installée ! La batterie interne ne peut être remplacée que par le même modèle ou par un modèle de remplacement recommandé (ID 17391) par une personne autorisée par Berthold Technologies. Les piles usagées doivent être éliminées conformément aux instructions du fabricant.

- ❑ Installez l'appareil de manière à pouvoir le mettre en marche et l'arrêter facilement.
- ❑ Si la sécurité de fonctionnement est compromise, les appareils doivent être éteints et débranchés.
- ❑ Si du liquide a pénétré à l'intérieur de l'appareil, débrancher la fiche secteur. Ouvrir l'appareil et le nettoyer ou le faire nettoyer par un centre de service agréé.
- ❑ Lors de l'ouverture de l'appareil, il faut éviter les charges électrostatiques (p. ex. dues à la moquette), car les décharges sur l'appareil peuvent endommager les pièces électroniques.
- ❑ Ne pas utiliser de liquides inflammables ou explosifs, ni de liquides dont le mélange est inflammable ou explosif.
- ❑ Respectez toutes les dispositions légales relatives à la manipulation des déchets biologiques, des réactifs et des échantillons de patients.
- ❑ Pour le nettoyage de l'appareil, veuillez consulter la partie correspondante de ces instructions d'utilisation.
- ❑ Le bon fonctionnement ne peut être garanti que si les pièces de rechange d'origine sont utilisées.
- ❑ Le degré de pollution est de 2.
- ❑ L'appareil ne doit pas être utilisé dans des zones à risque d'explosion.
- ❑ L'appareil ne peut être utilisé que dans des locaux présentant un degré de pollution maximal de 2.

Afin d'assurer la sécurité de l'utilisateur et le bon fonctionnement des appareils, effectuer les travaux d'inspection et d'entretien recommandés par le fabricant. Toutes les mesures d'entretien et de réparation allant au-delà de celles spécifiées dans ce manuel sont réservées aux techniciens autorisés.

2.2 Sicherheitshinweise



Die vorliegende Bedienungsanweisung enthält Informationen und Warnungen, die vom Benutzer befolgt werden müssen, um einen sicheren Betrieb der Geräte zu ermöglichen.



Dieses Zeichen weist den Benutzer auf wichtige Punkte hin, deren Beachtung unerlässlich ist.



Die folgenden Sicherheitshinweise sind sowohl vor der Inbetriebnahme als auch während des Betriebs des Gerätes unbedingt zu beachten. Vor Inbetriebnahme des Gerätes ist es zwingend erforderlich, die Bedienungsanleitung zu lesen, da ansonsten die Sicherheit des Gerätes und des Benutzers nicht gewährleistet wird.

Das Gerät wurde in Übereinstimmung mit den Sicherheitsanforderungen für elektronische und medizinische Messgeräte hergestellt. Bestehen für die Errichtung und/oder den Betrieb von Probenmessgeräten gesetzlich vorgeschriebene Regelungen, so ist es die Aufgabe des Errichters und Betreibers, diese einzuhalten.

Das Gerät emittiert keine gefährlichen Schallpegel. Sollten dennoch Geräusche wahrgenommen werden, die auf einen fehlerhaften Betrieb hinweisen, ist der Service zu rufen.

Der Hersteller hat alles unternommen, um ein sicheres Arbeiten der Geräte (bezüglich Elektrik, Elektronik und Mechanik) zu gewährleisten. Der Benutzer muss dafür sorgen, dass die Geräte so aufgestellt und installiert werden, dass ihr sicherer Gebrauch nicht beeinträchtigt wird.

Die Geräte sind werkgeprüft und wurden in betriebs sicherem Zustand ausgeliefert.



- ❑ Die Geräte dürfen nur von autorisierten Personen in Betrieb genommen und nur von eingewiesenem Personal bedient werden. Alle Benutzer, die mit den Geräten arbeiten, müssen zuerst diese Bedienungsanleitung lesen.
- ❑ Überspannungskategorie ist II.
- ❑ Die Geräte dürfen nur von dafür geschultem Personal betrieben werden. Es wird allen Anwendern empfohlen, diese Bedienungsanleitung vor Benutzung zu lesen.
- ❑ Transportsicherungen vor dem Einschalten entfernen.
- ❑ Die Geräte dürfen nur für den vorgesehenen Zweck eingesetzt werden.
- ❑ Berthold Technologies übernimmt keinerlei Gewährleistung, auch für Schäden gegenüber Dritten, die durch unsachgemäße Handhabung der Geräte hervorgerufen werden.
- ❑ Die Geräte dürfen nur innerhalb geschlossener Räume betrieben werden (trockene Umgebung).
- ❑ Die Geräte dürfen nicht in Höhen von mehr als 2000 m über dem Meeresspiegel betrieben werden.

- ❑ Die Geräte sind dafür ausgelegt, innerhalb des Temperaturbereiches von 10 bis 28 °C betrieben zu werden. Höhere Umgebungstemperaturen beeinträchtigen die Kühlung der Kamera (Sicherheitsabschaltung der Kamerakühlung).
- ❑ Die Geräte sind dafür ausgelegt, bei einer maximalen relativen Luftfeuchte von 80 % (bis zu 28 °C) betrieben zu werden (nicht kondensierend)
- ❑ Die Stromversorgung darf nicht mehr als ± 10 % des Nominalwertes aufweisen. Der Spannungsbereich beträgt AC 100 – 240 V ± 10 %
- ❑ Es liegt im Verantwortungsbereich des Anwenders, dass die Geräte nach den lokalen elektrischen Vorschriften installiert werden.
- ❑ Die Geräte sind mit einem 3-poligen Netzkabel ausgestattet. Dies ist eine Sicherheitsausstattung. Wenn die Steckdose keinen 3-poligen Anschluss unterstützt, muss ein Fachelektriker eine passende 3-polige Steckdose installieren oder einen passenden Adapter zur Erdung des Anschlusses bereitstellen. Zerstören Sie niemals die Sicherheitsvorkehrungen des geerdeten Anschlusses. Verwenden sie nur das beigelegte Netzkabel oder eines mit der gleichen Spezifikation und Zertifizierung.
- ❑ Die Geräte entsprechen den Vorschriften der IEC 61010-1 und EN 61010-1 für elektrische Messgeräte.
- ❑ Nicht öffnen, wenn das Gerät in Betrieb ist.
- ❑ Service- und Reparaturarbeiten dürfen nur von Fachleuten ausgeführt werden. Nach den Arbeiten muss der sichere Betrieb geprüft werden.
- ❑ Es dürfen nur die im Handbuch beschriebenen Wartungsarbeiten vom Anwender ausgeführt werden.
- ❑ Bei Wartungsarbeiten dürfen nur die angegebenen Teile verwendet werden.
- ❑ Vor dem Öffnen des Gerätes ist die Stromzufuhr zu unterbrechen.
- ❑ Um das Gerät vollkommen vom Netz zu trennen, kann das Netzkabel gezogen werden. Es ist auf ausreichend Abstand zur Wand zu achten, damit das Netzkabel problemlos abgezogen werden kann.
- ❑ Gerät ausschalten, bevor der Stecker gezogen wird.
- ❑ Wenn das Gerät geöffnet ist sind Sicherheitsmaßnahmen nicht mehr in Betrieb. Auf bewegliche Komponenten achten! Das Innere der Geräte kann Temperaturen erreichen, die Verbrennungen verursachen können. Einige Teile können heiß bleiben ohne sichtbare Zeichen, auch nachdem das Gerät abgeschaltet worden ist.
- ❑ Alle gelieferten Geräte und Zusatzgeräte sind geerdet ans Netz anzuschließen. **Schutzkontaktstecker** verwenden!
- ❑ **Achtung:** Die interne Batterie kann bei falschem Einbau explodieren! Die interne Batterie darf nur mit dem gleichen oder einem empfohlenen Ersatzmodell (ID 17391) von einer von Berthold Technologies autorisierten Person ausgetauscht werden. Verbrauchte Batterien müssen entsprechend den Herstelleranweisungen entsorgt werden.



- ❑ Stellen Sie das Gerät so auf, dass Sie es leicht ein- und ausschalten können.
- ❑ Bei Beeinträchtigung der Betriebssicherheit sind die Geräte abzuschalten und vom Netz zu trennen.
- ❑ Ist Flüssigkeit in das Innere des Gerätes gelangt, Netzstecker ziehen. Das Gerät öffnen und reinigen bzw. durch eine autorisierte Servicestelle reinigen lassen.
- ❑ Elektrostatische Aufladungen (z.B. durch Teppichböden) müssen beim Öffnen des Gerätes verhindert werden, da Entladungen am Gerät zur Beschädigung elektronischer Teile führen können.
-  ❑ **Es dürfen keine entzündlichen oder explosiven Flüssigkeiten oder keine Flüssigkeiten, deren Mischung entzündlich oder explosiv ist, verwendet werden.**
- ❑ Beachten Sie alle gesetzlichen Vorschriften für den Umgang mit biologischem Abfall, mit Reagenzien und Patientenproben.
- ❑ Zum Reinigen des Gerätes bitte den entsprechenden Teil dieser Bedienungsanleitungen beachten.
- ❑ Ordnungsgemäße Funktionalität kann nur bei Verwendung der Originalersatzteile garantiert werden.
- ❑ Verschmutzungsgrad ist 2.
- ❑ Das Gerät darf nicht in explosionsgefährdeten Bereichen verwendet werden.
- ❑ Das Gerät darf nur in Räumlichkeiten mit einem maximalen Verschmutzungsgrad von 2 betrieben werden.

Für die Sicherheit des Benutzers und die Funktionsfähigkeit der Geräte sind die vom Hersteller empfohlenen Überprüfungen und Wartungsmaßnahmen durchzuführen. Alle über die Betriebsanleitung hinausgehenden Wartungs- und Instandhaltungsmaßnahmen dürfen nur von autorisierten Technikern ausgeführt werden.

3 Warranty and Customer Service

3.1 Warranty Statement

The instrument is sold in accordance with the general conditions of sale of Berthold Technologies GmbH & Co KG and its affiliates and representatives.

Berthold Technologies warrants this product to be free of defects in material and workmanship for a period of 12 months from the date of delivery, ex works Bad Wildbad.

Berthold Technologies or its authorized representative will repair or replace, at its option and free of charge, any product that under proper and normal use proves to be defective during the warranty period.

Berthold Technologies shall in no event be liable or responsible for any incidental or consequential damage, either direct or indirect.

The above warranty shall not apply if:

- a) the product has not been operated in accordance with the operating manual
- b) the product has not been regularly and correctly maintained
- c) the product has not been repaired or modified by a Berthold Technologies authorized representative or user
- d) parts other than original Berthold Technologies parts are used
- e) the product and parts thereof have been altered without written authorization from Berthold Technologies GmbH & Co KG
- e) the product has not been returned properly packed in the original Berthold Technologies packaging

This warranty does not apply to any third-party product involved in the application.

Berthold Technologies reserves the right to refuse to accept the return of any product that has been used with radioactive or (micro)-biological substances, or any other material that may be deemed hazardous to employees of Berthold Technologies. Such products have to be properly decontaminated and marked.

Before returning products to Berthold Technologies ensure the devices are properly decontaminated.

A return material authorization (RMA) number must be obtained and clearly identified on the packing and documents. Please visit the Berthold Technologies website and fill in the required information to obtain the number.

<https://www.berthold.com/en/company/service/form-rma-return-material-authorization/>

Retain the original packaging for use if the instrument needs to be returned to Berthold Technologies.

3.2 Customer Service

Customer service will be provided in the first instance by the network of Berthold Technologies representatives. In the event of any problem experienced with your instrument, the first recourse should be **your local Berthold Technologies representative**. For further problems requiring hardware or software expertise, the Technical Support group at Berthold Technologies GmbH & Co KG will be available by phone, fax or email to deal with your queries.

Please use the following contact for your enquiries.

Berthold Technologies GmbH & Co.KG
Technical Support
Calmbacher Str. 22
75323 Bad Wildbad
Germany
Phone: +49 7081 177 171
Fax: +49 7081 177 301
Email: service@berthold.com
<https://www.berthold.com/en/company/service/form-more-services/>

Please visit the Berthold Technologies website for any service inquiries.

Please also make sure that you have the relevant information available before contacting Berthold Technologies. Helpful information would include:

- serial numbers, part number, revision:
see production label on instrument
- software and firmware versions
- monitor and log files (refer to the respective service manuals)

4 Quick Start

These quick start instructions provide a quick access to the program. You are guided step by step through the major program functions. The page numbers in the headers refer to the individual sections in the manual that provide more detailed information about the respective functions.

4.1	Installation	21
4.1.1	Requirements	21
4.1.2	Unpacking and Carrying	22
4.1.3	Camera Installation	23
4.1.3.1	Top view camera	23
4.1.3.2	Side view Camera	24
4.1.4	Connecting	25
4.1.5	Power On	25
4.1.6	Connecting to the PC	25
4.1.7	Software Installation	26
4.1.8	Instrument Controller First Time Initialization Setup	28
4.1.9	First Time Initialization Setup	31
4.1.10	User Management	34
4.1.11	Installing Fluorescence Filters	35
4.1.12	Filter Management	37
4.1.13	Filter Inventory	38
4.1.14	Filter Manager	39
4.1.15	Filter Manager – Backup and Restore	39
4.2	First Acquisitions	40
4.2.1	Quick Start Glossary	40
4.2.2	Launching the software	41
4.2.3	Defining a Measuring Template	41
4.2.4	Defining a Controller Template	47
4.2.5	Defining a Project	51
4.2.6	Starting an Acquisition	54
4.2.7	Image Sequences	56
4.3	Image Processing and Evaluation Functions	59
4.3.1	Load Images	59
4.3.2	Image Window	60
4.3.3	Line Plot	63
4.3.4	Surface Plot	64
4.3.5	Peaksearch	65
4.3.6	Define Areas Manually	66
4.4	Creating a Report	69
4.4.1	Creating a Rich-Text-Report	69
4.4.2	Excel Export	70
4.5	Using the Scheduler	73

4.1 Installation

4.1.1 Requirements

Please observe the following requirements when working with the NightSHADE evo:

- 2-5 mains outlets (depending on accessories). The maximum power consumption of the NightSHADE evo without PC and external devices is max. 600 W. The power consumption of the external cooling unit is 340 W. To power on/off the system with one switch we recommend connecting all devices to a common outlet strip.
- At least 160 cm width of workplace (the load-carrying capacity of the table or the desktop should be approx 100 kg). Make sure that the table or desktop stands firm and does not wobble.
- To use network resources (network printer, file server, etc.) a connection to a network is essential. A connection to the Internet is recommended for easier support, but it is not essential for operation of the NightSHADE evo.
- To allow ergonomic working, the NightSHADE evo should always be set up to the left of the PC, as the door hinges are on the left.
- To fill the cooling unit for the LED panels up to 4 liters of distilled water are required. It is recommended to place the external cooler on the floor (under the lab table etc.). Please note that a tubing connection to the NightSHADE evo is required for the coolant. In case the cooler is placed under the lab table a cut-out on the table area is required to get shortest possible tubing.



4.1.2 Unpacking and Carrying

The NightSHADE is delivered in a so-called flight case. The case consists of 3 parts, the lower base and two upper shells. The upper shells are locked to the base with butterfly locks and can be removed after releasing the locks.

Since the lower base of the flightcase is equipped with wheels, the instrument can be placed in front of the table to allow an easy placement on the table top.



The NightSHADE evo is heavy. To unpack the instrument from the flight case and lift it to the table at least 2 persons are required.

When the upper shells are removed the black handles at the side of the instrument can be popped out to carry the instrument.

One person on each side of the instrument must grab the handle to lift the instrument carefully.



The wheels must be secured using the built-in brake when the box is not used or stored to avoid uncontrolled movement of the flight case.

To activate the break just push the pedal (see image) at the corresponding wheel. Pull back the pedal to release the brake.

To lift up the instrument to the table both persons holding the handle must lift up the instrument simultaneously and lift it out of the lower base and up on the table. A third person that secures the instrument with both hands at the base plate after lifting it out of the lower base is recommended. Once the instrument is placed on the table the handles can be folded in.

- ❑ Unpack the boxes and make sure that the delivery is complete. If you should discover any sign of damage on the instruments or instrument parts, please notify the shipping agent and/or the distributor immediately.
- ❑ Provide adequate space for setting up the instruments. To set up the NightSHADE evo, the computer and the printer next to each other about 2 meters of desk space are required.
- ❑ The device feet of the NightSHADE evo are height-adjustable to compensate a possible unevenness of the table. By turning the device feet, their height can be adjusted to ensure a safe foothold of the NightSHADE evo.

4.1.3 Camera Installation

For safety reasons, the camera is delivered in a special case to prevent damage during transport. Therefore, the camera has to be installed first before it is taken into operation.
Please note: Do not turn on the NightSHADE evo as long a no camera has been installed.

The camera is connected to the instrument using two cables. One cable provides the supply voltages (2), the other USB-cable (3) is to control the camera and readout the data.



The cable (1) is used to supply the fan in the top cover to allow a proper air exchange. Make sure this cable is plugged in correctly as otherwise the fan will not work and the camera could get cooling problems during longer operations.

After switch on, the device is initialized.

If the device is transported over larger distances (e.g. in a car or truck), you have to dismantle the camera prior to the transport to rule out that the camera will get damaged. To prevent damage please remove or secure the base plate whenever you transport the device.

4.1.3.1 Top view camera

The NightSHADE evo camera has a USB port to control the camera and to transfer data. The camera is delivered in a separate case which also accommodates other accessories, such as the connection cable etc. Make sure no protective stickers are in front of the camera (protecting the shutter hole).



Install the camera that the red dot located on the bajonet lock is matching the red dot on the NightSHADE evo socket and fix it in the bayonet socket until it clicks into place. Connect the power supply cable (1) to the camera and as well as the outlet socket and screw it.

Make sure that the USB-connector (2) is plugged into the camera and outlet socket properly.

With this the installation of the camera is completed.

4.1.3.2 Side view Camera

The NightSHADE evo offers the special option that the camera can be mounted on the side, too. This feature enables the system to take images of plants from the side or to measure the sample dishes in an upright standing position. The upright standing position is used when the sample changer is assembled into the NightSHADE evo.

The power supply and the USB connection for the camera is located at the rear side of the instrument. Please use the supplied special cables to connect the camera to the instrument.

To assemble the camera at the side proceed as follows:

- Turn off the NightSHADE evo.
- Open the top cover and remove the USB cable and the power supply cable from the camera.
- Remove the blind plug from the side view camera holder. To remove the cover, turn it slightly.
- Dismantle the camera by turning it counterclockwise until it can be removed
- Place the camera in the same way at the camera holder at the left side of the instrument. Make sure that the red dot at the camera is matching the red dot at the instrument ba-jonett.
- Plug the power supply cable to the rear panel and then to the camera.
- Plug the USB control cable to the rear panel and then to the camera.
- Put the blind plug to the top position of the camera to avoid incoming light.
- Switch on the NightSHADE evo.



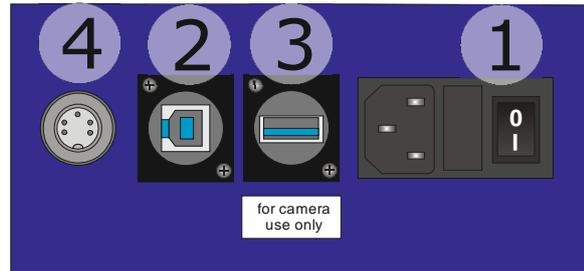
4.1.4 Connecting

The power supply and device communication ports are located on the rear panel of the NightSHADE evo. Plug the power cord into the mains inlet unit(1). The main power switch is located next to the mains inlet.

Connect the NightSHADE evo to the PC via the USB3 B-socket (2) on the rear panel using a USB3 cable (A to B).

The 7-pin Tuchel socket (4) is the power supply for the side mounted camera, the USB3 socket (3) is used for the control of the side view camera only.

To ensure secure operation only connect certified IT equipment to the instrument.



Check to make sure that the system voltage indicated on the instruments matches the line voltage.

If possible connect all devices via *one* outlet strip to the mains supply.

The NightSHADE evo is initialized after power on; a control lamp on the front panel is flashing.

The NightSHADE evo camera is ready for operation after it has cooled down to operating temperature; this takes approximately 10 minutes.

The cooling of the camera is initialized when the indiGO software is started (controller startup). Switching on the NightSHADE evo without starting the software will not cool down the camera.

The NightSHADE should only be connected to mains power supply systems that are secured by a 10A automatic circuit breaker.

4.1.5 Power On

After power on, the device starts an automatic initialization where all relevant components will be initialized. During this time, the LED on the front panel is flashing red. After the successful initialization the NightSHADE evo is ready for operation.

The indioGO software automatically recognizes the NightSHADE evo. No further steps are necessary.

4.1.6 Connecting to the PC

To assure the proper operation of the NightSHADE evo device drivers are necessary. All device drivers are automatically installed through the indiGO software setup.

When connecting the instrument to PC, new devices are detected through the windows device manager. Using the automatic installation (not using the Windows update driver repository) the drivers are installed automatically. This procedure can take a while.

Do not connect the instrument to the PC when the indiGO software has not been installed yet.

4.1.7 Software Installation

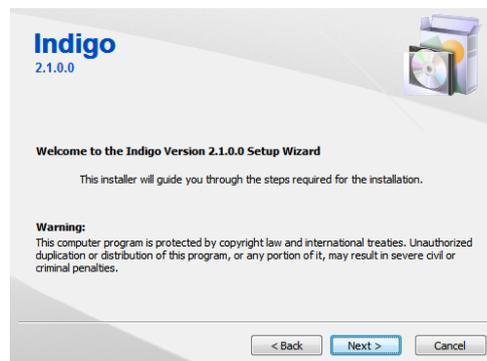
The indiGO software is included on the enclosed USB stick.

Use this media to install the software in order to use the correct individual calibration data which is automatically copied during installation. Keep this media in a secure place. You will need it again e.g. for a new installation on a different PC.

For the operation of the NightSHADE evo the indiGO software version 2.1 or higher is necessary.

We recommend using Windows 10 as operating system. Older operating systems than Windows 7 are not supported.

To start the installation program, click on the file **SETUP.EXE** on the data carrier. After launching the installation program a welcome screen appears:



After the welcome screen the folder for the software installation must be defined. If possible confirm the suggested folder. Click on **Next >** to get to the next page.



IndiGo can be installed in two different configurations. The evaluation or demo version will install a preconfigured system with instrument simulation to work without an instrument connected or to use it on office PCs for offline data evaluation. In this case select "Evaluation – Demo Version". The installation will start immediately.

To update an existing indiGO installation (Evaluation or full version) or new installation with full instrument control, select "Full/Update Version with Instrument control".

Depending on the hardware, a default selection of system filters might be available and can be installed through the setup. If you want to install a set of default filter slider and wheel tick the box "Install default filters". The default filters are empty filter sliders without any filters.



In this case the default systems filter management file will be copied to your computer. If you reinstall the software the current active systems filter management file will be overwritten.

To get to the next screen asking for the drivers to be installed on the system click on 



When working with the NightSHADE evo the PCI drivers are not required. Use the selections as suggested above.

Note: Due to the fact that the setup will install device drivers for the NightSHADE evo hardware on the computer, the Windows-system might show a warning that a signature of a Windows driver cannot be verified. In this case allow the setup to install these Windows drivers.

After the installation the controller will be initialized for the first time.

If the installation was performed using an instrument specific installation media (USB Stick) the configuration procedure can be slightly different since the installer has already copied the instrument configuration to disk from the USB Stick during installation.

4.1.8 Instrument Controller First Time Initialization Setup

The **Instrument Controller First Time Initialization Setup** starts via the setup. Please note that this Setup will only be executed when the software is installed on a computer for the first time without calibration data.

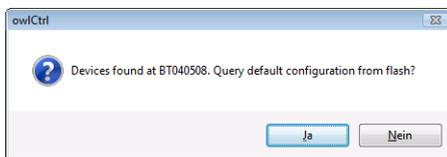


Select **Next >** to continue.



Please enter the instrument ident-number and the serial number. You can find both numbers on the back of the instrument.

You can check the instrument ident-number by starting „**Scan Devices**“. If the calibration for your device was stored on the NightSHADE evo by Berthold Technologies before it was delivered you can download the stored calibration parameters from the instruments flash memory now by confirming „**Yes**“.



If you want to recalibrate your instrument please select „**No**“.

You can load a default calibration for your instrument type via „**Load factory default settings**“. The factory default settings are chosen by the system via the given instrument ident-number.

Select **Next >** to continue.



This step of the NightSHADE evo Controller First Time Initialization Setup allows the import of further created configuration backups and calibration files.

Select **Next >** to continue.



To maintain the instrument the instruments controller needs to create an administrator account. Please enter a name and a password for the administrator account.

Select **Next >** to continue.



In the next screen log on in order to verify the administrator account.

Select **Next >** to continue.



Next, choose a start method for the controller. . With regular installations it is recommended to select the first option ("Start controller with indiGO"). The indiGO controller takes care of the hardware control and performs the measurements even when indiGO itself is not running. Choosing this option the controller will be started everytime indigo is launched (if not already running). After a reboot it will not restart automatically.

With instruments running unattended for a longer period it is recommended to choose the second option (Start controller with Windows) to have a working system even after a power failure when the PC is rebooting. In this case the controller is added to the startup group of Windows and starts automatically after login.

Select **Next >** to continue.



With this screen the NightSHADE evo Controller First Time Initialization Setup is completed. Select „**Finish**“, to restart and initialize the instrument controller now.

4.1.9 First Time Initialization Setup

To configure indiGO launch the software by double-clicking on the indiGO icon.



The first time initialization setup will start.



If the PC is directly connected to the NightSHADE evo via USB choose **“Server”** to install the instrument controller as well as the evaluation part.

In case of a network connection to a NightSHADE evo, select the **“Client”** installation.

Click  to continue.



If indiGo is installed for the first time the first time initialization setup will create a new and empty database.

If indiGO should be connected to an existing database, it has to be attached to this database. Use the browse command to allocate and define a suitable database.

Please note:

IndiGO manages all measuring data in a file-based database.

The indiGO software is able to manage the database on the local computer as well as on a network location. Therefore the database location can be set up to an individual folder.

The benefit of storage on the network is that one has access to the data from several evaluation stations, which is a major advantage in a multi-user environment.

Due to the huge amount of data indiGO does not use a server-based database like Microsoft SQL-server. Therefore the sharing of measurement data through a multi-user environment is not possible. For this reason measurement data should only be edited by one user at a time.

Old databases created with indiGO Version 1.x are not compatible. But data created with these indiGO versions or WinLight can be imported into the new indiGO database using the new indiGO software.

Click to continue.



This dialog box defines the connection address to the NightSHADE evo controller. If the controller is running on the same computer the default settings "LOCALHOST-1024" can be applied.

If the NightSHADE evo controller is running on a remote PC in the network, please set up the corresponding IP-address or the host name like "192.168.1.2-1024" or "IMAGING_PC-1024". The IP port number to contact the instrument controller is always 1024. Host and port number must be separated with a dash "-".



