

# SINGLE PLATE ELISA WALKAWAY AUTOMATION

The Crocodile miniWorkstation



**BERTHOLD**

# SMALL FOOTPRINT. RELIABLE RESULTS. CONVENIENT ELISA AUTOMATION.

## Automated unattended operation for low to medium throughput



Running your ELISA assays without automation requires an hour or more of your attention, making it difficult to plan other important activities while your assay is being processed.

The Crocodile 5-in-one miniWorkstation provides complete ELISA automation in a footprint similar to a stand-alone-reader, while ensuring reliable results.

Using the Crocodile will reduce assay times by eliminating the need to move plates between dispenser, washer, incubator, shaker and reader. Free-up precious time to perform other important activities.



8 well parallel washer, providing consistent washing performance

### Crocodile benefits at a glance

- **All-in-one ELISA automation:** Truly unattended operation – simply insert a microplate including samples and standards, choose the assay protocol and leave the work to the Crocodile.
- **Ultra-compact footprint:** With its compact width of only 26.5 cm, the system fits into any laboratory environment.
- **Plug & play setup:** The system is pre-installed and ready to start within minutes.
- **Open system software:** The user-friendly software easily adapts to any 96 well assay.
- **Wavelength flexibility:** In addition to the preinstalled filters for ELISA (405, 450, 492 and 620 nm), 4 positions are available for custom filters.
- **Robust design:** The Crocodile is very robust due to the small number of moving parts.
- **Simple maintenance:** Maintenance is also really simple: flushing of the system and routine replacement of the tubing is made easy by the intuitive design of the Crocodile.
- **Easy validation:** Installation Qualification and Operation Qualification steps are performed according to detailed instructions provided (optional Qualification package).

# ALL-IN-ONE ASSAY AUTOMATION

Requires the footprint of a reader. Offers the functionality of up to 5 different instruments.

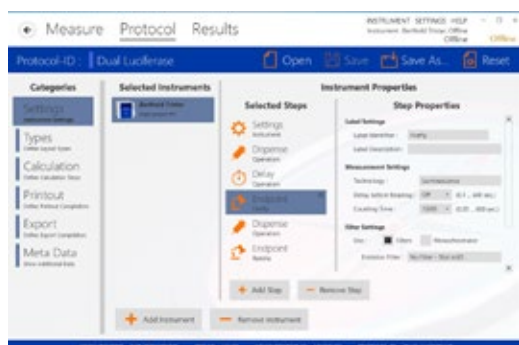
The Crocodile 5-in-one miniWorkstation provides the same functionality as five individual instruments in a footprint similar to a stand-alone reader.

1. Dispenser
2. Shaker
3. Incubator
4. Washer
5. Reader (5-in-one version only)



## TOTAL FLEXIBILITY. MAXIMUM PRODUCTIVITY.

Intuitive, easy-to-use software



### The LightCompass® Software

LightCompass® is a flexible software enabling easy data acquisition, analysis and reporting using Berthold Technologies' instruments. It has been optimized for different microplate and tube formats. LightCompass® features all the functions you can expect from an advanced scientific software, such as flexible creation of measurement protocols, curve fitting, qualitative analysis and compliance with FDA 21 CFR Part 11.

LightCompass® allows the easy import of your measurement data from the Crocodile Control Software to process it according to your assay needs. You can perform any desired complex calculation starting from a simple calculation of the CV up to more complex operations like curvefit and complete QC procedures all within a 21 CFR part 11 compliant environment. 3

# TECHNICAL SPECIFICATIONS

<b>General Specifications</b>	Sample format	1 × 96 well microplate (ANSI SLAS)
	Dimensions (H × W × D)	25.8 × 26.3 × 62.6 cm / 10.2 × 10.4 × 24.6 inch
	Weight	14.8 kg / 32.6 lbs (5-in-one); 13.9 kg / 30.6 lbs (4-in-one)
	Interface	USB interface
	Power Requirements	Desktop power supply and power cord. 100 – 240 V AC; 24 V DC / 3.3 A
	Operating Temperature	10 – 35 °C; 50 – 95 °F
	Humidity	Up to 75 % (non condensing)
<b>Incubator</b>	Temperature Range	From ambient +4 °C to 55 °C
	Temperature Stability	±1 °C across plate at 37 °C
	Incubation Time	Programmable
	Temperature Monitoring	Yes
<b>Reader (5-in-one version)</b>	Dynamic Range	0 – 3.0 OD at 450 nm
	Spectral Range	400 – 690 nm (pre-installed filters: 405, 450, 492, 620 nm)
	Filter Slots	8
	Reading Channels	8 independent photometric channels plus reference channel, mono and bichromatic reading
	Precision	<1.5 % CV (0.01 to 0.5 OD) <1 % CV (0.5 to 2.0 OD) <1.5 % CV (2.0 to 3.0 OD)
	Accuracy	±0.01 OD or 2.5 %, (whichever is greater)
<b>Dispenser</b>	Type	4 independent precision pumps
	Volume Range	10 – 2000 µl
	Precision	<2 % CV at 100 µl
	Accuracy	<5 % at 100 µl
	Reagent Support	Any bottle or vial type can be attached
<b>Washer</b>	Manifold Configuration	Autoclavable 8 way manifold with 2 parallel needles
	Dispense Volume	25 – 1000 µl
	Wash Modes	Standard, overflow, soak, two-point, sweep
	Wash Buffer Capacity	3 × 1 l bottles for different wash solutions
	Waste Liquid Container	2 l bottle
<b>Shaker</b>	Shaking	Independent linear motion 5 – 20 Hz
<b>PC Software</b>	PC Software	Crocodile Control Software for Windows allows creating and measuring assay protocols
<b>Ordering Infos</b>	84024-01	LB 925 Crocodile miniWorkstation 5-in-one incl. Crocodile Control Software licence
	84024-02	LB 925 Crocodile miniWorkstation 4-in-one incl. Crocodile Control Software licence
	37854-403	LightCompass® Professional
	37854-404	LightCompass® Plus
	67950	Crocodile Control PC-Software (CCS) (additional licence)
	84078	IQ/OQ/PQ Qualification Package for Crocodile miniWorkstation
	84083	Absorbance Test-plate for validation of 96 well microplate absorbance readers

## TRANSFORMING SCIENCE INTO SOLUTIONS

Berthold Technologies is a global technology leader in life sciences. Our extensive range of analytical system solutions made in Germany has been trusted by scientists since 1949. These range from small standalone readers, such as microvolume spectrophotometer and luminometers to various dedicated and multimode readers, microplate washers, microplate workstations, RIA and ELISA automation products to high-end imaging systems, HPLC radio detectors and gamma-counters. It is our mission to create a healthier world, a safer environment and more efficient manufacturing processes.

### Berthold Technologies GmbH & Co. KG

Calmbacher Straße 22 · 75323 Bad Wildbad · Germany  
+49 7081 1770 · bio@berthold.com · [www.berthold.com/bio](http://www.berthold.com/bio)

© Berthold Technologies. All rights reserved. All trademarks are the property of Berthold Technologies or their respective owners. Berthold Technologies reserves the right to implement technical improvements and/or design changes without prior notice.

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.