

ROD SOURCE SHIELDS LB 8300

Operating Manual



69560BA2 Rev.01 04/2023

BERTHOLD

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

industry@berthold.com

Table of contents

1	About this operating manual	5
1.1	Applicable documents.....	5
1.2	Some prior remarks	5
1.3	Storage	5
1.4	Target group.....	5
1.5	Validity of the manual	6
1.6	Structure of the manual.....	6
1.7	Copyrights	6
1.8	Representation	6
1.9	Symbols used.....	7
1.10	Structure of warnings.....	8
1.11	Symbols used on the device.....	9
2	Safety	10
2.1	Proper use	10
2.2	Ambient conditions during operation and storage.....	11
2.3	Qualification of the personnel	11
2.4	Radiation protection	13
2.4.1	Basic principles and regulations	13
2.5	Emergency procedure	15
2.7	Source type plate.....	16
3	System description	17
3.1	View.....	18
4	Transport.....	20
4.1	Safety instructions	20
4.2	Packaging.....	21
4.3	Temporary storage of the source	21
4.4	Scope of delivery	21
4.5	Transport to operation site.....	21
5	Installation	22
5.1	Safety instructions	22
5.2	Ambient conditions during installation.....	23
5.3	Temporary storage of the source	23
5.4	Preparing the installation	23
5.5	Checking the delivery.....	23
5.6	Installation of the shield	24
5.7	Marking.....	29
6	Usage	30
6.1	Safety instructions	30
6.2	Opening and closing the locking mechanism	31
7	Maintenance and repair.....	33
7.1	Safety instructions	33
7.2	Visual inspection.....	34
7.3	Leak test.....	34
7.4	Shutter check procedure	36
7.5	Removing and installing the source.....	37
7.5.1	Removing the source.....	37
7.5.2	Installing the source	40
7.6	Source replacement.....	42
7.6.1	Replacing the source	43

8	Uninstallation.....	44
8.1	Safety instructions	44
8.2	Dismounting the shield	45
8.3	Disposal of radioactive substances	46
8.4	Shipment of radioactive substances.....	46
9	Appendices.....	48
9.1	Check lists for tests	48

1

About this operating manual

1.1 Applicable documents

This manual contains the following document:

- Technical Information, Id. No. 69560TI (see appendix)

1.2 Some prior remarks

The product is handed over to you by the manufacturer BERTHOLD TECHNOLOGIES GmbH & Co. KG (designated as Berthold in the following) in a complete and functionally reliable condition.

This operating manual illustrates how to:

- set up/install the product
- operate the product
- carry out maintenance on the product
- dismantle the product
- dispose of the product

Read these instructions thoroughly and completely before working with the product. We have tried to compile all the information for safe and proper operation for you.

However, should questions arise which are not answered in this manual, please contact Berthold.

Store the instructions where they are accessible for all users at all times.

1.3 Storage

This operating manual as well as all product-related documentation relevant to the respective application must be accessible at all times during the life cycle near the device.

1.4 Target group

The product may only be installed, operated, maintained and repaired by trained personnel.

This manual is directed at qualified specialist personnel who are familiar with handling radioactive sources and heavy system components.

Specialist personnel refers to those who can assess the work assigned to them and recognise possible dangers through their specialist training, knowledge and experience as well as knowledge of the relevant regulations.

1.5 Validity of the manual

The manual is valid from the delivery of the Berthold product to the user until its disposal. Version and release date of this operating manual can be found in the bottom of each page. An alteration service is not provided by the manufacturer Berthold.

The manufacturer reserves the right to make changes to this operating manual at any time without stating reasons.

NOTICE



The current revision of this operating manual replaces all previous versions.

1.6 Structure of the manual

This manual has been divided into chapters. The order of the chapters should help you to familiarise yourself quickly and properly with the operation.

1.7 Copyrights

This manual contains copyright-protected information. None of the chapters may be copied or reproduced in any other form without prior authorisation from the manufacturer.

1.8 Representation

Identifier	Meaning	Example
Round brackets	Image reference	Connect the plug (fig. 1, item 1)
	Prohibited actions, procedures or processes within a figure.	
	Representation of the ionizing radiation within a figure.	

1.9 Symbols used

NOTICE



If this information is not observed, deterioration in the operation and/or property damage may occur.

IMPORTANT



Sections marked with this symbol point out important information on the product or on handling the product.

Tip



Provides tips on application and other useful information.



General warning symbol



Warning symbol danger of crushing



Warning symbol heavy loads



Warning symbol suspended load



Warning of radioactive substances



Wear protective helmet



Wear safety shoes



Do not touch the surface

1.10 Structure of warnings

Signal word



Source and consequence

Explanation, if required

- ▶ Measure

In case of emergency

-
- **Warning symbols:** (warning triangle) draws attention to the hazard
 - **Signal word:** states the severity of the hazard
 - **Source:** states the type and source of the hazard
 - **Consequence:** describes the consequences if warning is ignored
 - **Measure:** states how one can avoid the hazard.
 - **In case of emergency:** states how to react in case of direct danger.

Warning levels used

In this manual, warning instructions in front of instructions for action refer to risks of injury or damage to property. The hazard-prevention measures described must be observed.

DANGER



Indicates an **imminent**, major hazard, which will certainly result in serious injuries or even death if the hazard is not avoided.

WARNING



Indicates a **potential** hazard, which can result in serious injuries or even death if the hazard is not avoided.

CAUTION



Refers to a **potentially dangerous** situation, which can result in medium or minor physical injuries or damages to property, if it is not avoided.