IndiGO offers a user management. Three different modes are available:

- **<Disable User Management>** (not recommended)
- **<Set a Default User Account>**, user will be logged on automatically or
- User must log-in everytime the software starts.

In case no user management is required tick the box **<Disable User Management>**. No log-in is required and no user management is available. If the User Management has been disabled it can be re-activated via the system configuration at any time.



To set up an automatic log on to indigo using a default user account, tick the box **<Set a default user account>**.

By selecting **<Next>** you will be asked to set up a user name and a password. Please note that the created account will have administrator rights.

Click **Next >** to continue.

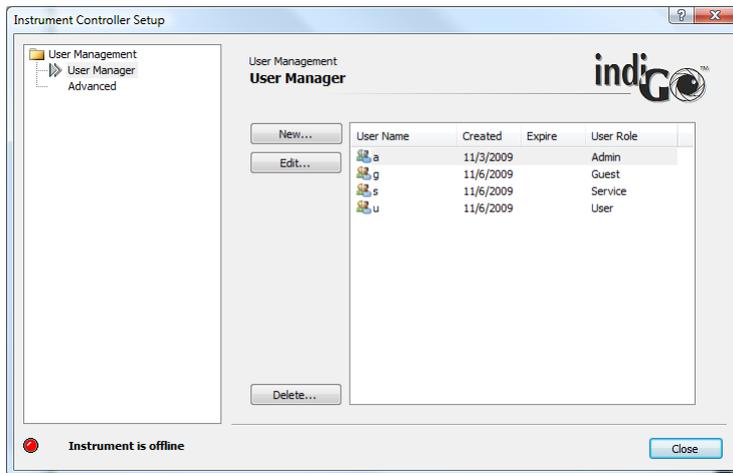
At the end of the configuration a summary page is displayed.



To end the first time initialization setup select **[Finish]**.

4.1.10 User Management

IndiGO supports an integrated user management.



The User Management can be enabled or disabled by the system administrator. Refer to the **<User Management> <Advanced>** settings to enable or disable the user management.

Only admins can access the user management.

Users are organized in groups, so called roles. You can assign one out of five roles for each user.

- Service
- Admin
- Supervisor
- User
- Guest

Service

Service users can change all settings of the system. New users can be created and existing users can be deleted. The access to measurements is restricted through the user management. Other users can be invited to own projects.

Admin

This is the system administrator role. Administrators can change all settings of the system. New users can be created and existing users can be deleted. Administrators have access to all measurements created. Administrators can change the ownership of measurements and manage invitations of all users.

Supervisor

This is the system supervisor role. Supervisors can change most settings of the system like heater temperature, filter management and system status. Supervisors have access to all measurements created by themselves or where they got invitations. Supervisors can change the ownership of measurements and manage invitations of all projects they have been given access to.

User

Users can change any settings necessary to run measurements. Applications and templates available can be created and changed. Users can access and change the filter management. The access to the user management and any controller service functions such as the system

calibration is denied. The access to measurements is restricted by the user management. Other users can be invited to own projects.

Guest

Guests can run measurements in projects where they have been invited to.

4.1.11 Installing Fluorescence Filters

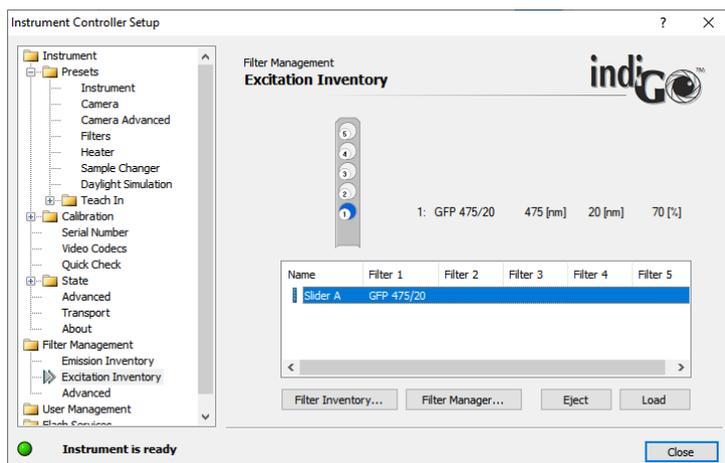
Excitation Filter Slider

The NightSHADE evo is equipped with automatic filter changers for the excitation and emission path. The excitation filter is located on a filter slider, which is accessible after opening the cover on the front panel of the NightSHADE evo. The filter slider can hold five excitation filters with a diameter of 25 mm (1 inch). The 6th position is reserved for luminescence measurements to avoid incoming light through the fibre optic.

Filters placed in the filter slider should be fixed with the snap rings provided with the device. Using excitation filters the inprinted arrow on the filter should match the arrow printed on the fluorescence filter housing. The arrow should face the sample, turning away from the light.



The filter slider can be ejected via the indiGO software (Filter Manager). The filter slider duct is protected by a light blocking plug, which must be mounted after inserting and loading the filter slider.



To eject the slider select .

When loading the filter please observe the correct installation direction (see above). To reload the slider, select . After that the slider is initialized by the NightSHADE evo.

Emission Filter Wheel

The emission filters are located on a filter wheel in front of the lens. A filter wheel can hold up to five filters with a diameter up to 50 mm using specific filter holder inserts.



The positions of the filters are numbered and identified by a respective number printed on the left side of the corresponding filter wheel position.

For luminescence and photographic measurements filter position 6 is used. Using emission filters the inprinted arrow on the filter should face the sample, turning away from the camera.

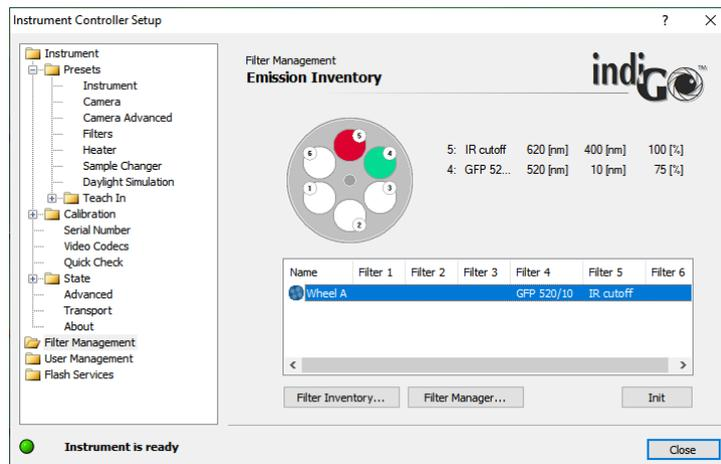
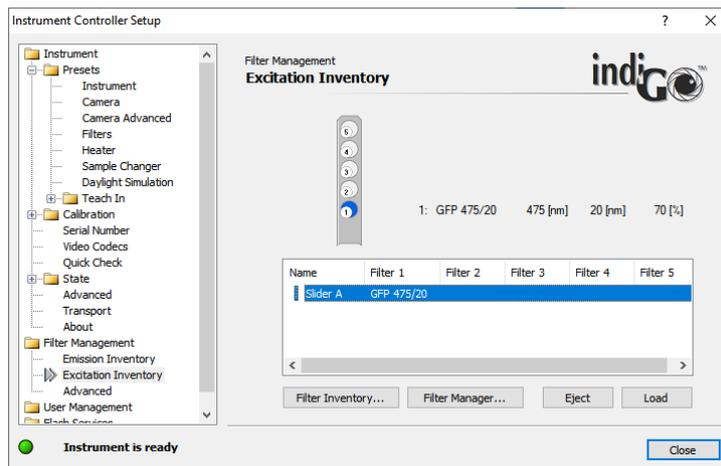
To insert the filter slides switch off the NightSHADE evo, open the filter flap at the front side of the instrument, move the wheel manually to the required position and slide in the filter holder. After moving the filter wheel manually it is required to initialize the wheel again. Otherwise the filter positioning might be wrong. It is recommended to change filters on the filter wheel while the instrument is turned off.

4.1.12 Filter Management

indiGO offers an integrated filter management. The filter management is used to describe the filter load on the excitation filter sliders as well as the emission filter wheels.

The system comes with a set of GFP filters. The recommended excitation position for the GFP exciter (478/20) is position 1.

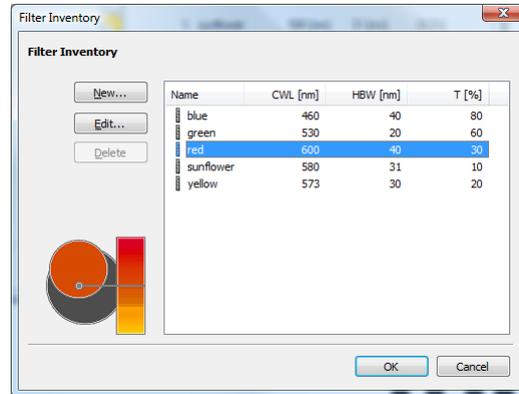
The GFP emission filter (520/10) is recommended to be placed in position 4. The IR-cutoff filter that comes with the NightSHADE evo is recommended to be placed in position 5



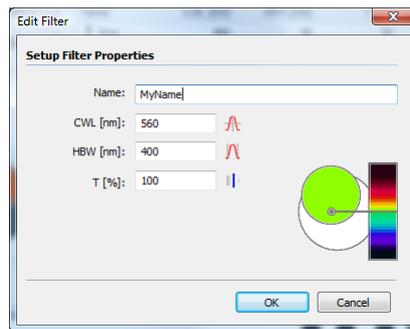
4.1.13 Filter Inventory

The filter inventory is used to collect the available filters. indiGO governs two filter inventory lists. One is used for excitation filters, the other for emission filters.

The inventory lists can be either reached via **<Excitation Inventory> <Filter Inventory>** for the excitations filters or via **<Emission Inventory> <Filter Inventory>** for the emission filters.



To add a new filter to the inventory list, select the **<New>** command.



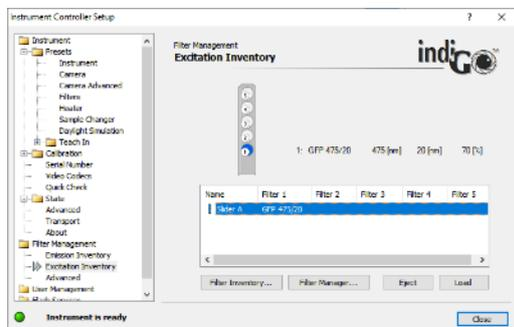
Each filter is described by the following properties:

- The name of the filter
- The center wave length of the filter
- The half band width of the filter (FWHM)
- The transmission rate

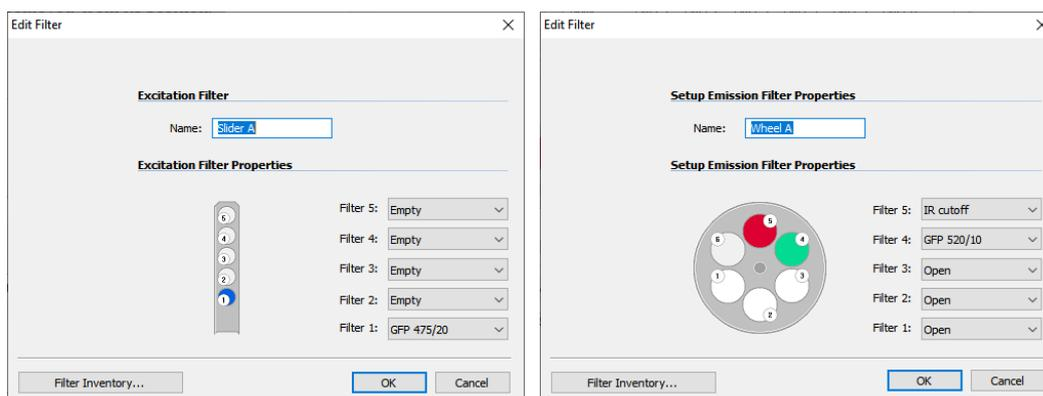
4.1.14 Filter Manager

Once a filter has been added to the filter inventory list it can be used for an excitation filter slider or an emission filter wheel.

The filter manager can be reached via **<Excitation Inventory> <Filter Manager>** for the excitations filters or via **<Emission Inventory> <Filter Manager>** for the emission filter.



To add a new excitation filter slider or an emission filter wheel to the manager select **<New>**. To change or edit the load of a slider or a wheel select **<Edit>**.



The loading can be changed using the filter dropdown list **<Filter N>**. Please keep in mind, that only available and unused filters are displayed in the droplist. Before you can assign a new filter to a position, either a new filter has to be added to the **<Filter Inventory>** or an existing filter has to be removed from its position.

4.1.15 Filter Manager – Backup and Restore

All data from the filter manager can be exported from the system and stored to an archive as a backup using the command **<Filter Management> <Advanced> <Export Filter Inventory>**.

A further created backup can be imported using the command **<Filter Management> <Advanced> <Import Filter Inventory>**.

4.2 First Acquisitions

4.2.1 Quick Start Glossary

Projects, Applications and Identities:

All parameters necessary to run a measurement are defined in an application.

Individual samples, called identities, can be organized in a project. Each project is connected with an application.

To perform a measurement, select an identity and the system automatically refers all settings of the project and the correlated application.

After the measurement is accomplished all data will be automatically transferred to the pre-selected identity of the project.

All acquired data is stored and organized as a structured tree. The main nodes are the projects containing identities, containing measurement data.

Example:

- Project 1
 - Identity 1
 - Measurement 1
 - Measurement n
 - Identity n
 - Measurement 1
 - Measurement n
- Project n

Projects and Applications

Applications consist of a pre-measurement, a measurement and a post-measurement.

When a project is created a mirror copy of the used application is saved with that project. This means that changing the original application template later on will not affect the project's parameters. Consequently changing parameter settings in the project (e.g. exposure time) has no influence on the original application template.

On the other hand before each measurement the application settings of a project can be modified. The last modification is always saved with the project. Settings used for each specific measurement performed can be found in the report.

Projects and Users

Projects can be associated to users. When created, a project belongs to a specified owner. The owner has the possibility to allow other users to access the project and to run measurements on identities associated with the project. To allow other users the access to a project, the owner has to create an invitation.

4.2.2 Launching the software

Depending on the system configuration it might be required to log-in after the indiGO has been started. When no other users are defined the standard log-in is as follows:

Log-in: Admin
Password: Admin

After log-in the main window of indiGO is shown:

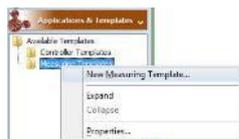


4.2.3 Defining a Measuring Template

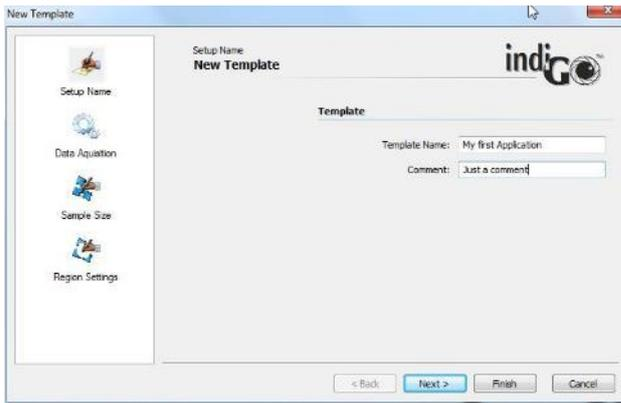
indiGO defines different measurement methods (e.g. luminescence measurement, fluorescence measurement) by an measuring template. In each template the measurement mode, the exposure time, the measurement sequence and the filter settings, (if available), are assigned.

Administrator access rights are necessary, to set up a measuring template. If this is not guaranteed the function to define a measuring template is not available to rule out faulty operations.

To set up a new template refer to **<Applications & Templates> <Measuring Templates>**. Select **<New Measuring Template>** from the **<Measuring Templates>** context menu.



First enter a name after indiGO has started the wizard. In addition a descriptive comment can be added. Due to Windows file and path name conventions special characters such as `< > ? " : | \ / * .` are not allowed.

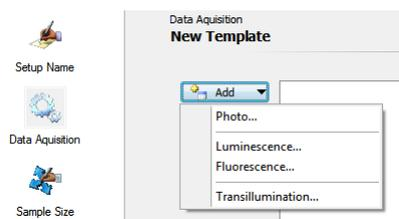


Select **Next >** to define the measurement workflow.

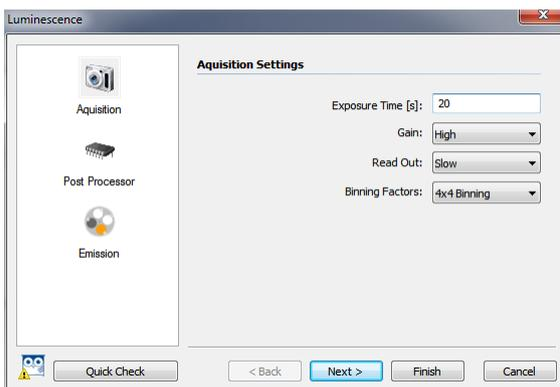
Four different kinds of measurements can be performed. Please note that depending on the system configuration some of the selections might not be available (e.g. due to a missing fluorescence illumination module).

Example:

To define a luminescence measurement, select "luminescence" from the dropdown list after clicking on **Add**



The Luminescence image wizard will pop up.

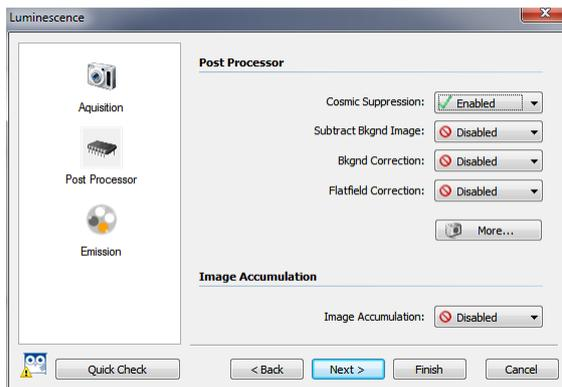


- **Exposure Time**
is the length of the time the CCD chip is exposed to the sample
- **Gain**
is an internal conversion factor. Low Gain offers high dynamic range while high gain gives higher intensity numbers but limited dynamic range. When using the NC100 camera the difference in signal between low and high gain is about factor 4.
- **Read out**
defines the speed to read out the camera. High speed is recommended for photographic images and high intensity measurements such as fluorescence since a higher

background noise will be obtained. Low speed is ideal for highest sensitivity luminescence measurements. Be aware that with low speed the camera readout can take up to 15 seconds.

- **Binning Factor**
is the combined readout of neighbouring CCD pixels. The result is an increased signal and thus an improved sensitivity and a better signal-to-noise ratio, however, at the cost of loss in spatial resolution.

Select **Next >** to get to the post processing definition.

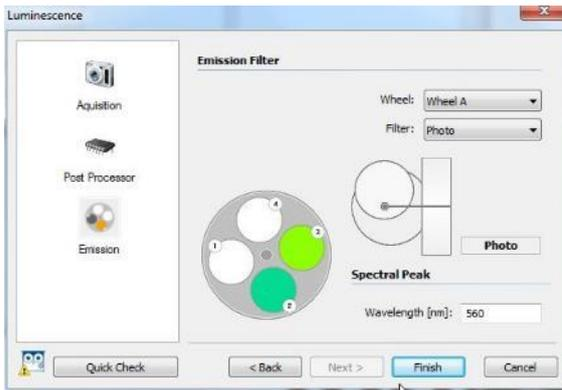


The Post Processor page offers additional image processing features which can be performed after the exposure to increase the image quality.

- **Cosmic suppression**
is an imaging filter to get rid of small high intensity pixels caused by cosmic ray during the exposure.
- **Subtract background image**
This function will subtract an individual background image from the taken image. The mode "Pre Image" or "Post Image" defines if the background image will be taken before or after the real image using the same settings. While taking the background image, the camera shutter is closed. The image background subtraction eliminates camera background pattern or memory images.
- **Background Correction**
subtracts a specific amount of counts from the current image. The background is calculated using the lowest intensities on the image.
- **Flatfield Correction**
corrects an uneven image caused by the vignetting of the lens
- **Image Accumulation**
With activated image accumulation more than one image is taken and added on top of the previous one. This allows a smoother background and offers a better cosmic suppression (if enabled). Up to 9 images can be added.

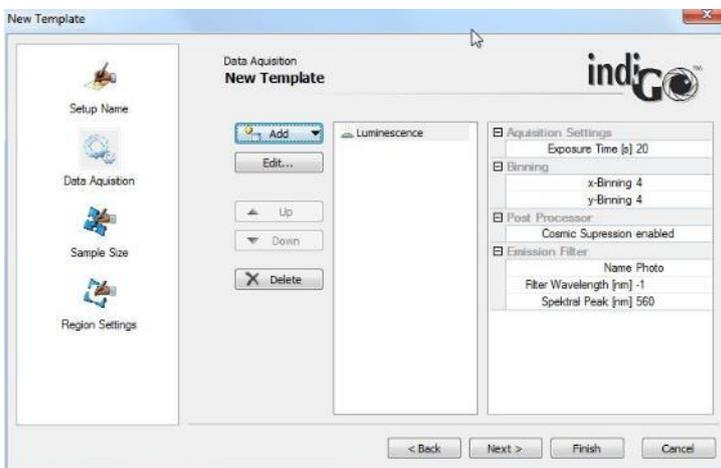
Enable the cosmic suppression to avoid spots caused by cosmic rays during longer exposures. Clicking on the **More...** button will open the parameter box to adjust the image processing filters. For save cosmic suppression use about 50% strength. Increasing this value will increase the risk that real small spots will be treated as cosmics.

Since luminescence is measured no filter is necessary. That requires the selection of the empty / photo position in the emission filter settings. To ensure a proper energy calibration check the wavelength mentioned in the input box (560nm for firefly luciferase).

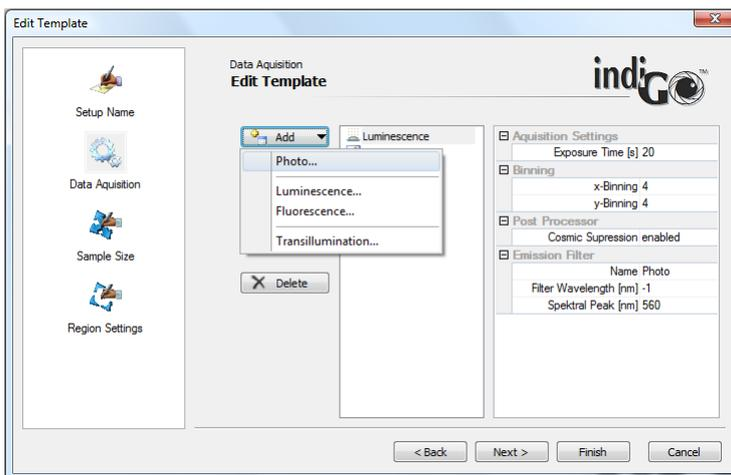


To close the luminescence parameter wizard select **Finish**.

The summary of the new created luminescence measurement is shown in the right column:

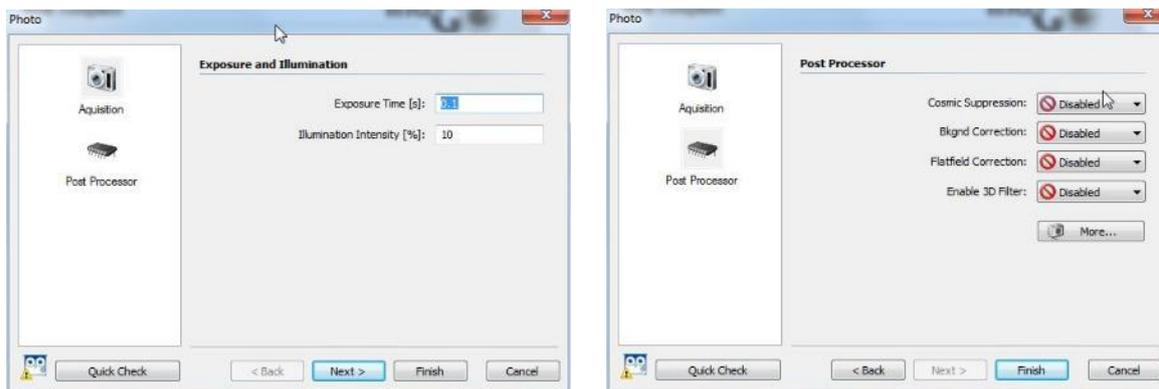


To add a photographic exposure to the list of measurements select **<Add>**, **<Photo>**.

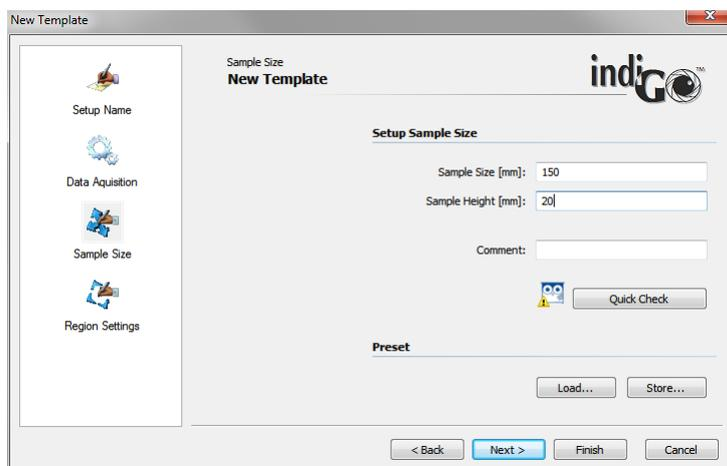


The default parameters (100ms exposure time, 10% illumination intensity) are suitable for most applications. Disable all post processors in the corresponding section. The “close iris” feature improves image quality (depth of focus). This feature is not available with in the NightSHADE evo

To close the photographic parameter wizard select **Finish**.



For the sample size definition dialog box select **Next >**.



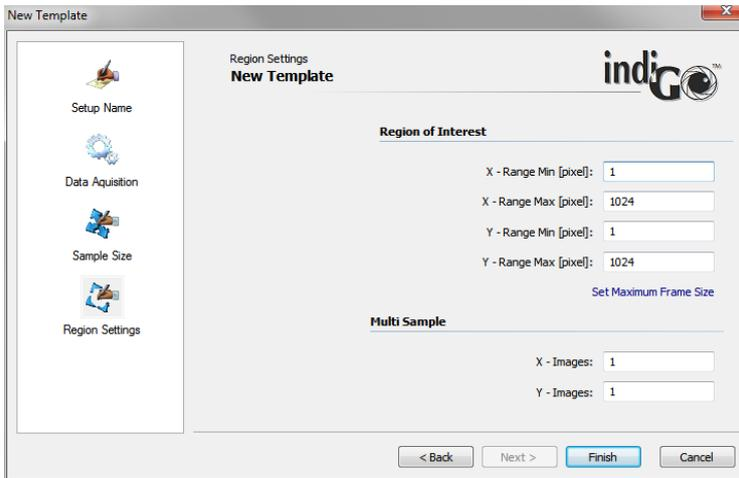
You can either select an already existing format using the load button or define a new sample format. The following parameters are required:

- Sample Size:**
 The sample size refers to the image size (width) in mm. To make sure that the sample will be measured completely it is recommended to define the image size a little bit larger than the actual size of the sample (hier ist bei der Shade doch egal, was man einträgt, sollen wir das nicht ändern?)
- Sample Height:**
 The sample heights refers to the focus height. If you are measuring flat samples (gels, blots) located on the base plate the sample height is 0 mm. If elevated samples are measured (e.g. animals or plants) the sample height has to be set to a relevant range (e.g. height of a special leaf). If you are analyzing animals, it is advisable to use half the height of the animal as sample height. If you are working with sample holders or other extra parts (e.g. transilluminators), you have to take their height into account when calculating the sample height.
- Comment:**
 You can enter a comment on the sample format which describes the format.

The **Quick Check** can be used to start an acquisition using the entered values. The result is shown in a preview where you can either check the focus or whether the entered sample size is adequate for the acquisition of the sample.

If the focus is not OK it can be adjusted by manually adjusting the focus ring at the lens.

To go to the next page select [Next >](#).



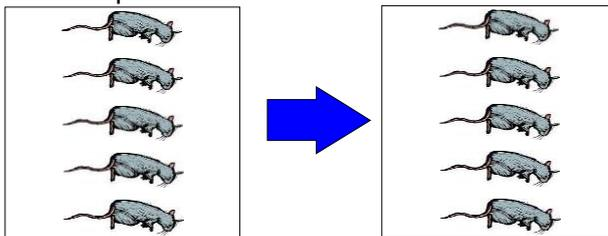
On this page you can define the setting for the “Region of Interest” as well as the “Multi Sample” parameters.

By using the “Region of Interest” a section of the CCD chip can be defined to get faster framerates. By default the whole chip area from 1 to 1024 for X and Y is used

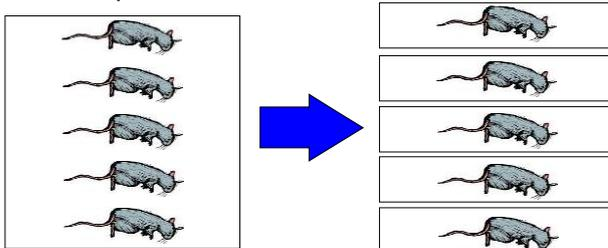
The command “Set maximum Frame Size” will adjust the frame size according to the selected binning factor since some cameras cannot handle full frame at different binning factors.

The MultiSample feature allows the possibility to split the image into image sections. Each section is treated as a single sample identity. This feature might be useful for animal imaging in case that more than one animal at a time is measured. The X-Images and Y-Images define the amount of samples in X and Y direction. This feature is limited to 5 images in each direction (max 25 sample-ID’s per measurement). Be aware that working with individual images will also require to run individual reports which is not useful for some applications.

Multisample X1 and Y1



Multisample X1 and Y5



In case of multi sample the required identities will be defined before starting a measurement.

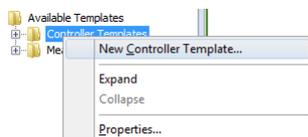
To close the application wizard select .

4.2.4 Defining a Controller Template

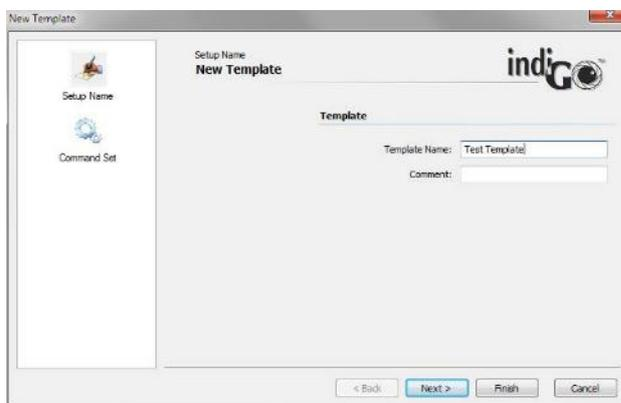
When creating a controller template the workflow is about the same as with a measuring-template. . But instead of defining measurement settings controller templates are used to control additional devices such as the LED panels in order to simulate environmental parameters (e.g. day & night cycle). Furthermore control templates are used to define pre- and postmeasurement settings such as “Daylight off and on” which allow to interrupt the light simulation and go back to its original setting after the measurement.

A controller template consist of a list of single commands. This list will be splitted in its single commands when executed using the so called scheduler. Therefore it is possible to perform measurements inbetween

To set up a new controller template refer to **<Applications & Templates> <Measuring Templates>**.Select **<New Controller Template>** from the **<Controller Templates>** context menu.

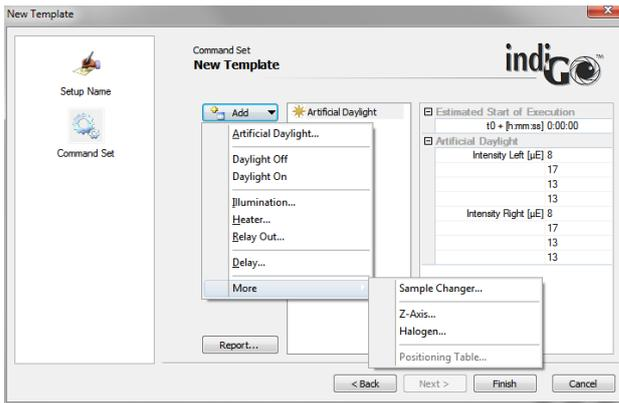


indiGO starts the wizard. In the first windowthe controller template can be named and a descriptive comment be added

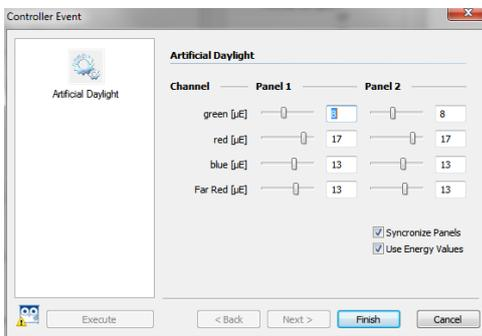


Select  to define the control workflow.

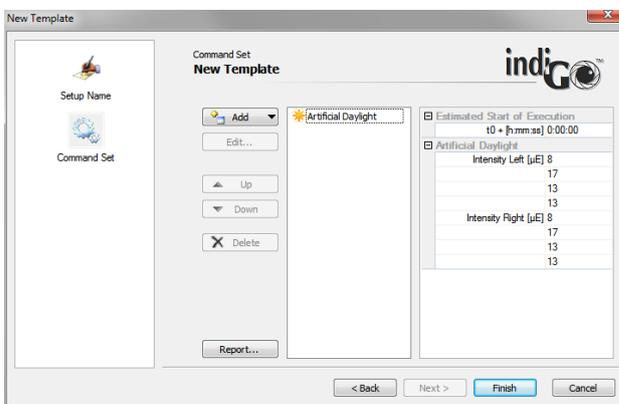
Different kinds of control actions can be performed. Please note that depending on the system configuration some of the selections might not be available (e.g. due to a missing LED illumination).



To turn on the LED illumination (Option) select **<Artificial Daylight>** from the list. Now the intensities of the different wavelength can be adjusted



Adjust the required intensities and click on **Finish** to accept the parameters. The control command shows up in the work flow list.



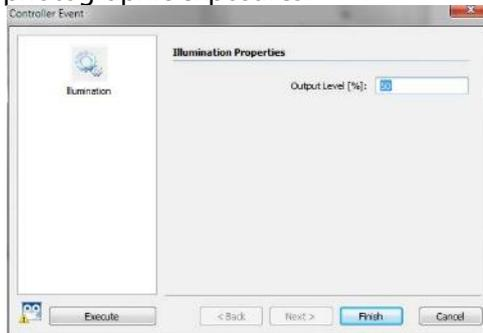
After setting all commands, close the dialog box and save the template by clicking on **Finish**

- **Daylight Intensity (option)**
This command sets the LED illumination panel to a defined intensity.
- **Daylight Off (option)**
Turns off the LED Panels. The previously used intensities are temporarily saved for future use (when using “Daylight On”). Should be added to a project as Pre-Measurement action to make sure the artificial daylight is off while measuring.
- **Daylight On (option)**
Turns on the LED panel using the last used intensities saved by the “Daylight Off” command. Should be added to a project as Post-Measurement action to make sure

the artificial daylight is turned back on to the previously used intensity after a measurement.

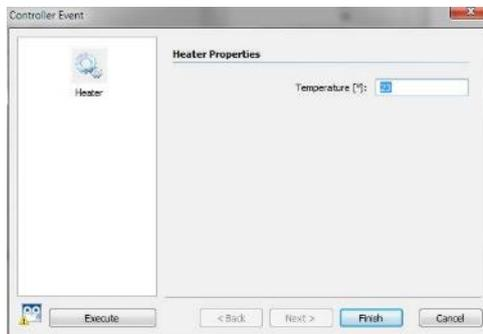
- **Illumination**

Controls the intensity of the white or infrared (option) LEDs in the dark room for photographic exposures.



- **Heater**

Sets the target temperature of the temperature controlled measuring table (Option).



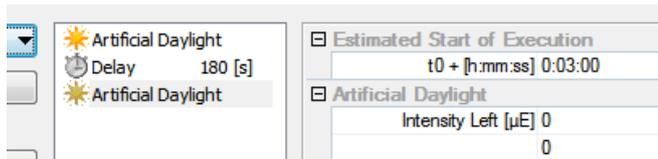
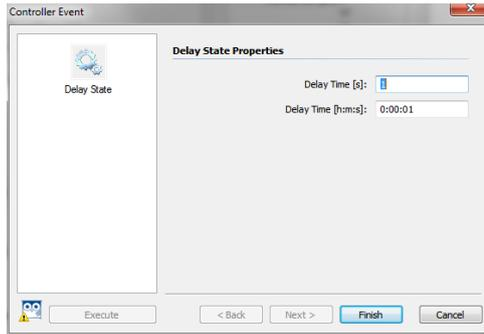
- **Relay Out**

Controls the power socket inside the dark room (Relay 1). Relay 2 is not used with this instrument.

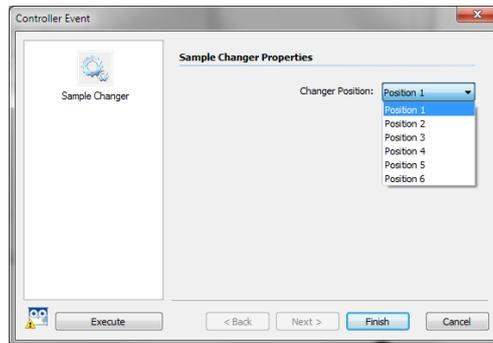


- **Delay**

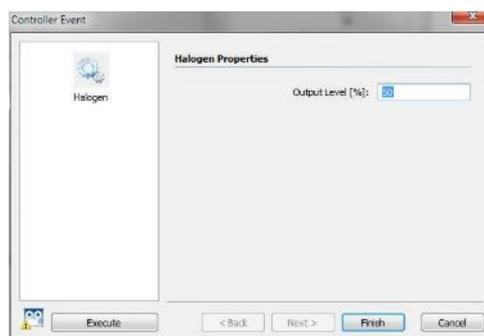
Defines a waiting time within the control sequence. When inserting a delay command the relative time of the control sequence (t0) will increase. The delay time can be inserted either in seconds or in h:m:s format (hour : minute :second).



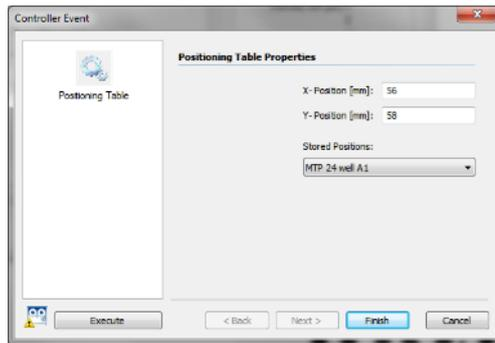
- **Sample Changer (option)**
Moves the sample changer to a defined position.



- **Halogen**
This command turns on the excitation light of the fluorescence illumination..

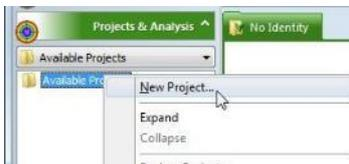


- **Positioning Plate (available with older NightSHADEs only)**
The positioning plate is an optional table to move the sample in X and Y direction. The required position can be entered in mm or a predefined position can be selected from the dropdown list.



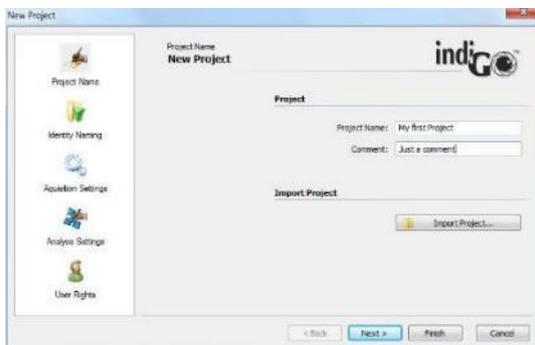
4.2.5 Defining a Project

To define a new project refer to **<Project & Analysis> <Available Projects>**.
Select **<New Project>** from the **<Available Projects>** context menu.



Projects are used to define non measurement-specific parameters, like user access rights, scale settings, etc.

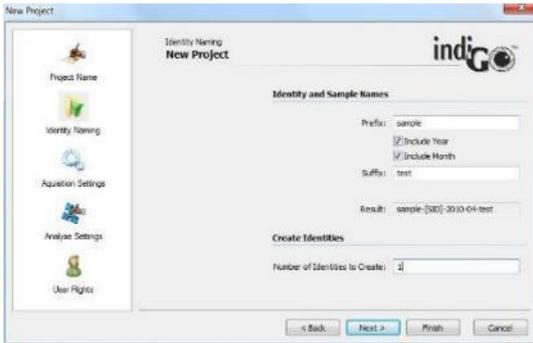
To create a new project indiGO starts the wizard:



The following parameters are necessary:

- **Project Name:**
Enter the project name under which the project is to be saved. Due to Windows file and path name conventions special characters such as < > ? " : | \ / * . are not allowed.
- **Comment:**
Enter a comment which describes the project.

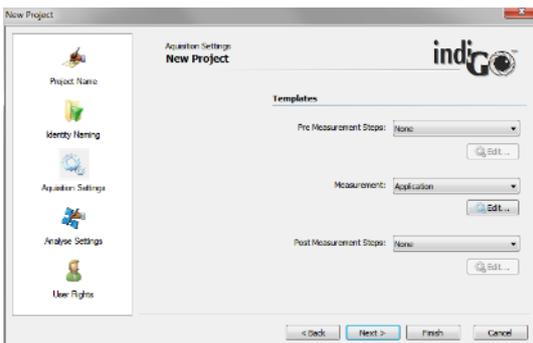
To go to the next page select .



On this page the number of sample identities and the sample names can be defined. A prefix, the current year and month, a 3 digit sample ID counter and a suffix can be specified if required. The resulting sample ID name can be seen under **“Result”**. The box **“Number of Identities to Create”** defines the amount of sample identities which will be created. More ID’s can be added to the project later, if necessary.

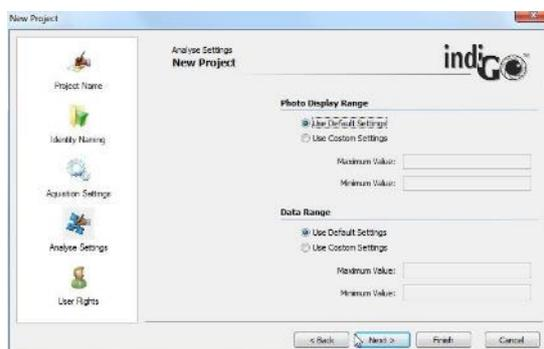
To go to the next page select [Next >](#).

Choose the **“Acquisition Settings”** used for the project.



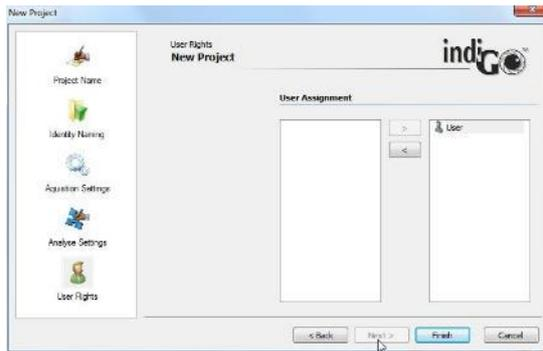
a **“Pre Measurement”** and a **“Post Measurement”** can be defined to be executed before and after the measurement. Under **“Measurement”** select the required measuring template.

To go to the next page select [Next >](#).



The display ranges for the b/w photographic and colored signal (data) image are defined on this page. If **“Use Default Setting”** is selected the system will calculate default values based on the photograph and the measured data. Individual settings for the minimum and maximum values can be selected in the **“Use Custom Settings”** mode.

To go to the next page select [Next >](#).



On the last page users can be assigned to the project. The projects creator automatically receives ownership of the project. Other users can be added to the project. To create new users refer to the indiGO user manager.

All defined users are listed in the left column and can be added to the project using the add command (). All users assigned to the project are listed in the right column; Users can be removed using the remove command ().

Users assigned to the project are authorized to accomplish measurements for the project. Users, defined as **“users”** can't change the measurement settings for the project. For users defined as **“guests”** the access to measure is denied.

To complete the project definition wizard select .

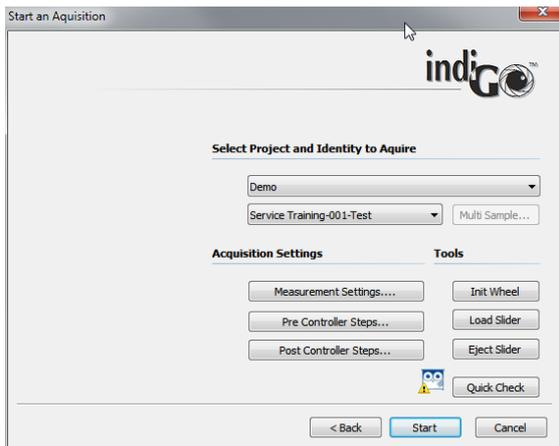
4.2.6 Starting an Acquisition

The system can carry out measurements on demand or scheduled using a time table.

First we will show here how to create an acquisitions on demand. Further details about how to use the scheduler can be found at the end of the chapter.

To start an acquisition refer to **<Acquisition & Scheduling> <Start an Acquisition>**.

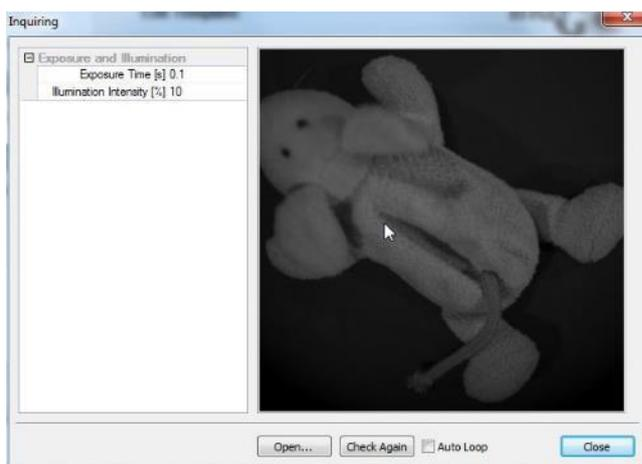
The acquisition wizard will be shown:



The project and identity might be pre-selected. Otherwise select the project and the identity from the provided drop down lists.

By selecting **Quick Check** a photographic preview of the sample is created using the project settings

This preview might look as follows.

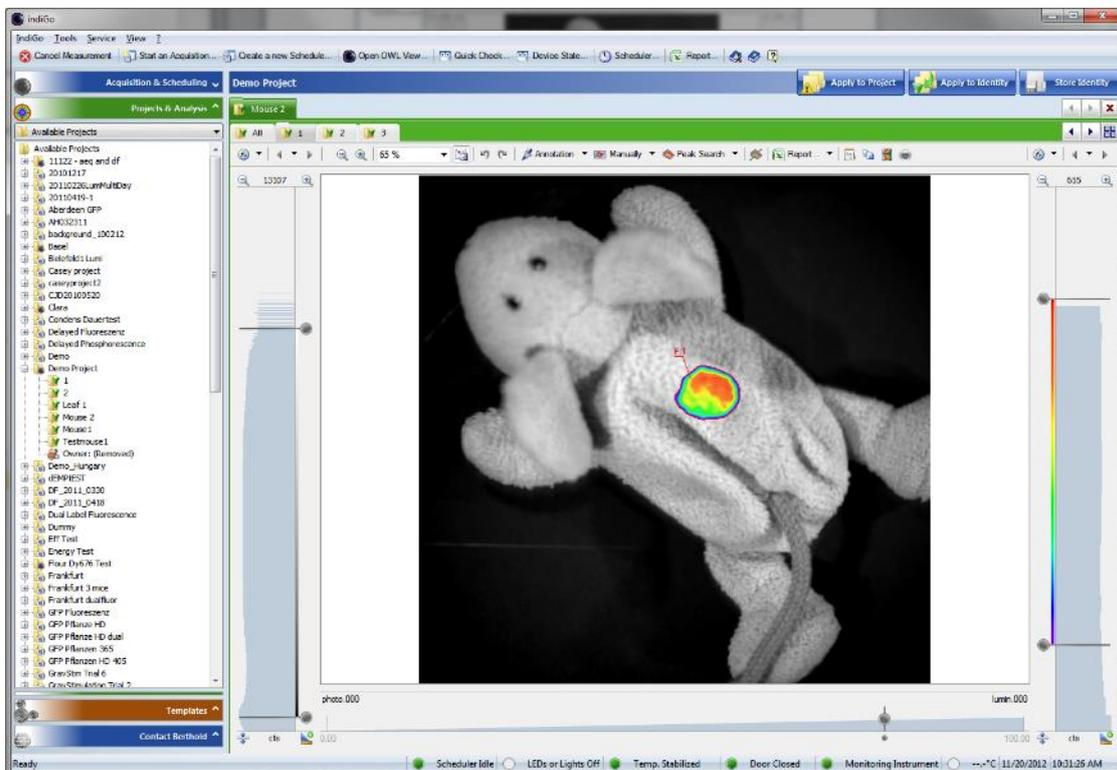


If the "Auto Loop" checkbox is activated the camera will take an image continuously. This might be useful for sample positioning or when adjusting the focus at the lens.

To focus the image just turn the manual focus wheel located next to the camera on the front left side. Note that it might take some time until the new image will show up on the screen.



To start a measurement select . Upon completion of the measurement, the data is automatically stored. The data is displayed by double clicking on the sampleID.



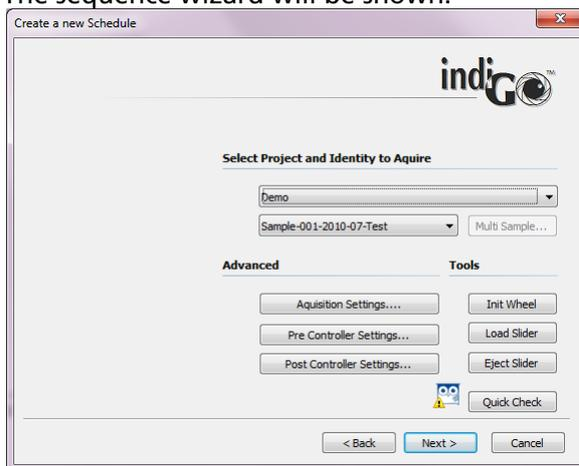
The measurement consists of a photographic image in grayscales and the luminescence image in pseudo color.

4.2.7 Image Sequences

In addition to single measurements indigo also offers automatic sequence measurements. These kind of measurements allow automatically triggered measurements in a predefined schedule.

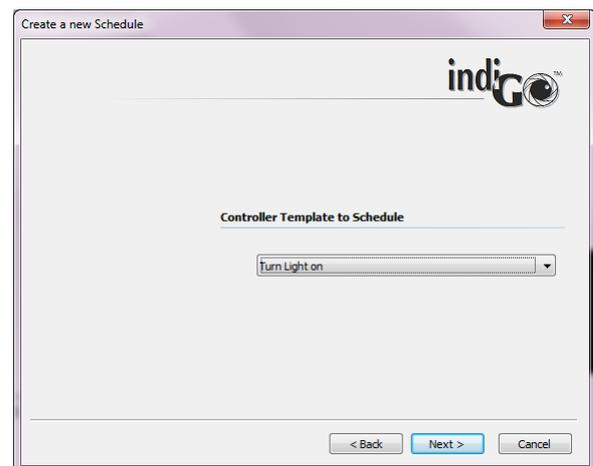
To start a sequence measurement refer to **<Acquisition & Scheduling>**, **<Create a new Schedule>**.

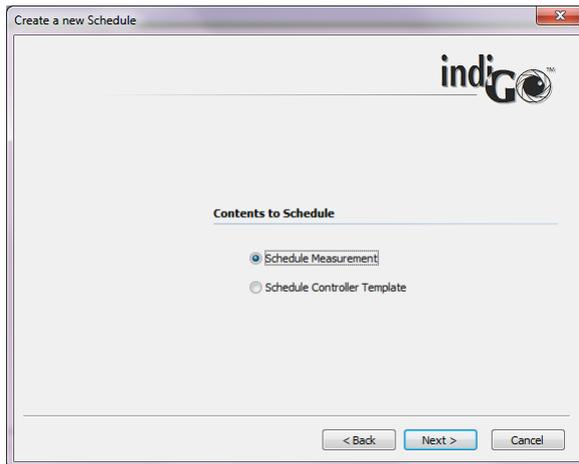
The sequence wizard will be shown:



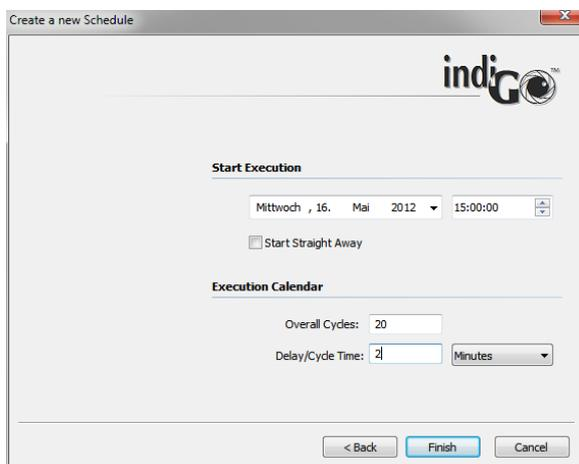
First of all specify the type of scheduled event. To set up a scheduler for measurements choose "Schedule Measurement", to set up an schedule for controll actions (lights on, mains switch, etc.) choose "Schedule Controller Template".

If choosing the measurement mode, first a project and sample ID has to be selected. In case of the controller template mode, the corresponding controller template has to be selected from the list.





Clicking on [Next >](#) leads to the following dialog box.

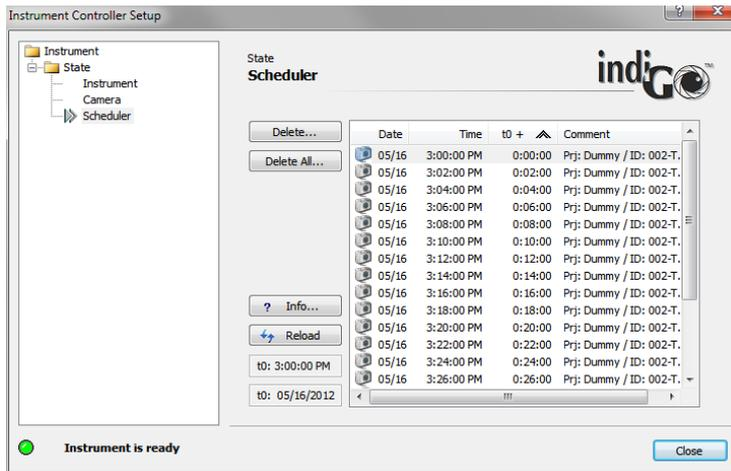


In this dialog box the starting time for the measurement- or control sequence can be defined. Alternatively "Start right away" can be selected by checking the box. This will enable an immediate start when [Finish](#) is clicked. All other events will be relative to this starting time.

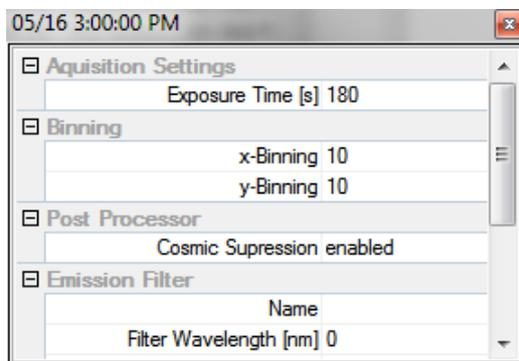
In the "Execution calendar" the total number of measurements / sequences as well as the interval time can be specified. Clicking on [Finish](#) will close the sequence wizard and the sequence is handed over to the scheduler.

Now the sequence is active and the instrument controller is taking care of the measurements. It's possible to evaluate other measurements or even exit indigo. Only the instrument controller must be running. This is the case as long as the indigo icon is visible in the taskbar.

To check the scheduler click on [Scheduler...](#). The scheduler will show up displaying all queued jobs with absolute and relative times.



Job details can be seen by doubleclicking the corresponding item in the list.



To delete a single job, click on the item to select it and click on . To delete the whole Schedule click on .

4.3 Image Processing and Evaluation Functions

4.3.1 Load Images

During the first acquisition images have been created and stored in the project folder. To re-open the previously taken images refer to the project folder first and select the appropriate identity. All identities associated to the project are located in the project folder. The measurements including the images are stored under the context of the identity and can be opened by either double-clicking on the identity or using the identity context menu.

Please note:

The contents of the project folder shown depend on the user account, based on the indiGO user access management.

The identity number is displayed on top of the data view. If several sample IDs are open, a cardfile tab for each identity is displayed, each cardfile tab has its own color.



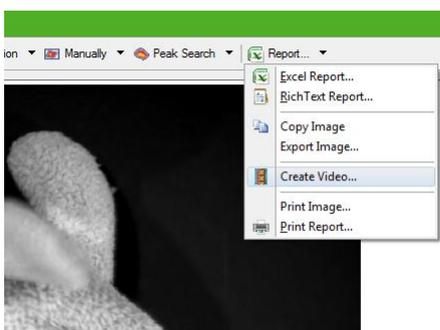
Below the identity number a row of index cards is shown containing the measurements acquired. The last acquired measurement is shown on the very right.

Using the   arrows you can scroll back and forth within the measurements or the identities.

To display a complete overview of the images, the so-called **"thumbnail view"** select .

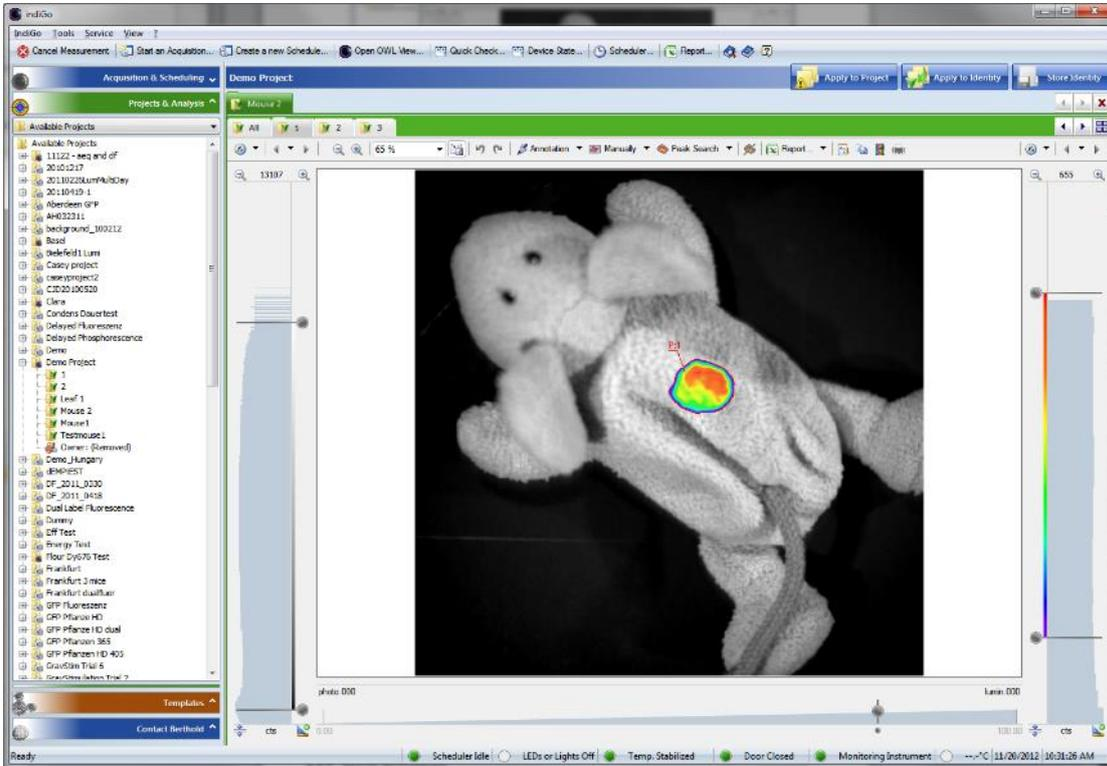
To close the identity select .

Using the **<create movie>** command in the **<Report>** menu a movie of all images in the current identity can be created.



4.3.2 Image Window

After an acquisition or the opening of an identity the measurement and the photographic image are displayed. By default the photographic image is overlaid by the measurement or signal image using pseudocolors.

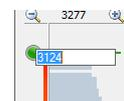


At the left margin of the screen you see the intensity scale and histogram of the photographic image, together with the associated sliders. These sliders can be used to change the brightness intensity of the photographic image. The top slider defines the upper brightness value, whereas the bottom slider defines the lower brightness limit.

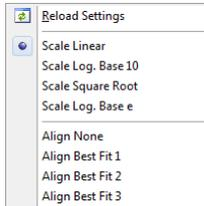
At the right margin of the screen you see the intensity scale and histogram of the measurement, together with the associated sliders. With these sliders the scaling of the pseudocolor signal image can be set. The top slider defines which intensity value has been assigned to the highest color value. Values above this threshold are not displayed (transparent area). The bottom slider defines which intensity in the signal image is displayed using the lowest color value. All values below this threshold are not displayed.

To modify the scale move the slider to the position required. Moving the mouse to the center of both sliders a center mark is displayed which enables you to move both sliders at once.

To enter a certain value rightclick on the slider and enter the new value in the input field.



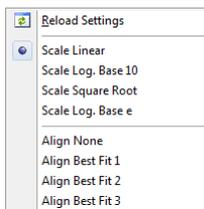
The system offers the possibility to zoom the histogram by using  3277 . The maximum range of the histogram is displayed on the zoom toolbar. By selecting  the histogram can be set to fullscale. With  the displayed unit can be changed from cts to %, nW or ph/s.



The scaling method can be changed via the context menu which can be invoked by clicking on the arrow next to the  icon.

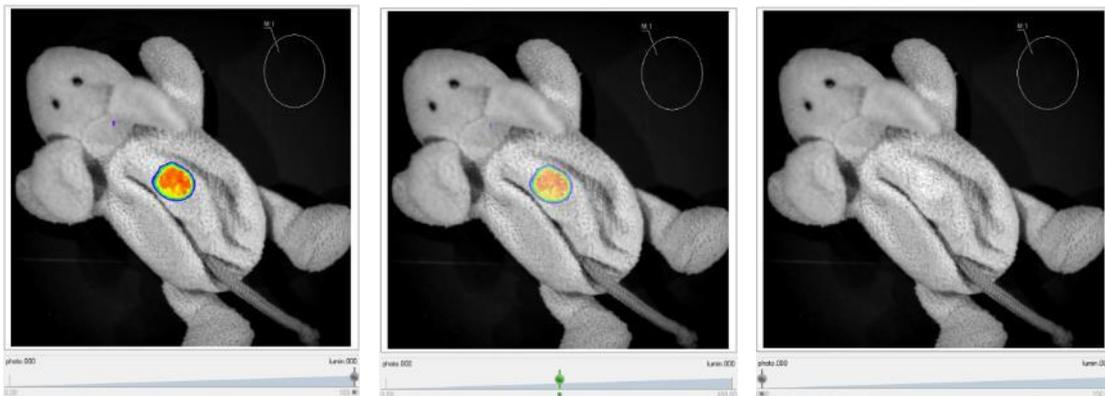
The following options for the scaling are available:

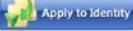
- Linear scaling
- Logarithmic scaling (base 10)
- Square root scaling
- Logarithmic scaling (base e)



In addition different autoalignment preset features (Best fit 1 to Best fit 3) are listed to adjust the min / max value based on the histogram.

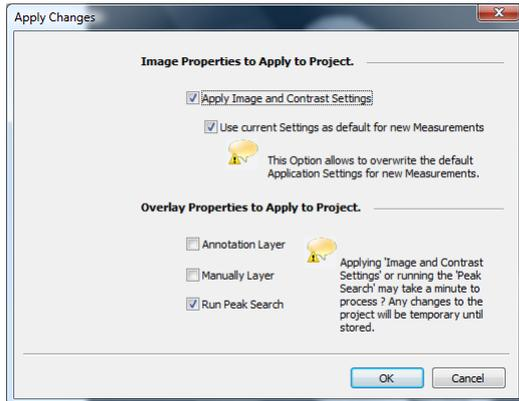
At the bottom margin of the screen a transparency slider is available. This slider can be used to define the transparency of the overlay. If the slider is in the 100% position, the overlay is completely opaque, whereas the overlay is fully invisible if the slider is set to the 0%.



Changes of the intensity scaling or the color threshold can be applied easily to all other measurements within the current identity using the  command. This simplifies the post run changes on measurements, as it is not necessary to modify every single image.

If the current changes should be applied to all identities measured within the same project, the  command can be used.

The command has to be confirmed and can be parameterized using the dialog shown below.



The following options are available:

<Apply Image and Contrast settings> will apply the image and contrast settings from the current image to all other measurements in the project.

If **<Use current Settings as default for new Measurements>** is selected, these settings are also valid for all future measurements in the project.

The selections **<Annotation Layer>** and **<Manually Layer>** will copy the areas and annotations to all other measurements within the project.

The **<Run Peak Search>** option will perform a peaksearch on all other measurements in the active project using the current settings.

The process will be started by selecting .

This opens all data of the current project (Apply to project) and applies the changes. To get back just close the data without saving.

Please note:

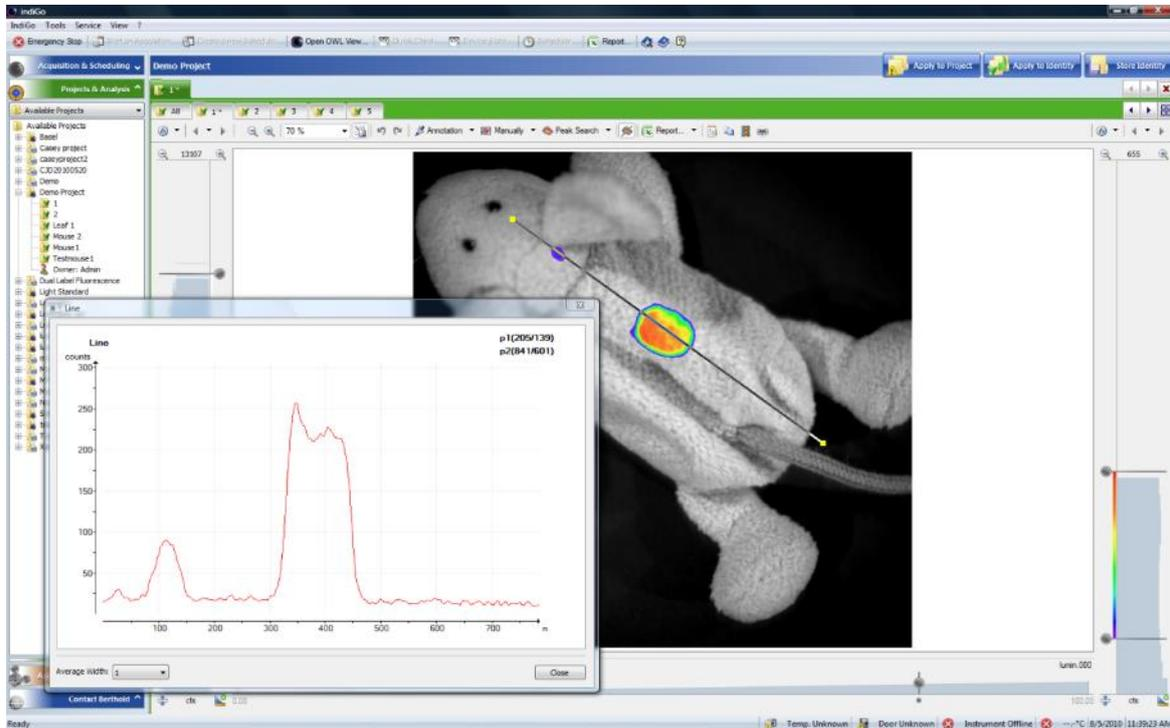
Depending on the selections and the amount of data this function may take some time to accomplish.

4.3.3 Line Plot

For further analysis of the measurements it might be helpful to create a 1D-line plot of the measured data. This function is recommended to evaluate the measuring data.

To use this function select .

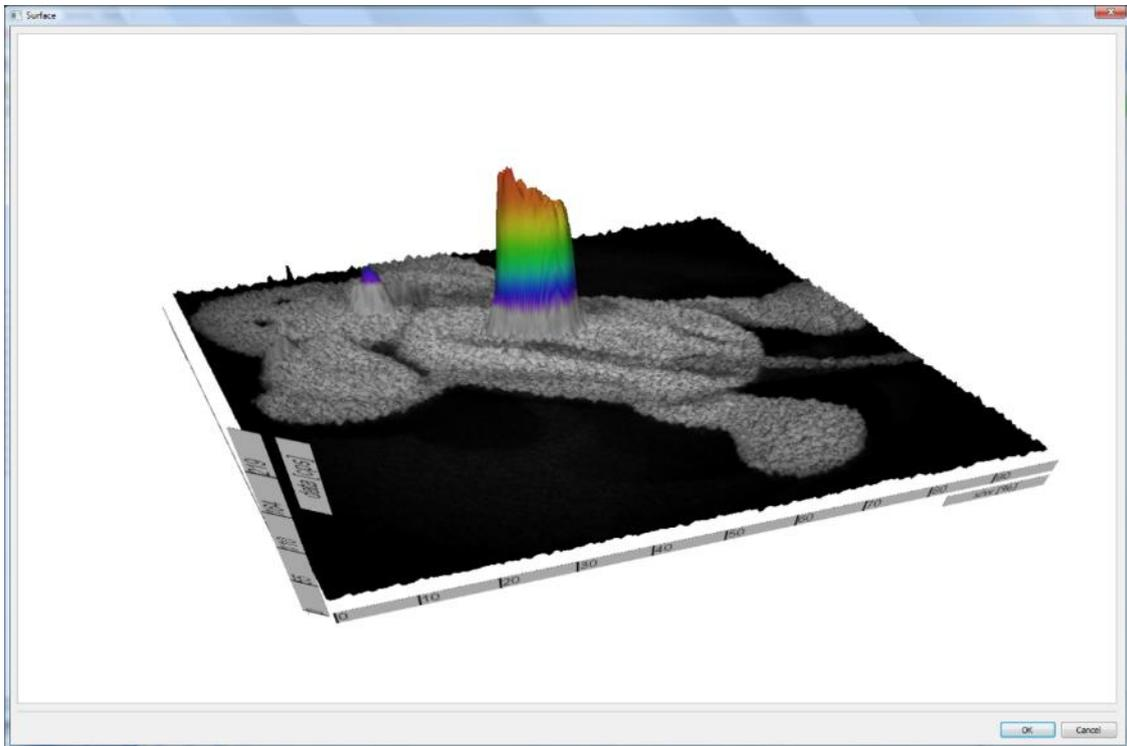
An additional window is opened where the line profile created from a given lead is displayed.



The pixel intensity along the selected path is displayed as a line diagram. The average width of the lead can be selected. Moving the lead updates the line profile.

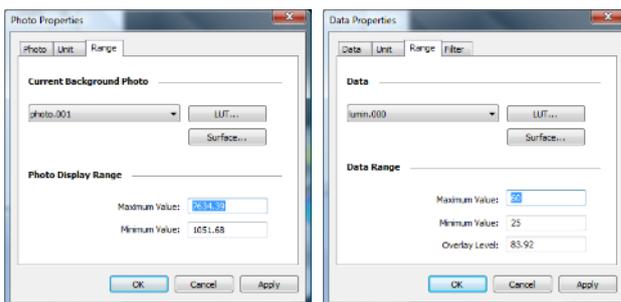
4.3.4 Surface Plot

To get a 3D surface plot select **<Surface>** from the **<Report>** menu. The surface can be used for 3D visualization of the measured data.



To create a 3D surface plot from the photographic image only please refer to associated  command. The same applies if you want to create a 3D surface plot from the measurement data only.

From the dialog **<Photo Properties>** or **<Data Properties >** select **<Range>** and start **<Surface>**.



To change the perspective of the 3D-view pick the graph and move the image to the position required. To zoom in use the mouse wheel.

To animate the surface double-click on the graph.

Please note:

The windows displaying the surface can be resized.

4.3.5 Peaksearch

Once the image has been loaded and the color view has been set up, you can start to analyse the evaluation.

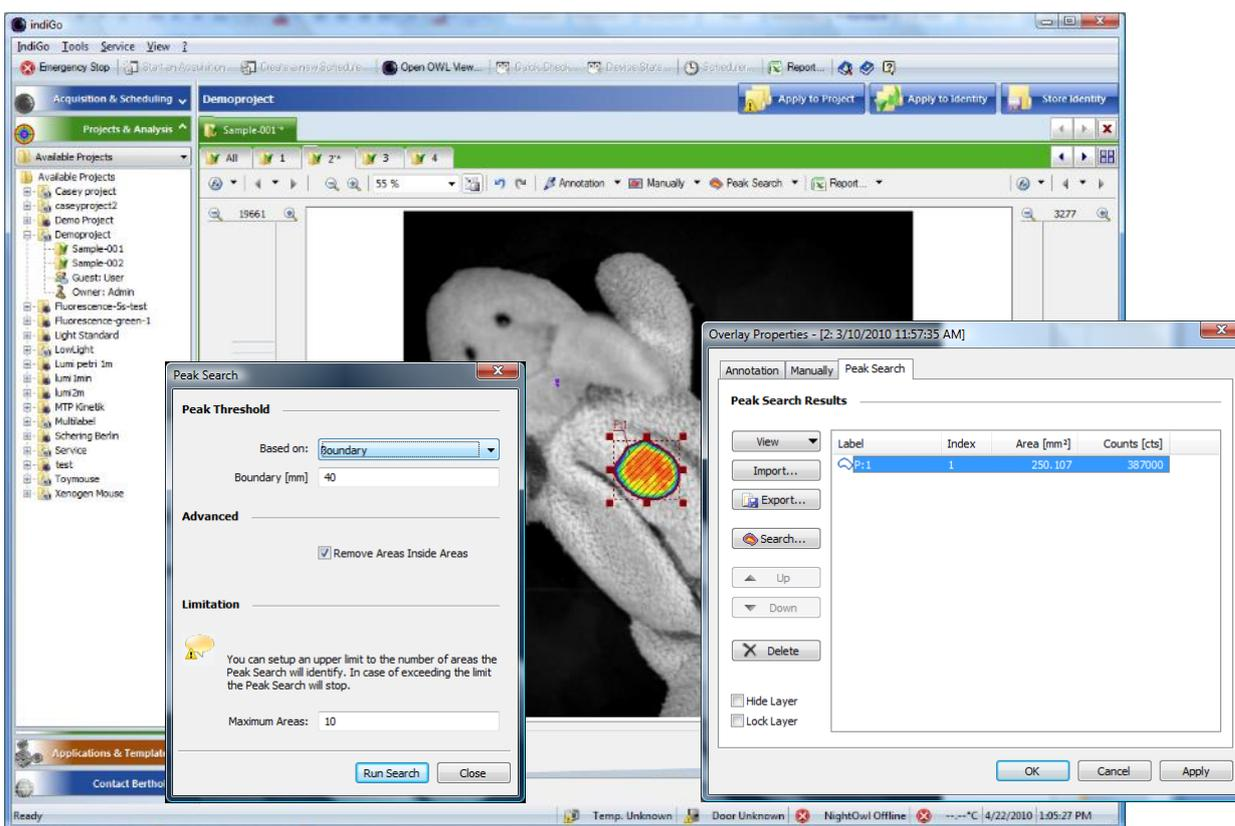
In the first step the desired areas has to be marked; now indiGO can calculate the intensity values.

The easiest method to find evaluation relevant data is to use the peaksearch.

Using the peaksearch all areas lying above a threshold value will be marked. With this method it is possible to evaluate different measurements under exactly identical evaluation conditions.

To start the peak search select  Peak Search .

The displayed overlay properties window gives an overview of all automatically found peaks in the current image.



To run a new search select  Search...

The following parameters can be edited:

- **Peak Threshold - Based on**

Using a droplist different search algorithms are available:

- **Area**
The area of every possible spot is analyzed and if the area is larger than a size specified in a parameter box below, the spot is validated as a peak. This is useful to get rid of smaller, unwanted spots.
- **Boundary**
The most common algorithm is the boundary discrimination. The boundary of every possible spot is calculated and if the boundary is larger than a specified size from a parameter box below, the spot is validated as peak. This is useful to get rid of smaller, unwanted spots.
- **Noise detection**
A special noise analysis filter is used to analyse the image and identify the peaks. For this method no parameters are required.

- **None**
No discrimination method is being used. Every spot which is above the given lower threshold will be treated as peak.
- **Advanced - Remove Areas inside Areas**
If this option is set areas which are located inside other areas will be declined.
- **Maximum Areas**
This parameter defines the maximum amount of peaks. If this amount of peaks is found the peak search stops despite of being completed. This will protect from long waiting times in case a very noisy image is used with hundreds of possible peaks.

To start the peaksearch select .

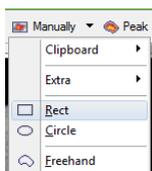
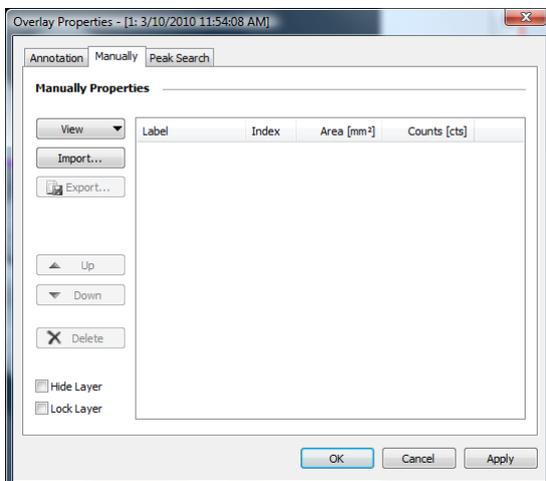
4.3.6 Define Areas Manually

The **"Manual Areas"** are available via the command **<Manually>**.

They are used for the manual definition of integration areas. Here you have the possibility to define free forms, such as circles, rectangles or polygons.

To start this function select .

A window is opened where an overview of all currently defined areas is displayed. If you have not defined any areas the list will be empty.

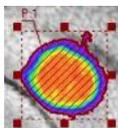


If you want to create new areas open the context menu by using the arrow on the button **<Manually>** and select the required shape (rectangle, circle or free-hand shape) from the context menu.

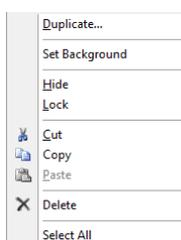
With the mouse cursor shown on the image the selected shape can be drawn by dragging into the required area of the image.

To define a freehand shape circle the desired area with several clicks on the left mouse button or hold it down completely (pencil mode). A double-click will end the procedure. The painted shape will be connected automatically and the freehand shape will be closed by a frame.

If you click on an area that has already been created, you will see "handles" around the edge of the area. The area can be resized by dragging the "handles" to the desired size.



The shape can be moved to another position. To do this you have to move the mouse to the boundary of the shape. The mouse pointer will be changed automatically to a hand cursor. If the hand cursor is displayed press the left mouse button and drag the shape to the required position. When pressing **<CTRL>** and **<Shift>** while moving the selected area this area will be duplicated. The copy mode is indicated with a little "+" next to the mouse cursor symbol. To select more than one area just press the **<CTRL>** key while selecting the area with a mouse click or drawing a box around the required areas.



A context menu for all shapes is available. To open the context menu select the shape and right-click inside the shape.

For editing the standard clipboard commands are available

To use the object for background subtraction select **<Set as background>**.

An object defined as background area is identified by a checkmark in front of the context menu entry. In this case the average intensity of the area is defined as background and subtracted from the intensities of the other areas in the report.

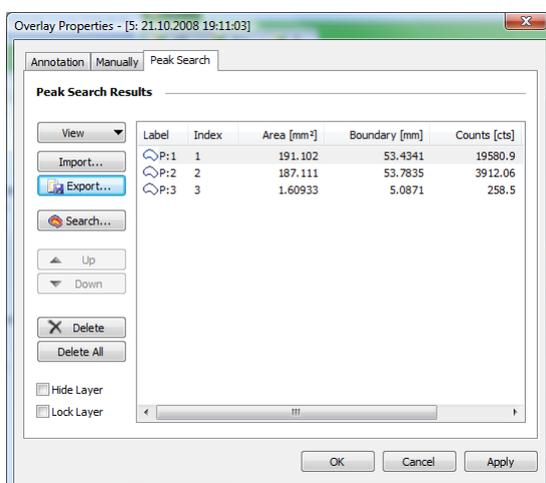
To create a copy of the current area select **<Duplicate>**.

To lock the whole layer, use the command **<Lock>**.

A locked area cannot be moved any longer unless it's being unlocked through the **"Overlay Properties" <Lock Layer>**.

The layer becomes invisible by selecting the command **<Hide>**.

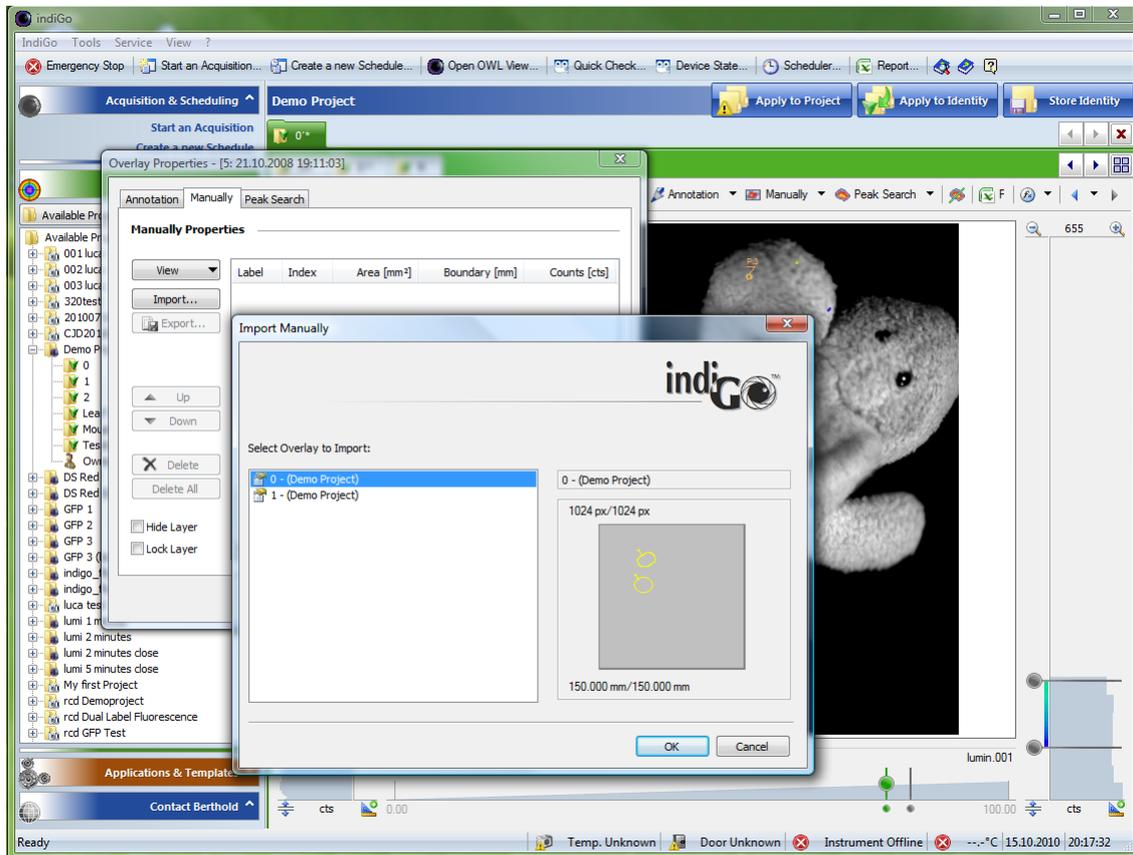
To delete all areas refer to **"Overlay Properties" <Delete All>**.



Further commands for the area management are available via the context menu of the area items list.

To export areas the **<Export>** command from the **"Overlay Properties"** can be used.

To import areas the **<Import>** command from the **"Overlay Properties"** can be used.

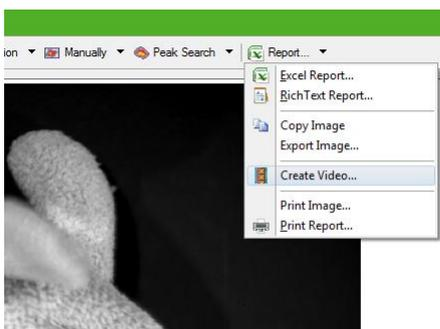


4.4 Creating a Report

4.4.1 Creating a Rich-Text-Report

indiGO offers the possibility to create and display measurement reports. The output can be adapted to the user's requirements.

Using the **<Rich-Text-Report>** command in the **<Report>** menu will start the Rich-Text-Report-Creation-Wizard.



To create a report, follow the instructions given by the wizard. In brief the following options can be set.

- **Create a Front Page**

This will create a report front page. The contents of the page, like your company name, address or logo can be setup through the indiGO options.



- **Export Images**

If this option is set, all overlay images will be included into the report. You can also select, whether the overlay areas will be included into the images. The meta data of each image will be displayed next to each image.

- **Units Area & Elongating**

Different units for the areas and the elongatings can be selected. Millimeter and centimeter can be selected for geometrical values (e.G. area or size of an area).

- **Measured Data**

Different units for the measured data can be selected.

The intensity data can be displayed in counts (raw values), the emitted photons (ph/s) or the

emitted power (nW). The photons as well as the power values are based on a calibration using a light standard. This calibration is only valid, when the sample is placed on the base plate and not higher than 50mm. In case of a different sample size height or position, the calibration is not valid and wrong results will be calculated using the default calibration. In this case a dedicated calibration must be performed using the light standard. This calibration can be applied to luminescence measurements only.

- **Rich Text Export Options**

The report can be opened to be printed or stored into a file for further usage.

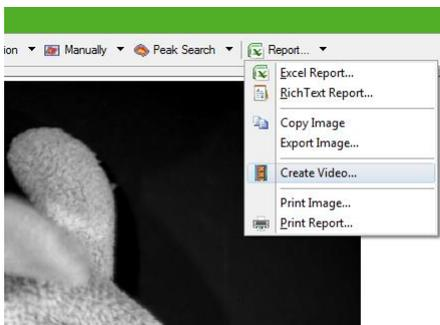


4.4.2 Excel Export

indiGO provides a very comfortable report function, the possibility to create Excel exports. In an Excel export, a complete measurement report, including all images, analysis data and in addition, diagrams containing a graphical view of the measurement results can be created.

To use this function Excel 2013 or later is recommended. Other spreadsheet programs are not supported.

Using the **<Excel Report>** command in the **<Report>** menu will start the Excel-Report-Creation-Wizard.



To create a report, follow the instructions given by the wizard. In brief the following options can be set.

- **Create a Front Page**

With this command you can create a report front page. The contents of the page, like your company name, address or logo can be setup through the indiGO options.



- **Export Area Details & Graphs**

You can select either bar graphs or line charts to be created.

- **Export Images**

If this option is set, all overlay images will be included into the report. You can also select, whether the overlay areas will be included into the images.

- **Units Area & Elongating**

Different units for the areas and the elongatings can be selected.

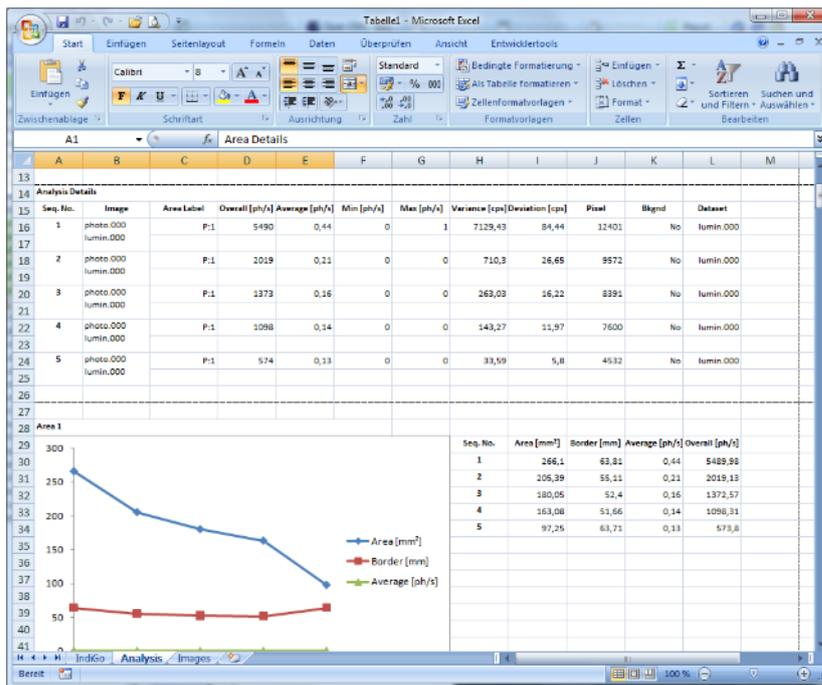
- **Measured Data**

Different units for the measured data can be selected.

- **Excel Export Options**

The report can be opened to be printed or stored into a file for further usage.





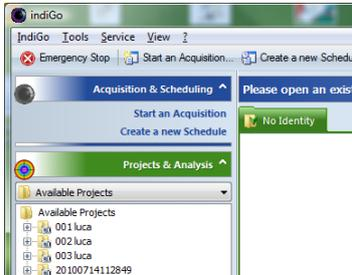
The spreadsheet will contain one to three worksheets depending on the selected options:

- **IndiGO**
The **"indiGO"** sheet contains the front page
- **Analysis**
The **"Analysis"** sheet contains all measurement results and the diagrams, depending on the settings.
- **Images**
The **"Images"** sheet includes the overlay images and all additional measurement information e.g. file name, date of acquisition, sample description and measurement parameters, to ensure reproducibility of the measurements.

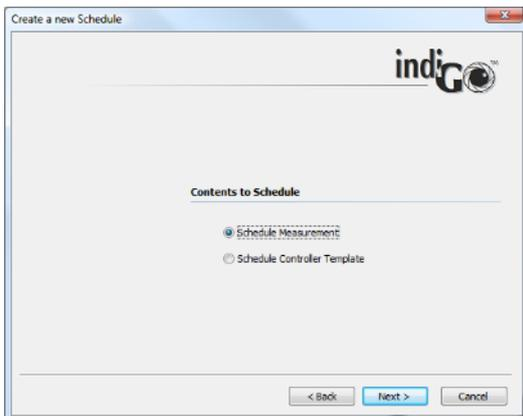
4.5 Using the Scheduler

Control sequences like sunset and sunrise or measurements for selected identities can be controlled via a scheduler. You can think of the scheduler as a simple list of time controlled events.

To start an acquisition, refer to **<Acquisition & Scheduling>** **<Create a new Schedule>**.



The system will start the **<Create a new Schedule>** wizard, which allows you to select the contents you want to schedule first.



To schedule measurements select **<Schedule Measurement>** and chose **Next >** to continue.

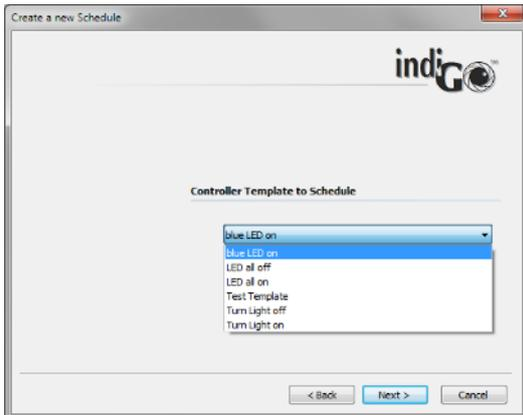


On the next page you can select the project and the identity you want to measure.

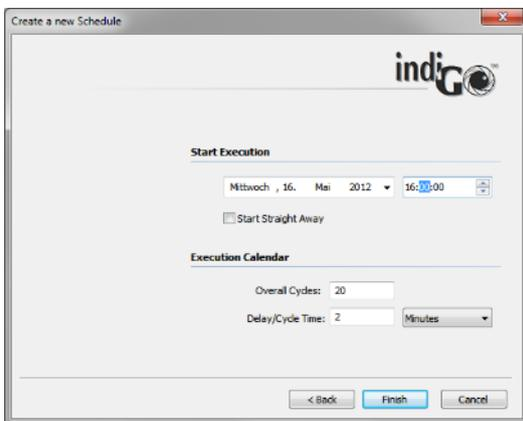
Using the **<Advanced>** or **<Tools>** commands you can validate the acquisition, pre- and post-controller settings or initiate a **<Quick Check>** to create a test image.

Chose **Next >** to confirm the selected options and continue.

When selecting **<Schedule Controller Template>** a dialog box pops up to select the desired control sequence.



Chose **Next >** to confirm the selected template and continue.

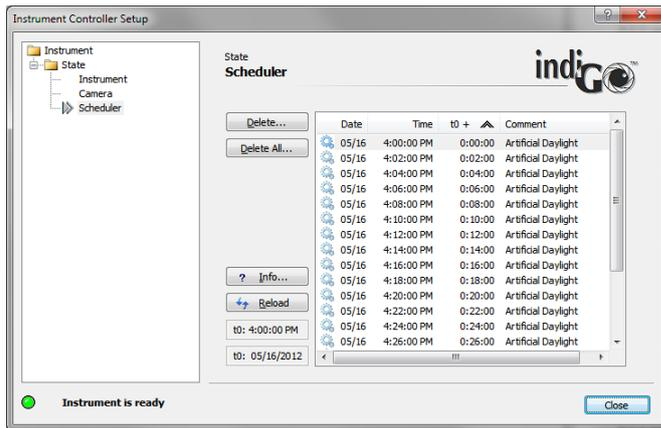


On the next page the **<Start Execution>** time can be selected. In the sample above, the first measurement will be executed on Wednesday the 16th of May at 16:00:00. When "Start Straight Away" is checked the start time cannot be entered. In this case the action starts right after clicking on **Finish**. All other repeats will start relative to this start time.

The **<Overall Measurements>** are set to 2 and the **<Delay/Cycle Time>** was set to 2, as well. Therefore the system will add two measurements into the scheduler. The first will be executed Wednesday the 16th of May at 16:00:00 and the second will be executed Wednesday the 16th of May at 16:02:00.

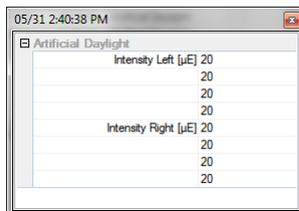
The measurements will be executed in the background by the scheduler and will be sorted automatically into the indiGO structured tree.

The contents of the scheduler can be verified or deleted via the **<View> <Scheduler>** command from the main menu of indiGO.



The scheduler shows the list of scheduled measurements and control commands with absolute starting time and relative time. The time of the first event in the list (when the scheduler was empty before) is treated as time zero.

To highlight an item just click on it. To highlight more than one item press the CTRL key while clicking on the items. A doubleclick on a single, highlighted item or clicking on will show detailed information about the event.



Highlighted items can be deleted by clicking on . To delete the whole list of items click on .

To update the list, click on . This will remove past events.

5 System Description

5.1	Use and Function	78
5.2	Design of the Acquisition Device	79
5.3	Acquisition Chamber	81
5.3.1	Light-tight Door with Door Button	81
5.3.2	Illumination of the Acquisition Chamber (Dark Box)	81
5.3.3	Connector Field	81
5.3.4	Camera and Lens	82
5.3.5	Filter Wheel	82
5.3.6	Filter Slider	83
5.3.7	Fluorescence Illumination	84
5.3.8	Bottom Plate	84
5.3.9	Top Part of Acquisition Device	85
5.3.10	Connections at the Rear Panel	85

5.1 Intended Use

NightSHADE evo LB 983 is a sensitive Imaging System for luminescence and fluorescence measurements to be used on the benchtop. It is used to acquire and evaluate images in the macro imaging area in a large intensity range.

Not for use in in-vitro diagnostic (IVD) procedures.

This instrument is not designed or intended for use with installations or equipment in hazardous environments. Servicing of the instrument must only be performed by Berthold Technologies Field Service Engineers or service staff authorized by Berthold Technologies.

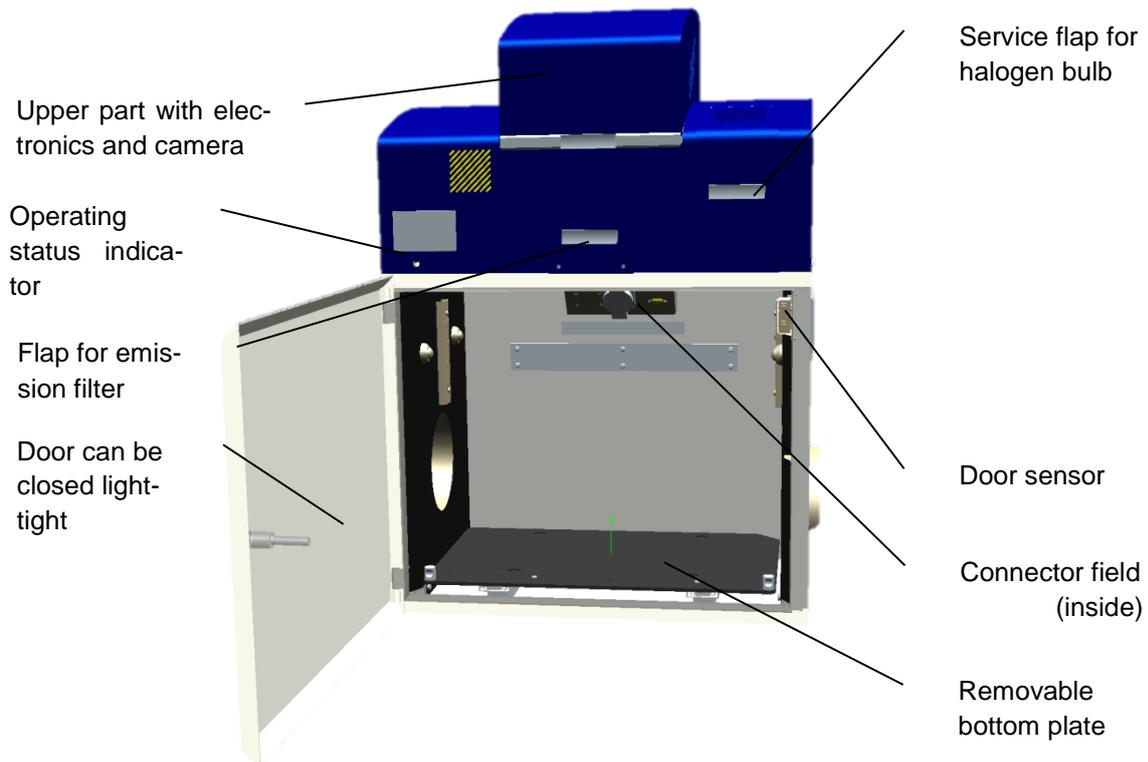
NightSHADE evo LB 985N includes the following components:

- **Acquisition device** including the transport mechanism with camera and lens and the electronics. The high-resolution CCD camera and the special wide open lens guarantee excellent sensitivity and image quality.
- The control of the acquisition system and the image processing and evaluation take place under control of the effective **software system *indiGO***. The easy-to-use and clearly structured Windows software covers the entire spectrum of image processing functions.
- The **PC system** including a standard computer (Pentium 4® or other processor) with high-resolution monitor. We recommend a resolution of at least 1024x768 pixels.
- **Image output** to any Windows printer (laser or ink-jet printer) or video printer.



5.2 Design of the Acquisition Device

The acquisition device includes the hardware and electronics required for image acquisition. It comprises the acquisition chamber (dark box) including the camera lift with camera, lens and filter wheel. The electronics and excitation light source for fluorescence measurements are accommodated in the upper part of the device. The ventilation cap is designed for drawing off the waste heat caused by the camera cooling.



Front side

- Light-tight door with handle to close the acquisition chamber
- Door for loading and unloading the filter slider for fluorescence emission
- Service flap for halogen lamp and excitation filter slider loading and unloading.
- Operating status indicator

- Inside the *acquisition chamber*
- Inductive sensor detecting the door position
- Bottom plate (removable) to position the acquisition objects
- 4 LED's illuminating the acquisition chamber
- Connector field
- Liquid cooling connectors (for LED panel)
- LED panel control connector
- Outlet of the fluorescence light for connection of the illumination light guide with optical switch
- 4 fibre optic illumination light guides for fluorescence illumination (with -40 version only)

Rear panel/Top part

- Connection ports
- On/off switch
- Mains fuse
- Mains connector
- USB B port (in)
- USB A port (out – for side view camera)
- Camera power outlet (for side view camera)
- Liquid cooling in/outlet (for LED panels)
- Ventilation slits

At the side

- Carrying handles
- Camera bayonet for side view (left)
- Flange port for external feed through

Bottom

The NightSHADE evo stands on four feet; the height of these feet can be adjusted to balance uneven setup locations.

5.3 Acquisition Chamber

5.3.1 Light-tight Door with Door Button

To close the acquisition chamber light-tight, the door frame is provided with an elastic sealing and a handle which, when turned by 180°, will pull the door into the frame and squeeze it against the sealing. An inductive sensor indicates if the door is closed correctly.

Always make sure that the door is properly closed. If the door is not closed and you try to take an image, you will be alerted by the program.

Images can be taken even while the door is open. The camera is insensitive to daylight, i.e. the camera will not be destroyed by overexposure.

5.3.2 Illumination of the Acquisition Chamber (Dark Box)

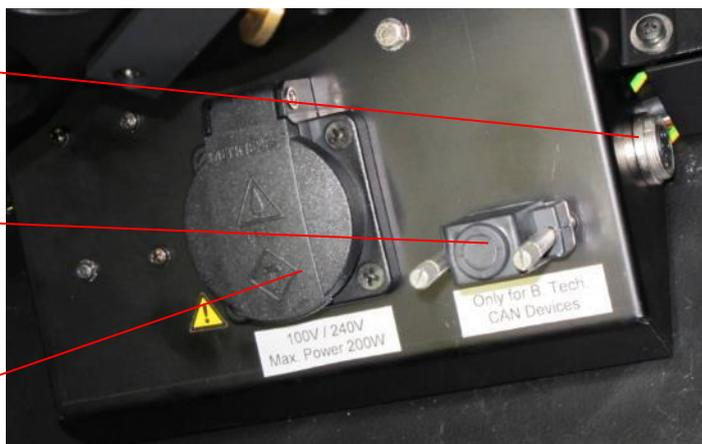
Four LED's installed on the walls of the dark box ensure constant illumination of the acquisition chamber. However, the illumination depends on the camera setting and with very low camera position it can be affected adversely. If necessary, an additional light source has to be installed (power is supplied via the outlet strip in the acquisition chamber).

The illumination of the acquisition chamber is automatically controlled by the software.

5.3.3 Connector Field

The connector field is located in the right rear corner of the acquisition chamber; it includes the following ports:

- Round sockets:* Control connector for LED panels (left and right)
- CAN socket:* System connectors to control instrument options (e.g. turn table)
- Mains socket:* Mains socket enabled via software



The mains socket is controlled through the software (see chapter 4.2.4) To prevent exposure to dangerous UV light when using e.g. a UV transilluminator the socket will switch off while the door is open.

Max. power outlet is 200VA (AC 100 – 240V +/- 10% according to the supply voltage).

The CAN socket controls external system options such as the turn table (see chapter 10.1.2) or the temperature controlled measurement table (see chapter 10.1.2).



When no external equipment is connected the dummy plug (terminator) must be connected to this socket to ensure a proper termination of the internal bus.

The two round connectors at both sides of the connection box are the power supply and control outlet for the optional LED panels (see chapter 10.1.1).



Connect the LED panels only while the instrument is powered off to prevent electrical damage of the LED panel.

5.3.4 Camera and Lens

NightSHADE evo LB 985N uses a slow scan CCD camera, hermetically vacuum encapsulated and cooled down by a Peltier element to -70°C (the operating temperature is reached after about 10 minutes).

To ensure the proper operation of the camera cooling the ambient temperature should not exceed 28°C .

The camera is connected with the electronics via two cables (voltage supply and USB communication) on its top side.

Focusing of the sample is done by turning the manual focus wheel of the lens, located at the front left side. The lenses are changeable using a quick change bajonet system. To mount the lens both red dots of lens and carrier plate must match.



Maximum sample size for standard resolution images:

Ikon (with 25mm lens) : 266 x 266mm at baseplate level up to 127 x 127mm (sample must be raised up).

With Macro lens: 130 x 130mm at baseplate level up to 60 x 60mm (sample must be raised up).

Discharge electrostatics before taking out the camera

Clean the camera and lens only with special cleaning products from your photo shop (see page 115).

5.3.5 Filter Wheel

The filter wheel is installed below the camera. The filters are located on specific filter holders. These holders can be slid into the filter wheel.



Please observe the correct installation direction (arrow on the filter wheel) when loading the filter. Please refer to the filter manufacturer's description. The filters should be secured from above against falling out by the snap rings which are included with the device.

In the filter manager you can specify filter properties. Make sure the filter numbers printed on the filter holder matches the filter position indicated on the filter wheel.

Filter position 5 should remain empty for the acquisition of photographic images.



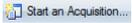
Do not reach into the filter wheel while the wheel is moving → Risk of injury.

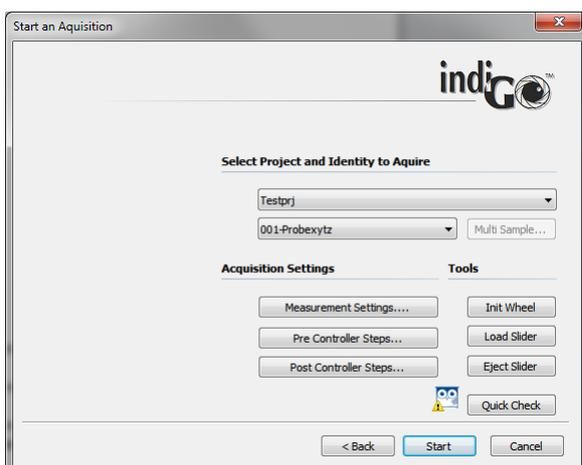
Filters can be exchanged by manually moving the wheel while the instrument is turned off.

5.3.6 Filter Slider



The filter slider can accommodate five excitation filters with a diameter of 25 mm (or 1 inch). It is located under the flap on the right top side of the instrument. A light blocking knob keeps out light to ensure best performance. To get access to the slider just remove the knob.

The filter slider can be ejected via the indiGO software. The eject command can be found under  Start an Acquisition...



To eject the slider select  Eject Slider .

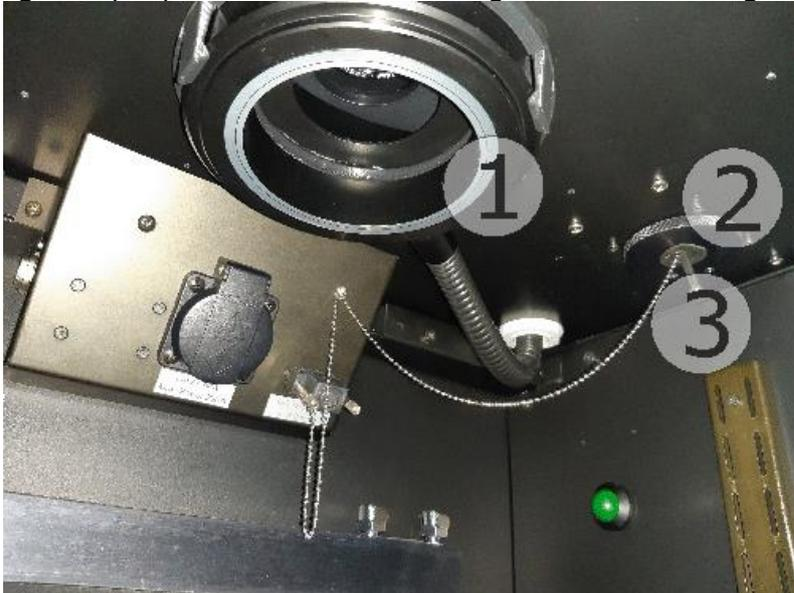
When loading the filter please observe the correct installation direction (arrow on the filter wheel).

To reload the slider select  Load Slider . After that the slider will be initialized by the instrument.

The Berthold Application Support is available to recommend filters when you are selecting the optimum filter combination for your fluorescence application.

5.3.7 Fluorescence Illumination

The built-in halogen light source together with the filter slider create the excitation light for fluorescence measurements. The light is guided into the dark box to the ring light (1). For alternative illumination devices using standard 15mm fibre optic connectors an auxiliary light output port (2) is located at the right side at the ceiling of the dark box.

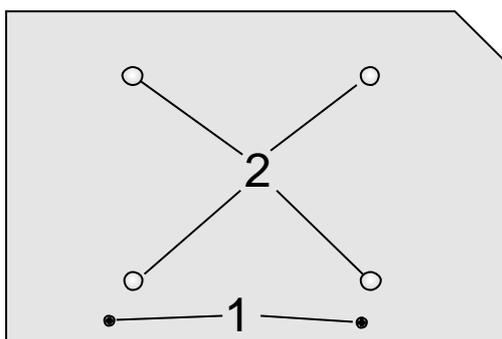


When the ring (2) is in the lower position, the light is guided to the built-in ring light. When pulling the ring to the upper position an optical switch guides the light to the auxiliary light output port. The light guide is fixed in the socket using internal spring loaded balls. To remove it pull the light guide slightly.

A blind plug (3) is placed inside the light output to prevent outside light entering the dark box. Please always keep this rubber plug in place if the auxiliary port is not required.

5.3.8 Bottom Plate

The bottom plate helps you to position the acquisition objects. It includes two bolt heads (1) on the front panel to help you position and fix the samples or sample masks accurately. In addition, there are four wells (2) for accurate positioning of transilluminators.



The bottom plate can be removed for cleaning, if liquid was spilled.

In addition, the bottom plate can be pulled out from the front to facilitate the loading of samples. When you push the bottom plate back in again, make sure that the plate snaps into place with an audible click.

5.3.9 Top Part of Acquisition Device

The upper part of the instrument contains the camera, the fluorescence light source (option) and the complete electronics.

To open lift up the lid.

Inside 2 connectors are visible. One is for the camera power supply, the other is the USB dataport to control the camera. Remove the cables before removing the camera.

The camera can be taken out by turning it counter clockwise until the two red dots are matching each other. Now the camera is unlocked and can be removed by lifting up the camera.



Please note:

The ventilation slits of the cooler must always be open!

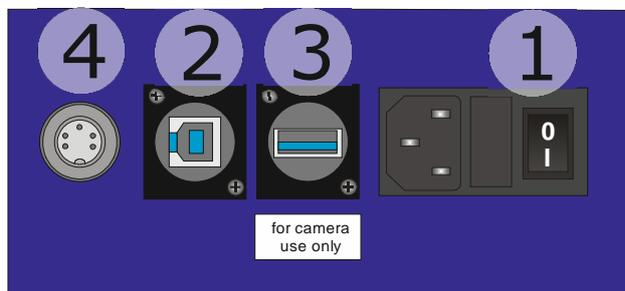
5.3.10 Connections at the Rear Panel

The power supply and device communication ports are located on the rear panel of the NightSHADE evo.

Plug the power cord into the mains inlet (1). The main power switch is located next to the mains inlet.

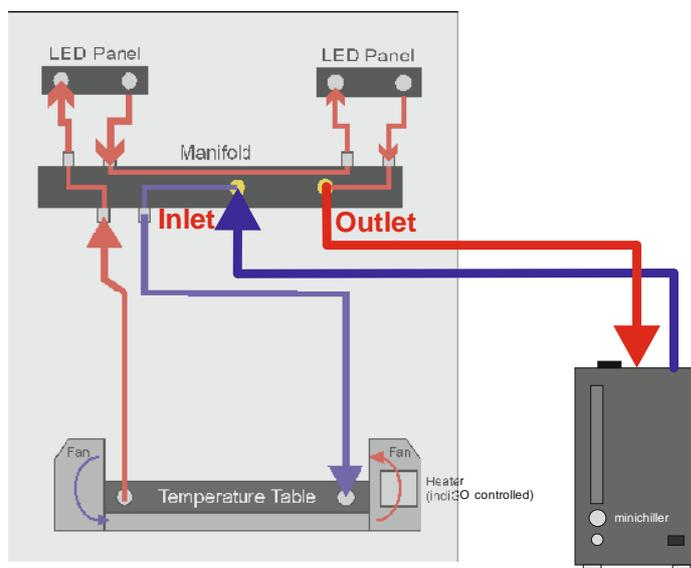
Connect the NightSHADE evo to the PC via the USB B-socket (2) on the rear panel using a USB cable (A to B).

The 7-pin Tuchel socket (4) is the power supply for the side mounted camera, the USB A socket (3) is used for the control of the side view camera.



The cooling liquid connectors are located on the lower side of the rear panel. These quick tube connectors connect the chiller to cool down the LED panels the inlet connector is located in the center of the rear panel.

Darkroom (rear view)



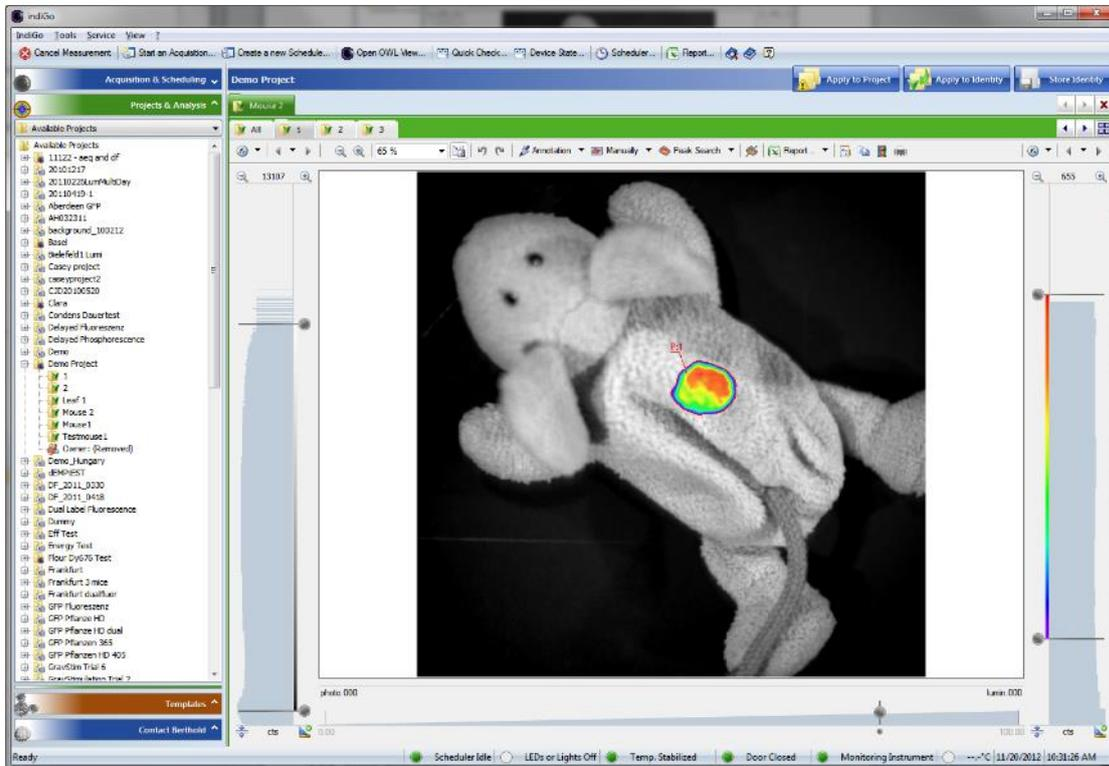
6 indiGO Software Documentation

6.1	Graphical User Interface	88
6.2	Application Toolbar	88
6.3	Identity Context	89
6.4	Identities and Measurements	91
6.5	Contrast Slider	91
6.5.1	Photo Properties Quick Menu	92
6.5.2	Photo Properties Dialog	92
6.6	Data Slider	94
6.6.1	Data Properties Quick Menu	94
6.6.2	Data Properties Dialog	94
6.7	Toolbar	97
6.8	IndiGO Menu	102
6.8.1	New	102
6.8.2	Start a new Acquisition	102
6.8.3	Create a new Schedule	102
6.8.4	Report	102
6.8.5	Print Image	102
6.8.6	Print Report	103
6.8.7	Logoff	103
6.8.8	Exit	103
6.9	Tools Menu	104
6.9.1	Backup Projects	104
6.9.2	External Tools	104
6.9.3	Import indiGO 1.x Project	104
6.9.4	Import Winlight Data	105
6.10	Service Menu	106
6.10.1	Filter Manager	106
6.10.2	User Manager	106
6.10.3	Controller Service	106
6.10.4	Device State	107
6.10.5	Transport	107
6.10.6	Quick Check	108
6.10.7	Controller Version	108
6.10.8	Messages	108
6.10.9	Options	109

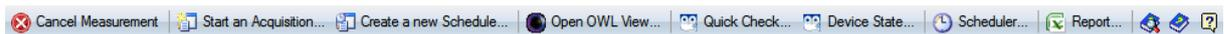
6.1 Graphical User Interface

The indiGO graphical user interface is clearly structured and offers a variety of setup options.

Measurements are automatically displayed as overlay.

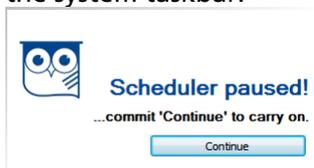


6.2 Application Toolbar



On the toolbar you find the following commands:

- **Cancel Measurement**
Will cancel the currently running measurement. The following message will show up in the system taskbar:



To continue the operations in the scheduler click on . Otherwise all actions are paused.

- **Start an Acquisition**
Start a new measurement for the current selected or any other project.
- **Create a new Schedule**
Create a new schedule for the current selected or any other project.
- **Open OWL View**
Start the indiGO image viewer.

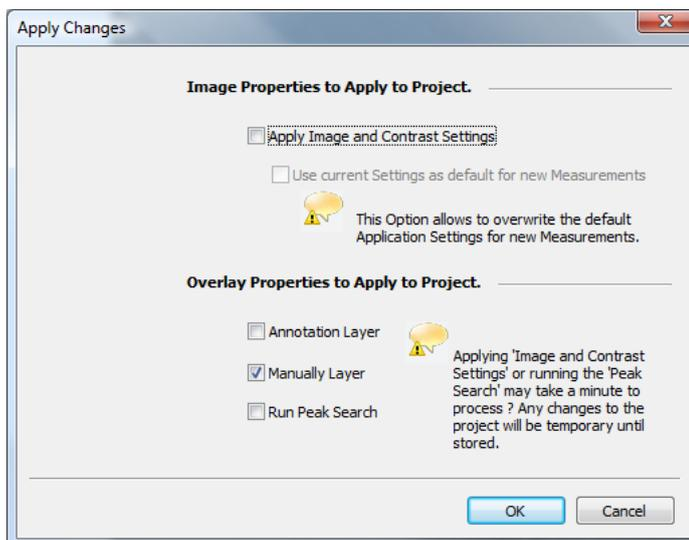
- **Quick Check**
Start the <Quick Check> utility to create a photo or run a measurement for test circumstances.
- **Device State**
Open a window to display the current state of the NightSHADE evo.
- **Scheduler**
Open the editor for the scheduler
- **Report**
Create a report for the current active project
- **Help Search**
Search the online help
- **Help**
Start the online help
- **About**
Display program version and copyright information

6.3 Identity Context



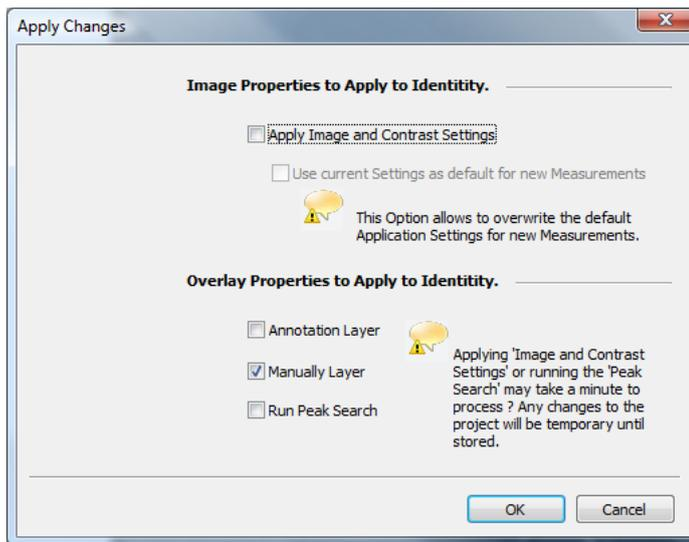
On the left of the context bar the actual selected identity name is displayed. On the right are three quick commands available.

- **Apply to Project**
With this command the current image and contrast settings, as well as the overlay properties of the annotation and the manual layers can be applied to the current project. You also have the possibility to start the peak search for all measurements of the project. .
To perform this operation, all images assigned to the selected project will be opened. According to the amount of images the operation can take some seconds up to minutes.



- **Apply to Identity**

With this command the current image and contrast settings, as well as the overlay properties of the annotation and the manual layers can be applied to all measurements of the current identity. You also have the possibility to start the peak search for all measurements of the identity.



- **Store Identity**

Save all changes made to the identity to disk.

6.4 Identities and Measurements



- **Top Row**
Currently opened sample identities
- **Center Row**
List of all measurements belonging to the current selected identity
- **Bottom Row**
Measurement or data toolbar

6.5 Contrast Slider



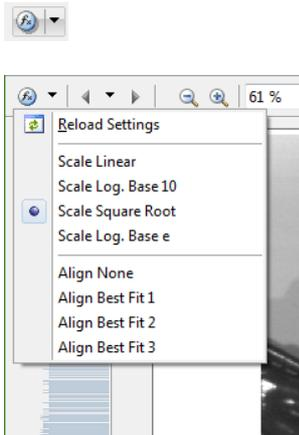
With the contrast slider on the left side of the image the brightness of the photographic display is defined. The top handle sets the upper brightness value, whereas the bottom handle is for the definition of the lower brightness limit.

The brightness can be scaled using several scaling algorithms and can be aligned automatically. These operations can be reached via the **<Photo Properties>**.

The **<Photo Properties>** can be reached through a dialog or the **<Photo Properties Quick Menu>**.

6.5.1 Photo Properties Quick Menu

To access the drop down menu simply click on the arrow next to the icon shown below.



- **Reload Settings**
Reload the last stored settings for the scaling and the alignment.
- **Scale**
Scale the contrast of the photo using the selected algorithm.
- **Align**
Align the contrast sliders using the selected algorithm.

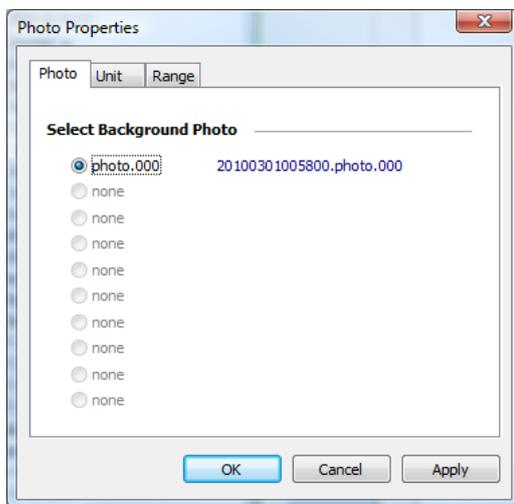
6.5.2 Photo Properties Dialog

To access the photo properties dialog simply click the icon shown below



- **Photo**
The Photo page is used to select the active photo.

If a measurement contains more than one photo, the photo displayed with the measurement can be selected.

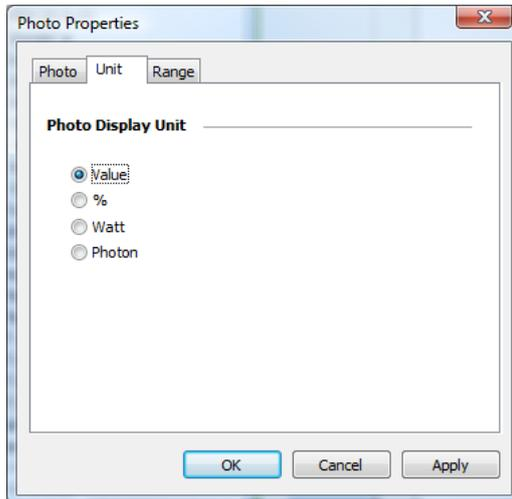


Reference:

The real name of the photo is displayed as a link. In our example: 20100301005800.photo.000, where *201003* is the date the photo was taken. In our example: 03. March 2010. 005800 is the recording time. In our example: 12:58:00 AM. By following the link, the indiGO image viewer will be opened and the photo will be displayed.

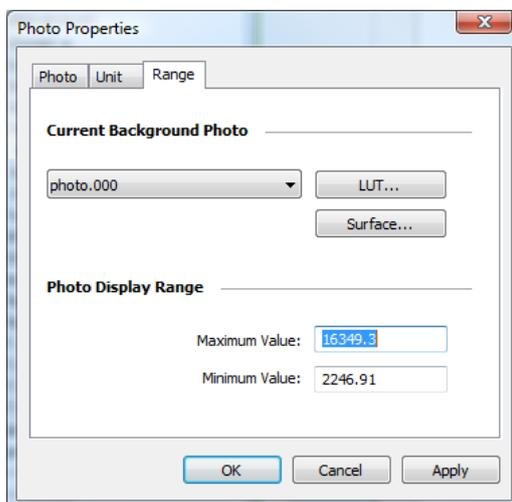
- **Unit**

The Unit page can be used to setup the unit of the contrast slider

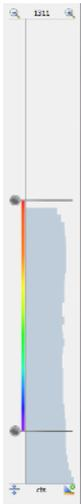


- **Range**

The Range page can be used to setup the contrast slider. For further analysis a LUT and a surface is supported.



6.6 Data Slider



On the right margin of the screen you see the intensity scale of the measurement, together with the associated slider. With these two sliders you can set the scaling of the pseudocolor signal image and the lower threshold of the color view. The top slider defines the intensity value.

The top slider defines which intensity value has been assigned to the highest valued color. Values above this threshold are displayed in white.

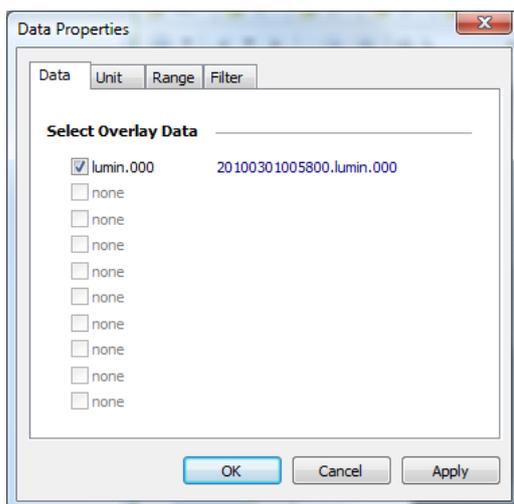
The bottom slider defines which intensity in the signal image is displayed with the lowest valued color. All values below this threshold are not displayed.

6.6.1 Data Properties Quick Menu

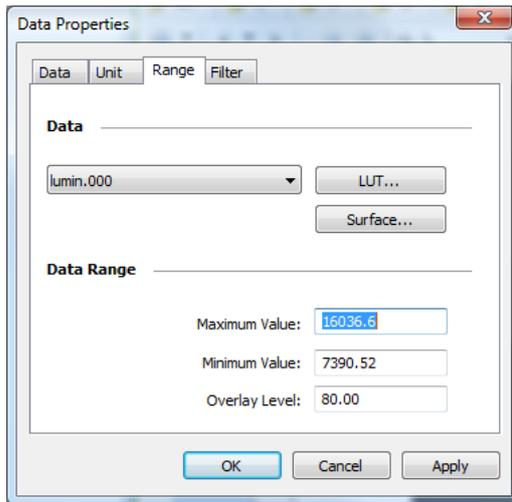
The <Data Properties Quick Menu> supports the the same functions as the <Photo Properties Quick Menu>.

6.6.2 Data Properties Dialog

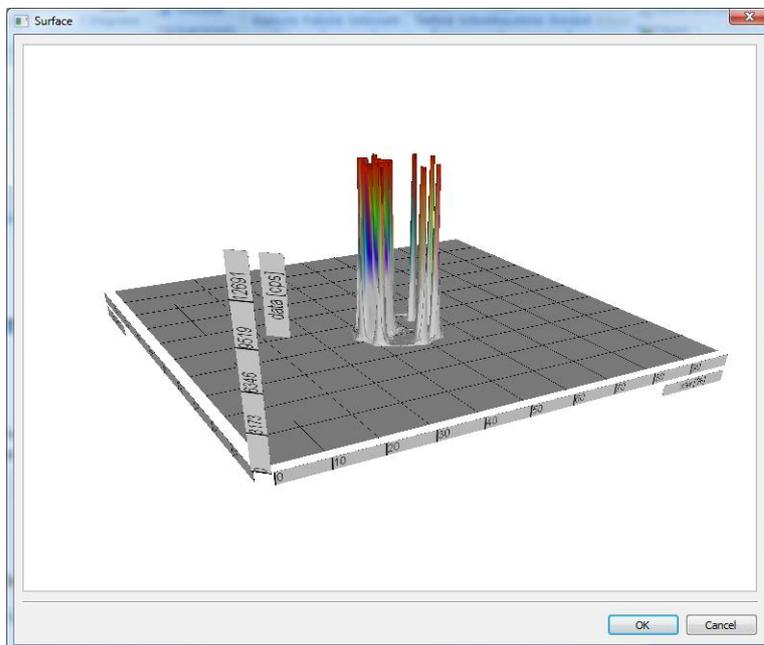
- Data**
 The Data page is used to select the measurements displayed. Each data image recorded can be tagged to be relevant and displayed along with the photo.



- Range**
 The Range page can be used to setup the data slider.

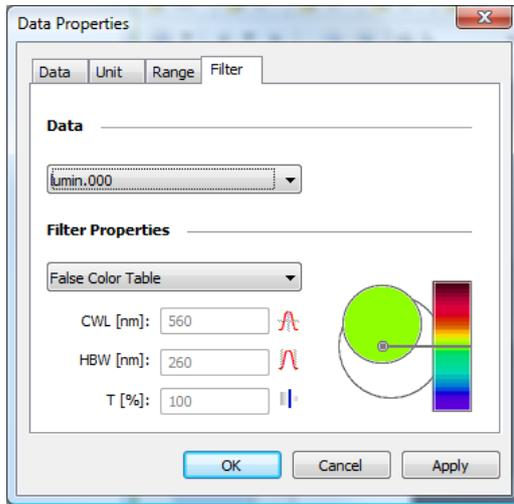


For further analysis a LUT and a surface is supported. In addition the overlay intensity can be set through **<Overlay level>**.



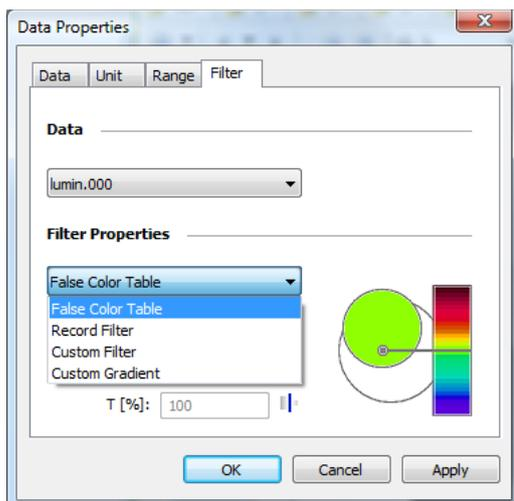
- **Filter**

The Filter page can be used to change or setup the filter properties used to render the overlay. This might be useful when more than one signal layer must be displayed in order to assign different properties to each signal layer.



The following options are available:

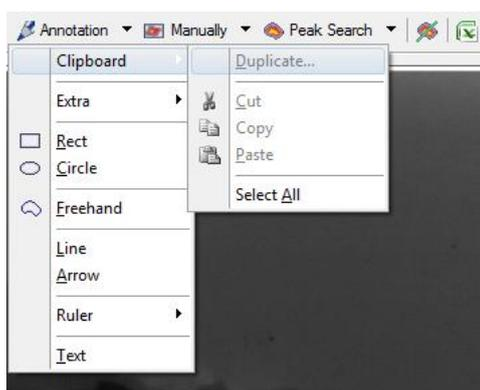
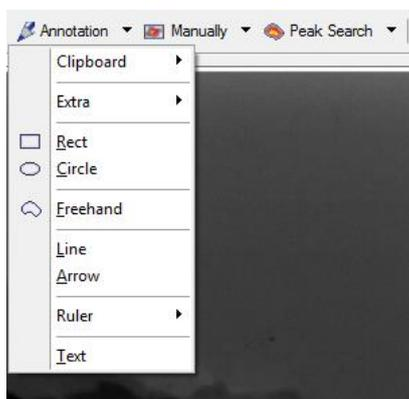
- **False Color Table**
Activates the system default false color table
- **Record Filter**
Activates the filter properties used for recording
- **Custom Filter**
Enables the user to setup a custom filter, using a center wave length and a bandwidth.
- **Custom Gradient**
Enables the user to define a gradient using a center wave length.

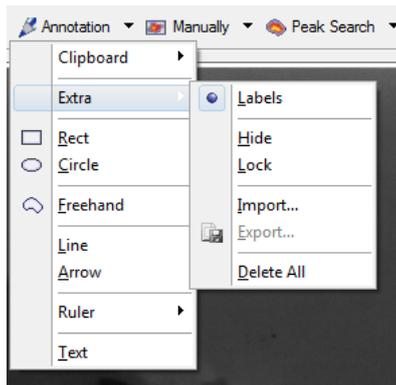


6.7 Toolbar

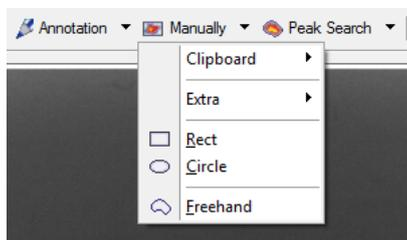


- **Decrease the zoom**
Used to decrement the zoom
- **Increase the zoom**
Used to increment the zoom
- **Current zoom**
Used to setup the zoom manually, displays current zoom factor
- **Align**
Align image into the view; used to calculate a best fit zoom factor
- **Undo (area editing)**
- **Redo (area editing)**
Used for undo and redo while area editing is active
- **Annotation (Menu)**
Create annotation areas

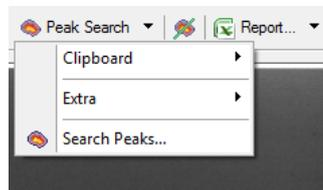




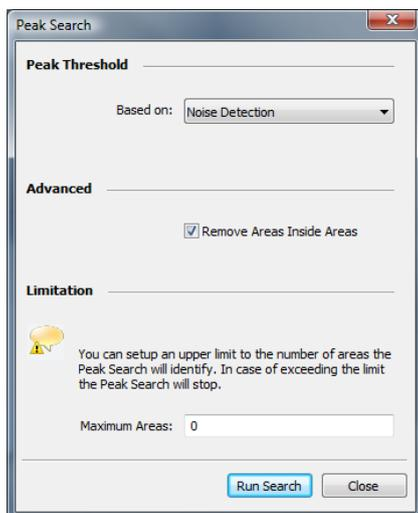
- **Manually (Menu)**
Create manual areas



- **Peak Search (Menu)**
Starts the peak search

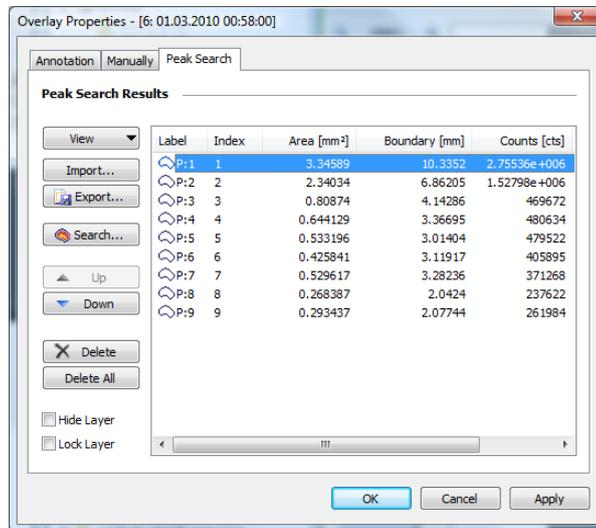


- **Peak Search (Command)**



- **Overlay Properties (Dialog)**

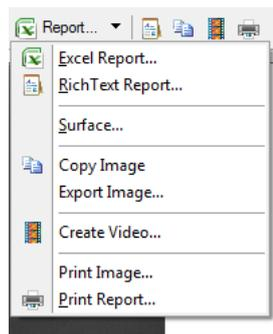
The <Overlay Properties> dialog can be opened via the <Annotation>, <Manually> or <Peak Search> commands



- **Line Plot**

Starts the Line Plot Tool

- **Report (Menu)**



- **Report (Command)**

Starts the wizard to create an Excel report. Same function as <Excel Report>.

- **Export Rich Text Report**

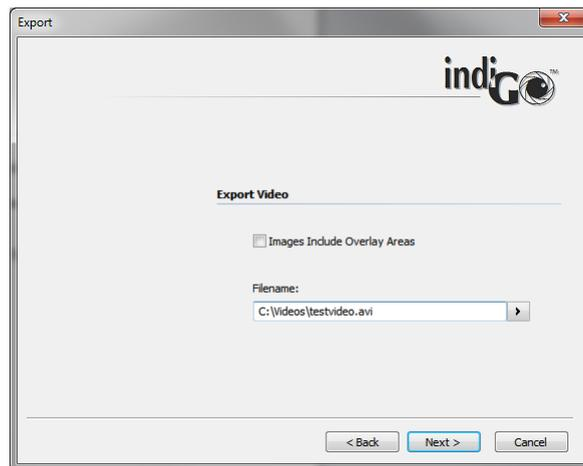
Starts the wizard to export a Rich Text report. Same function as <Rich Text Report>.

- **Copy current view to clipboard**

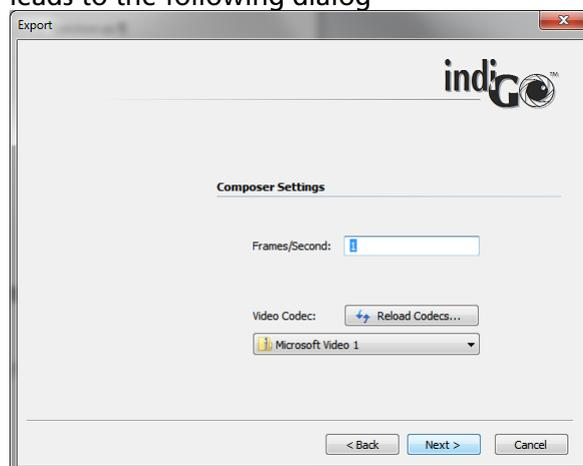
The current presentation of the measurement will be copied to the clipboard. Same function as <Copy Image>.

- **Render Video**

Starts the wizard to create a video of the measurements. Same function as <Create Video>.



The output name and destination of the video must be entered. Clicking on [Next >](#) leads to the following dialog



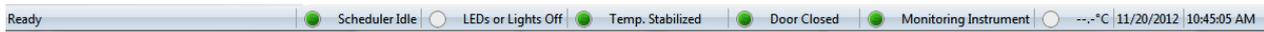
The frames per seconds for the video can be entered here. Depending on the amount of sequence images this will define the length of the video. A codec must be selected. When running this function the first time, all available video codecs must be identified by clicking on **<Reload Codecs..>** The "Microsoft Video 1" codec is recommended since it is available on almost any PC.

- **Print Report**

Starts the wizard to print a report. Same function as **<Print Report >**.

In the area below the image is a separate slider, which can be used to define the transparency of the data images.

6.8 Status Bar



In the status bar the current status of the hardware and software is shown.

Ready shows the current status of the system. When a measurement is running a progress bar is shown.

Scheduler Idle displays the status of the Scheduler. If the scheduler is busy (measurement / control jobs pending) the icon is yellow.

LEDs or Lights Off indicates if the optional LED panels inside the instrument are turned on or off.

Temp. Stabilized shows if the camera is initialized and the cooling status.

Door Closed displays the status of the dark room door. When the door is open an additional warning in the Windows status bar pops up.

Monitoring Instrument indicates if the instrument is initialized and properly connected to the PC.

---°C shows the current temperature of the heated sample tray (if installed)

11/20/2012 | 10:45:05 AM displays the current date and time.

6.9 IndiGO Menu

6.9.1 New

Project

To create a new project, start the wizard. For further details please refer to section 4.2.5.

Controller Template

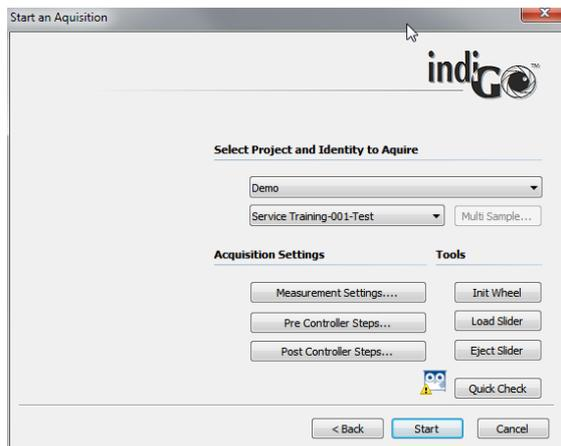
To create a new controller template, start the wizard. For further details please refer to section 4.2.4.

Measuring Template

To create a new measuring template, start the wizard. For further details please refer to section 4.2.3.

6.9.2 Start a new Acquisition

This command will start the acquisition wizard for on demand measurements.



Several quick check commands are available to verify the settings made to the application. For further details please refer to chapter 4.2.6.

6.9.3 Create a new Schedule

The system will start the **<Create a new Schedule>** wizard, which allows you to select the contents you want to schedule first.

For further details please refer to chapter 4.2.7

6.9.4 Report

indiGO offers the possibility to create and display measurement reports. The output can be adapted to the user's requirements. For further details please refer to chapter 4.4

6.9.5 Print Image

indiGO offers the possibility to create and display measurement reports. The output can be adapted to the user's requirements. For further details please refer to chapter 4.4

6.9.6 Print Report

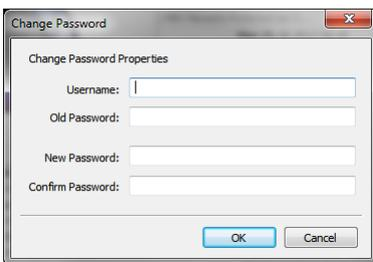
indiGO offers the possibility to create and display measurement reports. The output can be adapted to the user's requirements. For further details please refer to chapter 4.4

6.9.7 Logoff

This command will change the current user account.



When clicking on "Change password" a new password can be assigned. The old and new password must be entered as well as the user name.



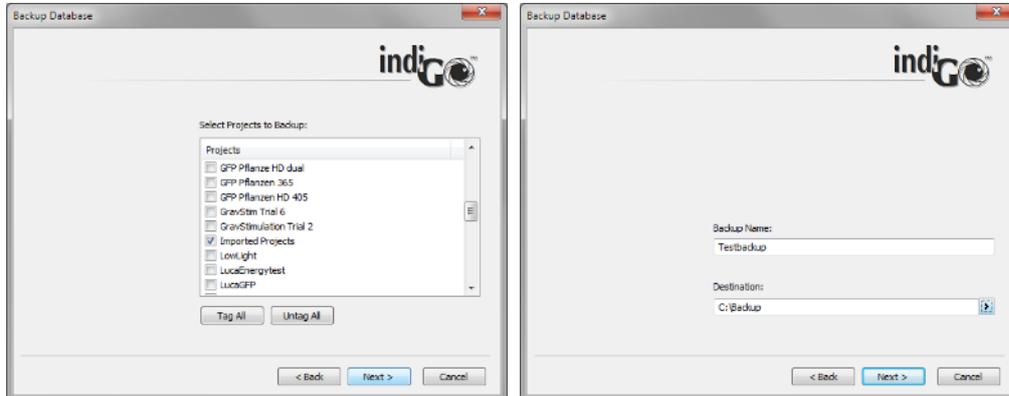
6.9.8 Exit

Quit indiGO to come back later.

6.10 Tools Menu

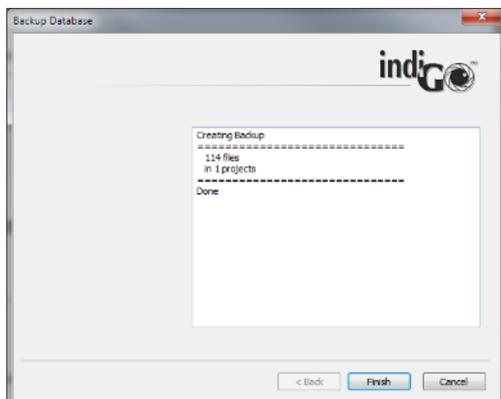
6.10.1 Backup Projects

indiGO can create backups of the database or selected projects. Backups are created using the **backup smart wizard**.



The wizard will guide you through the steps necessary to successfully create a backup. When selecting the Backup first define which projects have to be backed up. Clicking on [Next >](#) leads to the dialog box to select the backup target directory. This should be a remote network drive or removable storage to assure a proper backup. A Backup name can be assigned. This is the name of the file created by the backup (extension zip.dba). Depending on the amount of data and the computer- / transfer performance the backup may take a while.

After a successful backup a summary page will pop up.

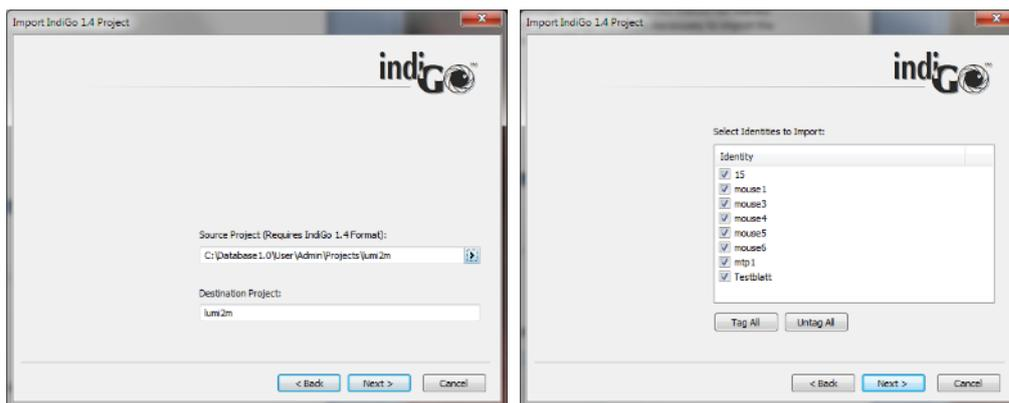


6.10.2 External Tools

indiGO supports some external tools which can be accessed through this menu.

6.10.3 Import indiGO 1.x Project

Projects and measurement data created with indiGO 1.4 can be imported into indiGO for reanalysis via this command. A smart wizard will guide you through the steps necessary to import the data. Please follow the instructions given by the wizard.



After launching the wizard just select the project file (.proj) of the indiGO 1.x project to import.

The "Destination Project" will define the Project name where the imported files will be copied to. An import project name cannot be used for data acquisitions.

Clicking on **Next >** will lead to a selection dialog box. Here you can select the Sample IDs to import. Uncheck the boxes next to the SampleID which should not be imported. The data import will start when clicking on **Next >**. After the successful import a summary page will pop up.

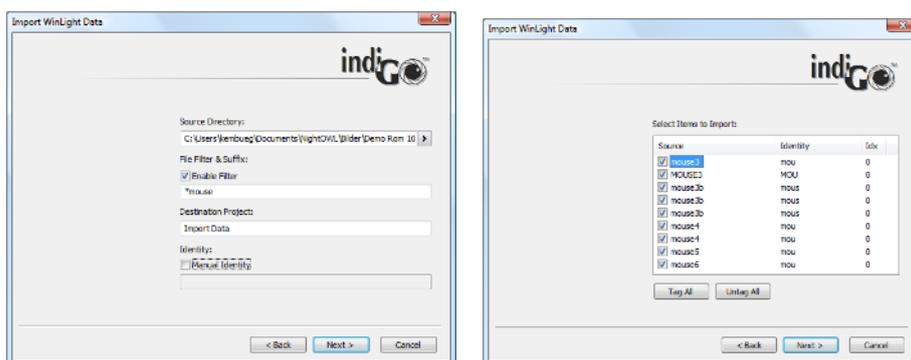
6.10.4 Import Winlight Data

Projects and measurement data created with WinLight can be imported into indiGO for re-analysis via this command. A smart wizard will guide you through the steps necessary to import the data. Please follow the instructions given by the wizard.

The file structure of the files to be imported must fit the file structure of images created by the "ReporterGen" application in WinLight.

Photographic image xxx_yyy-p.fts
Signal Image xxx_yyy-F.fts

Only image pairs (signal image plus photo image) can be imported.

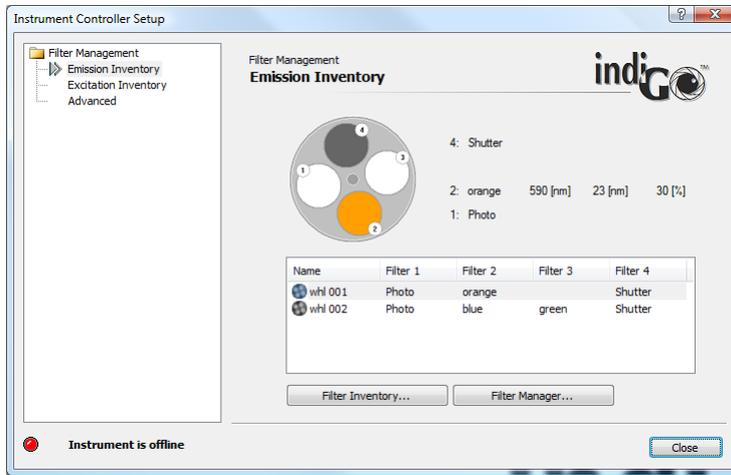


Using this structure xxx will be taken as SampleID (Identity). The 3 digit number yyy represents the sequence image number. The suffix defines the image type (photo -p; signal -f).

6.11 Service Menu

6.11.1 Filter Manager

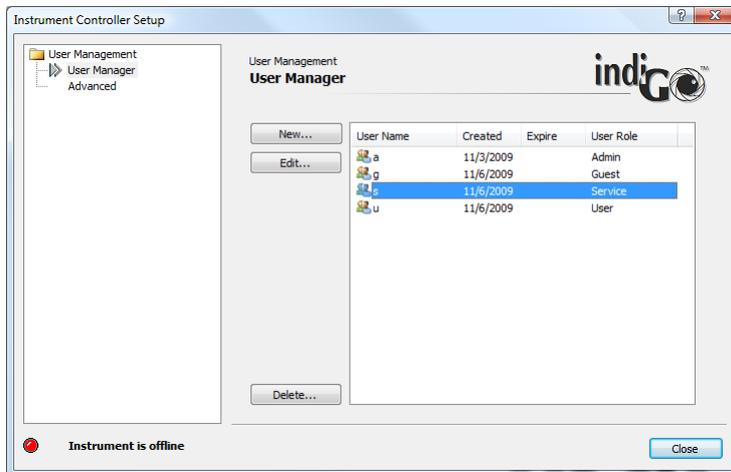
This command will start the indiGO filter manager.



For further details please refer to section 4.1.14

6.11.2 User Manager

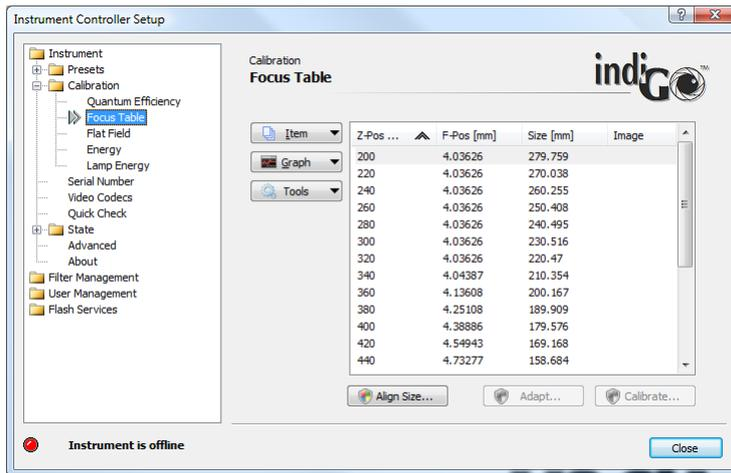
This command will start the indiGO user manager.



For further details please refer to section 4.1.12.

6.11.3 Controller Service

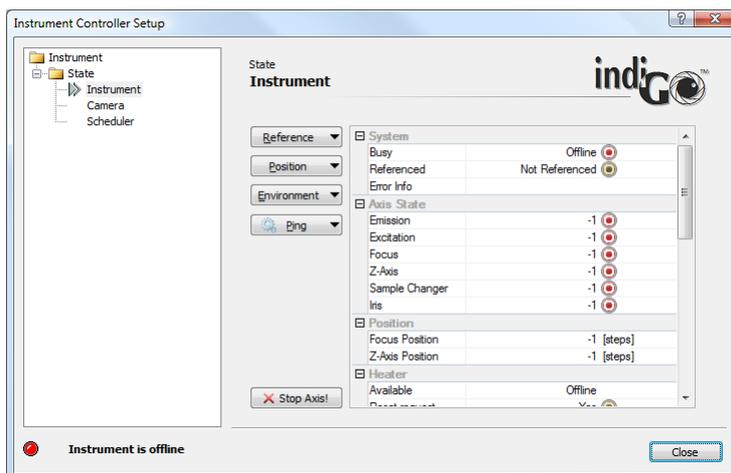
The <Controller Service> command will start the Instrument Controller Setup.



The Instrument Controller Setup is used to setup your NightSHADE evo or NightOwl device and to perform all system calibration tasks.

6.11.4 Device State

The **<Device State>** command can be used to display the current states of the instrument, the camera and the scheduler.



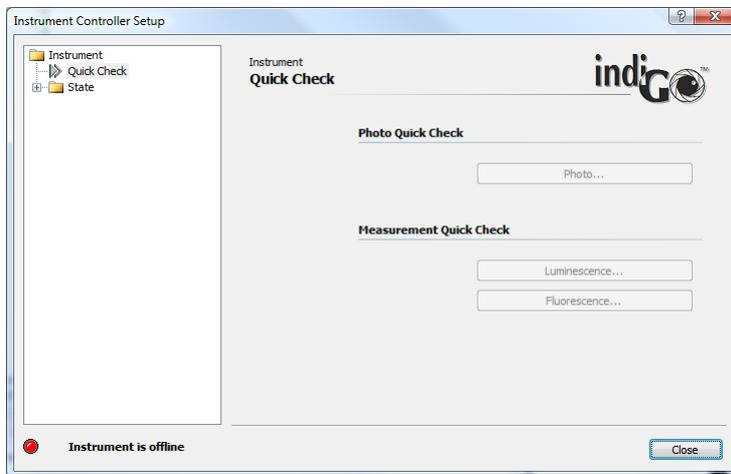
Raw device commands and test operations can be executed using the context command buttons.

6.11.5 Transport

This menu item has no functionality using the NightSHADE evo.

6.11.6 Quick Check

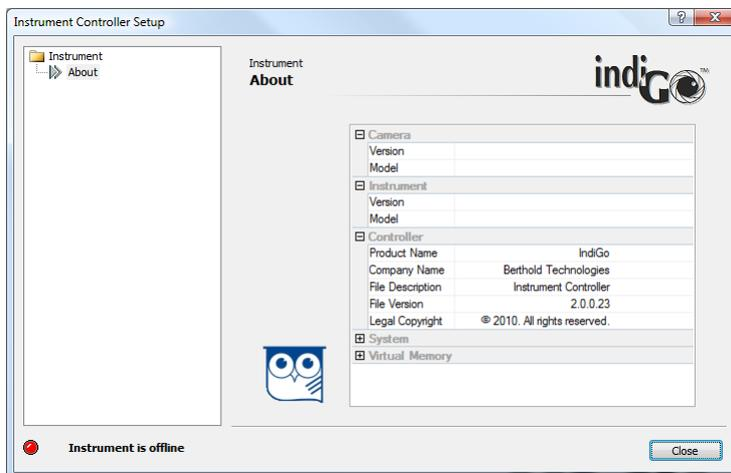
The **<Quick Check>** command can be used to get access to highlevel quickcheck commands for photo, luminescence and fluorescence.



Each highlevel quickcheck command will start a smart wizard which will guide you through the necessary steps to perform the quickcheck.

6.11.7 Controller Version

The command will display the current controller and instrument version information.

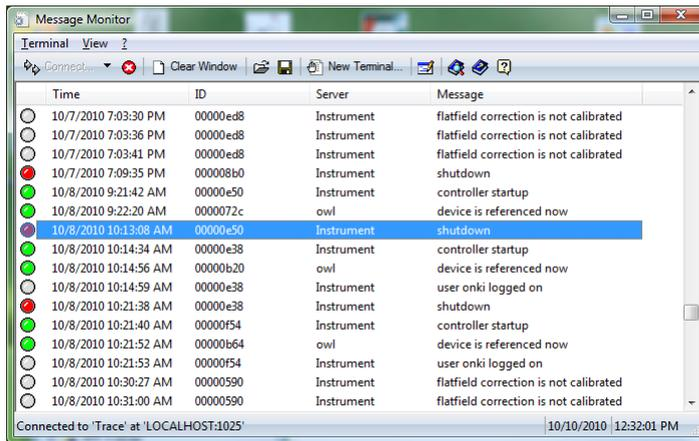


Please note:

Watch the owl while moving the mouse.

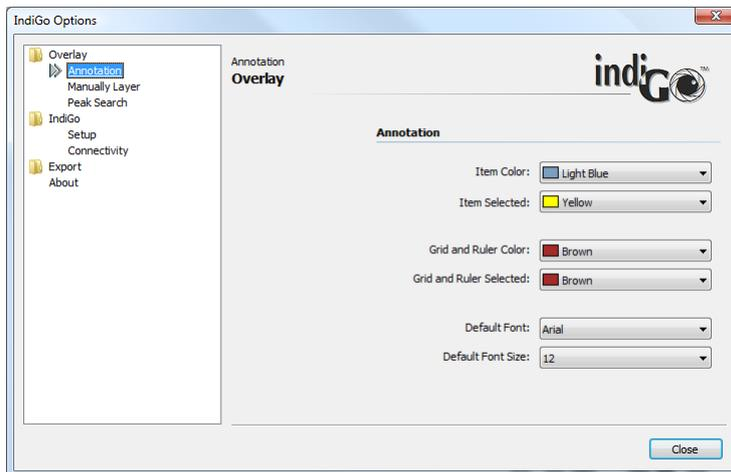
6.11.8 Messages

The **<Messages>** command will start the systems **Message Monitor** which can be used to review internal indiGO trace and error messages.



6.11.9 Options

The Options command allows you to setup further indiGO options.

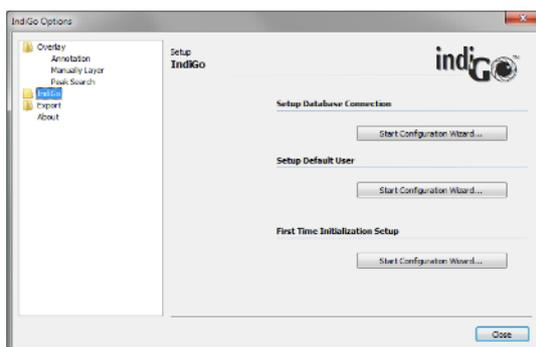


Overlay

Access to overlay colour options. Here user specific colors for the different data types (peak search areas, manual areas etc.) can be defined.

IndiGO

Access to setup configuration wizards and connectivity settings necessary for update and support mail access.



Setup Database Connection

Using this wizard the connection type can be changed. This might be required when the computer was moved away from the system or in case the database was changed from the local disk to a network drive.

Enter the type of connection first (usually direct connection to the instrument) and select the location of the database. If the database is not present at the selected location, a new, empty one will be created.



Finally the connection must be entered. In case of a local connection, the default "LOCALHOST-1024" is OK.



Setup Default User

With this wizard a default user can be assigned who logs in automatically after launching indiGO.



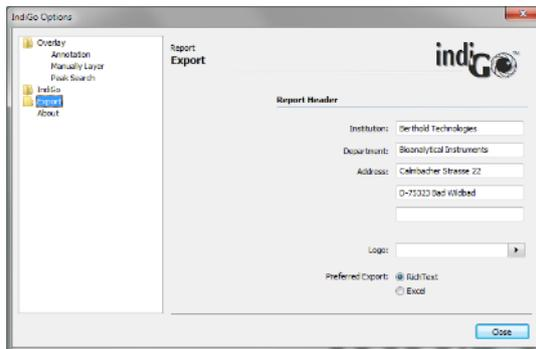
The user management can be disabled completely or a default user can be assigned.

First Time Initialisation Wizard

This button will start the initial initialization. See chapter 0 for details

Export

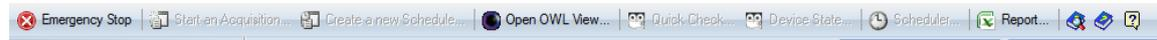
Export report settings like address, company logo and preferred format.



6.12 View Menu

6.12.1 Toolbar

This command shows or hides the toolbar.



This option is not persistent and the toolbar will be shown on the next start of indiGO.

6.12.2 Status Bar

This command shows or hides the status bar.



This option is not persistent and the statusbar will be shown on the next start of indiGO.

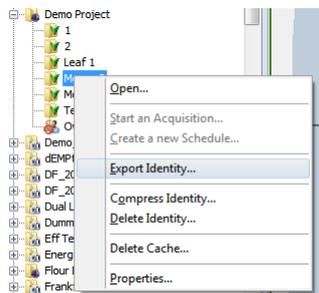
6.12.3 Scheduler

Open the editor for the scheduler. For further details please refer to section 0.

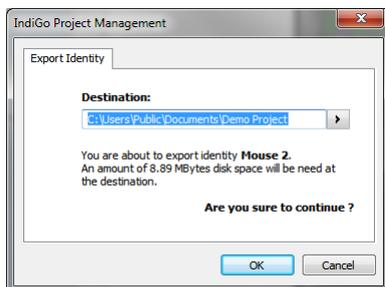
6.13 Batch Image Export

IndiGo allows the batch export of images within a sample identity. The images are exported as 16 bit gray scale TIFF files with meta data attached as XML file.

To start the export, select the required sample ID from the tree and right click on it.



Select "Export Identity" and a dialog box to select the destination folder will pop up.



Click on the arrow to select the desired folder and on **OK** to start the image export. A subfolder with the sample ID name will be generated. Depending on the amount of images, the export could take a while and requires a large amount of disk space.

The file name structure is as follows:

Sequence number (3 digits)_original filename (date coded 20120523123456).image type (photo; lumin;fluor).imagenumber (000-009).tif

7 Trouble Shooting

7.1 Hardware

7.1.1 Door status always open

If the door status is always open, the reason could be a disaligned door sensor. The door sensor is located at the upper right front side of the dark box. To adjust the door sensor refer to 8.2.3

7.1.2 Fluorescence lamp not working

The built in halogen fluorescence excitation light source has a lifetime of approx. 30h. When the lamp is no longer working replace the lamp. Refer to 8.2.4

7.1.3 No Fluorescence light output

When using the 4 light guides for fluorescence illumination it could happen, that the fluorescence light is not present at the light guides. In this case check the position of the optical switch (see chapter 5.3.7) the optical switch must be in the lower position to get light output at the 4 fibre optic light guides.

7.2 Software

7.2.1 Surface plot not visible

When running the surface plot from the Report menu a box with an error message (D3D device error) pops up instead of the surface plot. This indicates that the installed graphics card or the corresponding driver does not support Microsoft Windows DirectX 9 features.

In this case a driver update will solve the problem since many older PC do not have the latest drivers installed. In case the graphics card does not support DirectX 9 a new graphics card is required.

8 Cleaning and Maintenance

8.1	Camera and Lens	115
8.1.1	Removing the bottom plate	115
8.1.2	Cleaning the Dark Box	115
8.1.3	Cleaning the Camera	115
8.1.4	Cleaning the Lens	115
8.2	Darkbox	116
8.2.1	Replacing the Air Filter	116
8.2.2	Replacing the Fuse	116
8.2.3	Setting the Door Sensor	116
8.2.4	Replacing the halogen lamp	117

8.1 Camera and Lens

8.1.1 Removing the bottom plate

The retractable bottom plate is removable. The bottom plate is locked by a latch underneath the telescopic rail. To remove the plate open the latch on both rails. The black plastic latch is located inside the rail. After unlocking the plate it can be carefully slid out to the front.

To reinstall it carefully slide in both rails at plate to the counterpart located at the bottom of the instrument until both latches lock with an audible click. Make sure the plate is inserted properly with the knobs facing the front.

8.1.2 Cleaning the Dark Box

The dark box should only be cleaned using a mild detergent and a lint-free tissue or a vacuum cleaner. To this end, the bottom plate which is locked by a latch underneath the telescopic rail should be removed (see above).

8.1.3 Cleaning the Camera

☒ *Clean the camera only with pressurized air spray or optical paper!*

☒ *Remove the camera for cleaning.*

Dismantling and cleaning the camera

- Turn the camera to unlock the bayonet lock.
- Carefully lift the camera off from above.
- Do **not** detach the cable from the camera.
- If the camera has a shutter, start an acquisition with long exposure time via the software. This opens the camera shutter at the bottom and allows you to remove dust particles.

Caution: Do not damage the shutter by touching it when it's closed!

- Clean the camera using pressurized air spray.

Installing the camera

- Install the camera with both screws to the transport unit. The port must face the camera lift.

8.1.4 Cleaning the Lens

The lens can be cleaned when it is removed from the instrument. To do so, carefully unscrew the lens from the C-Mount thread. Hold the lens with both hands since it could be damaged when falling down on the bottom plate when the thread is completely unscrewed.

It should only be cleaned using pressurized air spray or cleaning agents available in photo shops!

To reinstall the lens hold it with both hands and carefully screw it to the C-Mount. Make sure the F-stop is set to 0,95 and locked to avoid loss of light.

8.2 Darkbox

8.2.1 Replacing the Air Filter

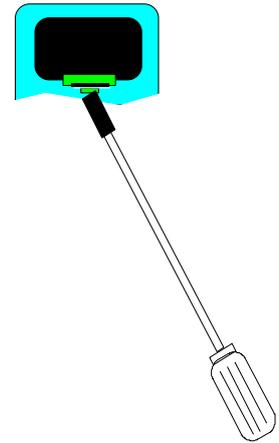
The air filter should either be cleaned or replaced, depending on the level of contamination. Remove the black ventilator cover on the rear of the instrument and clean the filter (blow it out or wash it) or replace it.

8.2.2 Replacing the Fuse

The mains fuses are located on the rear panel of the *NightSHADE evo*. The fuse holder and the mains connection socket are one unit.

Proceed as follows:

- Pull the power cord.
- Insert a screwdriver from below into the opening of the fuse inset.
- Pull out the fuse inset with the screwdriver like a drawer.
- Replace the fuse.
- Insert the fuse holder again until it clicks into place.



- + ***Use only certified fuses (original spare part) that match the value and the line voltage indicated on the device. (250V T 6,3AL)***

8.2.3 Setting the Door Sensor

In case the sensor on the door frame should be displaced so that there is no contact although the door is closed, you can easily re-adjust the sensor as follows:

- Open the locking screw at the bottom of the sensor using a 1.5mm Allen key.
- Push the sensor out inserting the Allen key at the rear side of the sensor housing, so much that it protrudes by about 1 cm. Do **not** tighten the locking screw.
- Carefully close the door and lock it so that the sensor is pushed through the door into its correct final position.
- Open the door carefully and tighten the locking screw.

8.2.4 Replacing the halogen lamp

The halogen lamp is located close to the excitation filters. It is protected by a black cover plate.



To replace the lamp proceed as follows. Make sure that the same type of lamp is used to guarantee proper light output.



Be careful when touching the lamp since it could be hot when the lamp was in use before.

- Switch off the instrument and unplug the mains connector. If the lamp was in use before allow some time to cool down the lamp.
- Remove the screw that closes the flap to open the top flap to get access to the lamp module.
- Pull out the black protection cover to get access to the lamp.
- Remove the lamp from the snap in holder.
- Unplug the lamp socket and plug in the replacement lamp.
Do not touch the reflector of the lamp to avoid damage of the lamp.
- Place the lamp back into the snap in holder. Make sure the lamp is fitted properly in the holder (round edge fit with lamp reflector).
- Insert the protection cover. Make sure the cover is properly fixed and both pins are securely fastened in the mounting holes.
- Close the top flap and fix the screw.

9 Technical Data

Camera	Peltier/air-cooled CCD camera C-Mount adapter
Resolution	1024 x 1024 Pixel x 13 μ
Quantum efficiency (max.)	>90%.
Noise	Readout: Σ 7 e ⁻ RMS
Exposure time	from milliseconds to hours
Binning of pixels	Variable in X- and Y-direction to increase the signal/noise ratio
Temperature	Ambient temperature during operation: 10° to 28° C Cooling temperature - XX 0°C Ambient temperature during storage: -20° to 70° C No condensation
Lens	standard: 25mm f = 0,95 macro: 50mm f = 0,95
Interface	USB 3 (B-type connector)
Illumination	internal illumination can be adjusted via the software optional white light fluorescence illumination (400-800 nm)
Max. sample size	266 mm x 266 mm
Min. sample size	130 x 130 mm with standard lens (65 x 65mm with macro lens)
Dimensions	Acquisition device: 600 mm x 900 mm x 400 mm (WxHxD) Camera head: 110 mm x 175 mm x 110 mm Sample chamber (inside): 400 mm x 380 mm x 450 mm
Weight	approx. 40 kg (depending on version)
Operating voltage	AC 100-240 V (+/-10%); 50-60Hz
Power rating	600VA, (max. 200VA through internal power socket)
Fuse	250V T 6,3AL
Computer	Minimum requirements: Standard PC with 3 GHz Pentium processor (MultiCore recommended) 4 GB RAM. Free USB2 port Graphics card with 1280x1024 pixels resolution 32 bit colors, DirectX 9c Support 17" LCD monitor Windows 10 (32 or 64 bit)

10 Appendix

10.1 Accessories

10.1.1 LED Panels



The LED panels are used to simulate environmental conditions like day / night changes while the sample is placed inside the dark box and no measurement is running.

The LED panels are available in 2 different versions with the following wavelength configuration:

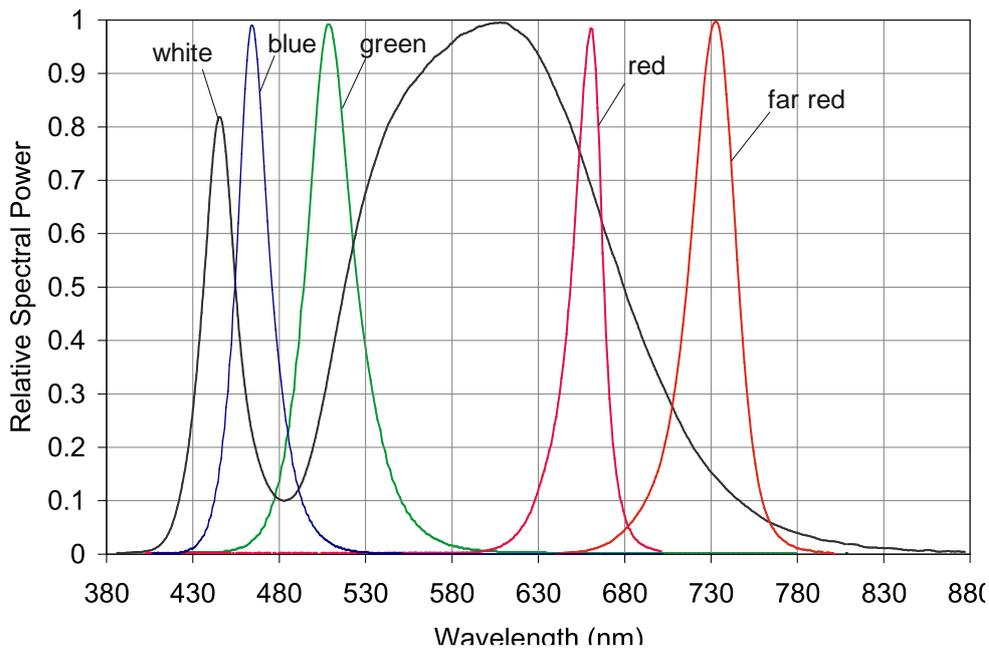
- Green panel (ID 56589-01)
 - Blue (470nm)
 - Green (520nm)
 - Red (660nm)
 - Far red (730nm)
- White panel (ID 56589-10)
 - Blue (470nm)
 - White (cold white)
 - Red (660nm)
 - Far red (730nm)



Make sure to wear standard safety glasses when using the LED panels due to the possible high intensity of the light source (CE EN 166, EN 172). These safety glasses ensure that bright, intense light intensity will not harm the beholders eyes.

The panels are coded for the left and the right hand side. They are fixed on the upper left and right side above the sample using the hooked fixing.

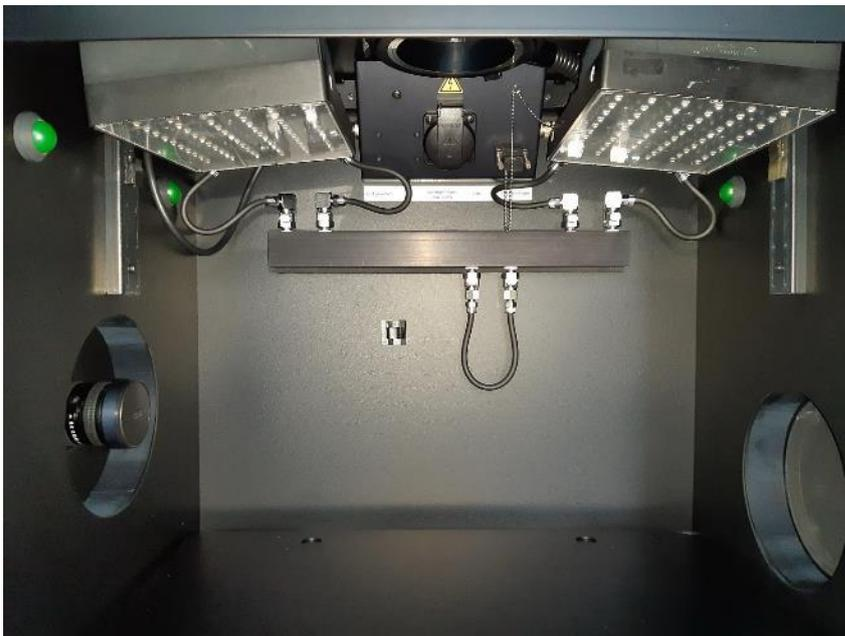
The spectra of the single LED channels are shown in the diagram below



The electrical and cooling liquid tubing connections inside of the dark box are shown in the image below:



Connect the LED panels only while the instrument is powered off to prevent electrical damage of the LED panel.



The control cable of each panel is plugged into the socket at the corresponding side of the connection box (see 5.3.3).

Water cooling connects to the water supply bar located at the rear side of the dark box. The plug at the lower side of the supply bar is a bypass plug (instead of the temperature controlled sample table) and must be present to ensure a flow of cooling water.

External connectors for the cooling liquid are located at the rear side of the miniChiller and the rear panel of the NightSHADE evo:



Tubing connector at the miniChiller and at the rear panel of the NightSHADE evo



When the miniChiller or any other device using liquids to cool the LED panel is connected to the NightSHADE evo it is not allowed to operate 100/230V devices such as transilluminators etc. inside the dark room due to electrical safety regulations.



The preset temperature of the cooling water in the miniChiller must not be set below ambient temperature since this could lead to condensation on the LED panel.

The "Artificial daylight" control must be enabled in the indiGO service backend in order to control the intensity of the LED panels.



To operate the cooling device please refer to the dedicated manual of the minChiller.

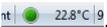
10.1.2 Temperature Controlled Measurement Table

The temperature controlled measurement table is used for measurements which require a temperature controlled environment.

If you are working with the heated macro table, plug the connector into the 9-pole Sub-D-connector on the power box at the center top inside the darkbox (see 5.3.3.) The device must be switched off, since after the NightSHADE evo is switched on the heating system is detected and may get initialized.

If the heating system is plugged in during ongoing operation, it will work only after the device has booted up again. This is done by clicking on  and selecting **<Boot Device>** from the **<Reference>** dropdown list.

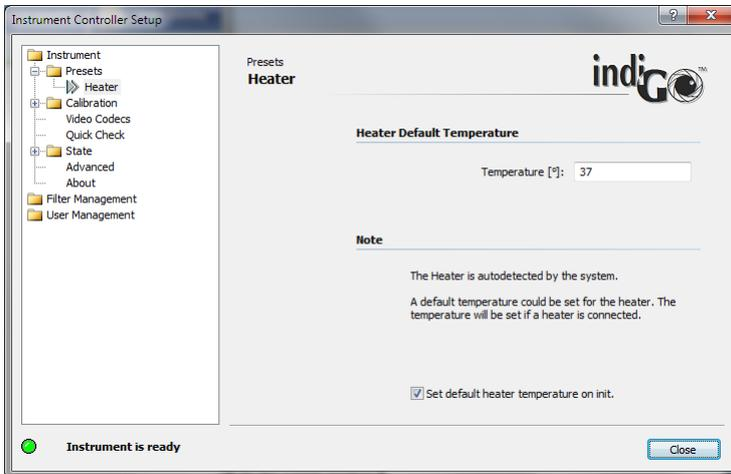


After approx. 5 to 10 minutes the heated measurement table has reached its working temperature (preset 37°C). The current temperature is displayed in the bottom status bar of the indiGO user interface .



The default temperature can be defined in the configuration backend of the software accessible through the instrument controller. Right click on the controller icon and log-in with a valid user account information selecting Setup from the context menu.

In the **<Presets>** **<Heater>** section the default temperature can be set. When the checkbox at the bottom is checked indiGO will set the default temperature at system initialization.



10.1.3 Temperature controlled tray with anti condensation



The temperature controlled tray can be placed inside the NightSHADE evo to allow a temperature controlled base plate when using petri dishes. The temperature is regulated by the water cycle used by the LED panels as well. Since the miniChiller can only cool down liquids it might be necessary to use the miniStat 125 or the miniChiller Plus when temperatures above or around ambient are required. The tube connections from chiller to NightSHADE evo are described in chapter 10.1.1.

A water cycle flows through the base plate of the tray which stabilizes the temperature. To avoid condensation of the petri dish lids an airflow is established over the dishes in a circular flow direction. This temperature should be about 3-5 degrees above the table temperature and can be set using indiGO (see 10.1.2).



The temperature controlled tray must be connected to the water cycle using the lower two connectors at the water supply bar. When the tray is not used, both outlets must be shorted (bypass) as shown below. If this is not done, the LED panels are not cooled properly due to an interruption in the water cycle. This could lead to an error state of the LED panels due to overtemperature.

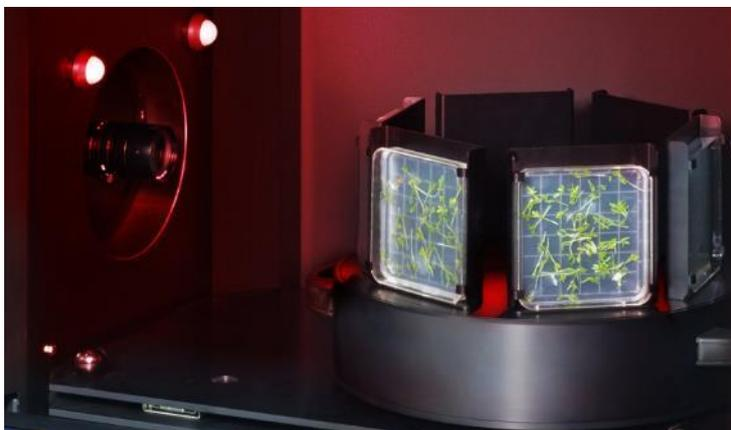


The temperature table itself is positioned inside the darkbox by the 4 feet. Make sure the 4 feet fit inside the corresponding wells located at the base plate.



Avoid temperatures who are significantly below ambient as this could lead to condensation.

10.1.4 Turntable



The turntable is used together with the sideview option. Using the turntable up to 6 samples can be measured in a automatic sequence. 3 different sample holders can be selected:

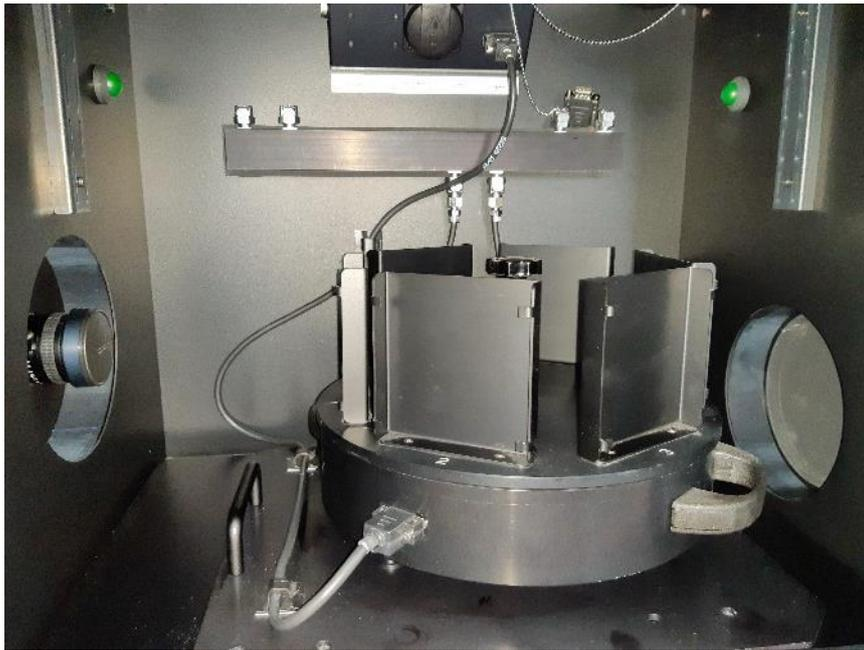
- 100mm square petri dishes
- 130mm square petri dishes
- DeWit culture tubes

Installation:

- Turn off the NightSHADE evo
- Place the turn table on the base plate that the rubber feet of the table will fit the positioning holes and markers on the positioning plate. The orientation of the turn table can be seen on the image below. Depending on the sample size the turn table can be lifted up to change the image center. In addition the positioning plate offers 2 different positions to adjust the sample size.

When using the 130mm petri dishes the turn table must be placed at the outer right position to ensure the whole plate can be imaged using the standard lens.

The 100mm petri dishes can be completely measured at the left position of the position plate. In this case the adjustable feet of the turn table must be pulled out to elevate the turn table (bajonet lock) to place the sample center in the center of the optical axis.



- Connect the 9 pin cable supplied with the turn table to the connector at the connecting panel at the upper position inside the dark box. Make sure the cable is not blocking the table or gets squeezed by the retractable base plate. Remove the dummy plug and keep it in a safe place. When not using the 9 pin connector, the dummy plug must be present
- Switch on the NightSHADE evo to initialize the turn table

Before removing the turntable make sure, the instrument is switched off.
The sample changer feature must be activated in the indiGO service backend in order to control the turn table



The position of the turn table is defined by the number of the sampleID. 001 represents the sample position 1, 006 moves the turntable to position 6. 007 will use position 1 again.

If further movement of the table is required a different position can be selected using a control sequence (see. Chapter 4.2.4)

10.1.5 Test Light Source

The Test Light Source is used to calibrate and test the System.



It contains a low light source which will glow for about 15 min. without external supply. To charge the unit, use the included mains adapter.

Charging is indicated by a red LED on the test plate. As soon as the plate is fully charged, the green LED will light. It takes about 1-2 minutes to charge the plate.

Now the plate is ready to be used in the NightSHADE evo. The spot with the white reflector (microplate position E9) emits light with an energy of approx 20nW (see individual certificate). This spot can be used to check the NightSHADE evo calibration. The measuring time should not exceed 1 second since the camera could be saturated otherwise.

Position E6 gives a lower signal intensity. This position can be used to check the detection limit with additional neutral density filters on top of the light source.

A special center plate is optionally available to place the E9 light spot at the center of the image.

The required light position is controlled through the sliding switch at the side of the plate (position E9 -> overl.; E6 -> 96). When the plate is not in use put the slider switch to position „OFF“.

10.2 Index

1D-line plot	63	filter wheel	82
acquisition	54	first time initialization	31
Admin	34	Flatfield Correction	43
Air Filter	116	Fluorescence Filters	35
Anti condensation	124	fuse	116
Application	41	Fuse	116
Apply to		Gain	42
Identity	61	Guest	35
Project	61	halogen lamp	117
Areas		Identities	51
Manual	66	Import	
Peaksearch	65	indiGO 1.x Data	104
Artificial daylight	122	WinLight Data	105
Background Correction	43	LED Panels	120
Backup Projects	104	Messages	108
Binning Factor	43	miniChiller	122
bottom plate	84	Noise	119
Camera Installation	23	Peaksearch	65
Cameras		Photo Properties	92
Ikon	82	Project	50
CAN socket	81	Quantum efficiency	119
C-Mount thread	115	Quick Check	54, 108
connector field	81	quick start	20
Connector Field	81	Range	93
Consignes de Sécurité	12	Read out	42
Contrast Slider	91	rear panel	25
Controller Template	47	Rear Panel	85
Controller Version	108	Resolution	119
Cosmic suppression	43	safety instructions	9
Customer service	19	Sample ID	51
Data Slider	94	Sicherheitshinweise	15
Database Connection	110	Side view camera	24
Device State	107	Software Installation	26
Door Sensor	116	Technical Data	119
Emission Filter	36	temperature controlled measurement	
Excitation Filter Slider	35	table	123
Exit	103	Test Light Source	127
Export		Toolbar	97
Batch Image Export	112	Top view camera	23
Excel	71	Transport	107
Report Settings	111	Turntable	125
Rich Text	70	typographical conventions	7
Exposure Time	42	User	34
Filter Manager	39	User Management	34
filter slider	83	Warranty	18

Modifications due to technical advancement reserved.