1.11 Symbols used on the device



Ionising radiation

The shield is equipped with a radioactive source. Please note the handling instructions. Please observe the transport instructions in this operating manual.

2 Safety

2.1 Proper use

The source with shield is used in connection with a detector and a suitable evaluation unit provided by Berthold to measure the radiation intensity occurring during a radiometric measurement.

The shield was developed as shield/protective container for radioactive sources and may be used exclusively for this purpose.

Usually, the shield contains a radioactive source. The notes on radiation protection contained in the present manual as well as any statutory requirements in this respect are to be strictly adhered to.

The following constitutes proper use

- Adhering strictly to the instructions and operation sequences and not undertaking any different, unauthorised practices which could put your safety and the operational reliability of the shield at risk.
- Observing the provided safety instructions.
- Carrying out the prescribed maintenance measures or having them carried out for you.

The following constitutes improper use and is to be avoided

- Any non-compliance with the present operating manual for the supplied products
- Applying conditions and requirements which do not conform to those stated in the technical documents, data sheets, operation and assembly instructions and other specific guidelines of the manufacturer.
- Using the product after any repair carried out by employees who have not been authorized by Berthold.
- Using the product in a damaged or corroded condition.
- Dismounting the unit while the radiation beam outlet is in the open position (except for situations in which the locking mechanism is defective and the beam outlet can no longer be closed).
- Operation without the safety precautions provided by the manufacturer.
- Any modification to design and function, except for any activities provided for and described in the present manual.
- Restructuring or changing the system components.
- Manipulation or avoidance of existing safety equipment.

Berthold shall only accept liability for / guarantee the correspondence of the product to its publicised specifications.

If the product is used in a way which is not described in this manual, the product's protection is compromised and the warranty becomes void.

2.2 Ambient conditions during operation and storage

The shield was specifically designed for use in rough ambient conditions. The compliance with the operating conditions specified below contributes to guaranteeing the permanent functionality of the shield and the prevention of damage.

Shields containing radioactive substances and sources are to be stored in a lockable storage room complying with the national requirements for the storage of radioactive substances.

The higher the dust and dirt content of the environment, the more likely the rotational motion of the shutter mechanism can be impaired or entirely blocked. For this reason, the functional test intervals (see chapter 7) should be adjusted for the ambient conditions.

Highly combustible or explosive substances must not be kept in the vicinity of shields in order to prevent a fire from spreading to the radioactive substances. Furthermore, the ambient condition requirements in the document "Technical Information" are to be observed.

2.3 Qualification of the personnel

NOTICE



A minimum requirement for all work on or with the product would be employees with general knowledge who are instructed by an expert or authorised person.

At different parts in this manual, reference is made to personnel with certain qualifications who can be entrusted with different tasks during the installation, usage and maintenance.

The four groups this refers to are:

- Employees with general knowledge
- Experts
- Authorised persons
- Radiation Safety Officer

Employees with general knowledge

NOTICE



Employees with general knowledge must always be guided by one expert at the very least. When dealing with radioactive substances, a radiation safety officer must also be consulted.

Employees with general knowledge are e.g. technicians or welders who can undertake different tasks during the transportation, assembly and installation of the product under the guidance of an authorised person. This may also refer to construction site personnel. The persons in question must have experience in handling the product.

Experts

Experts are people who have sufficient knowledge in the required area due to their specialized training and who are familiar with the relevant national health and safety regulations, accident prevention regulations, guidelines and recognised technical rules.

Expert personnel must be capable of safely assessing the results of their work and they must be familiar with the content of this manual.

Authorised personnel

Authorised personnel are those who are either designated for the corresponding task due to legal regulations or those who have been authorised by Berthold for particular tasks. When dealing with radioactive materials, a radiation safety officer must also be consulted.

Radiation safety officer

In order to ensure proper handling and compliance with the statutory requirements, the company has to appoint a radiation safety officer in accordance with the applicable national law (in Germany: Strahlenschutzverordnung [German radiation protection regulation]). The radiation safety officer must implement the statutory radiation protection requirements in order to protect employees against damage to their health caused by handling radioactive materials.

NOTICE



Dangerous goods officers must not perform any activities as radiation protection officers, unless they underwent a special training as radiation safety officer!

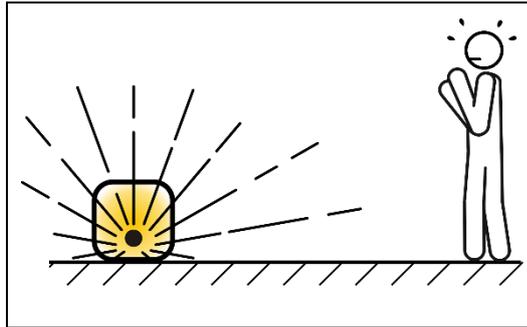
Radiation safety officers must have a special training with attendance of an officially recognized course and appropriate professional experience.

2.4 Radiation protection

2.4.1 Basic principles and regulations

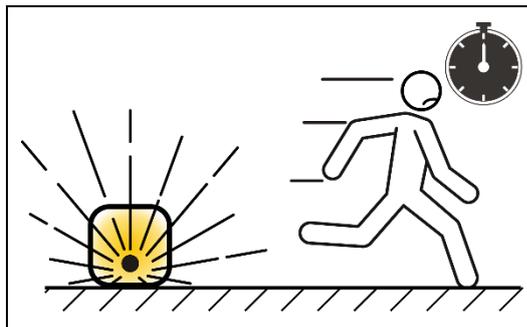
The amount of radiation absorbed by the body (exposure to radiation) is determined by three parameters from which the basic radiation protection regulations can be derived:

Distance



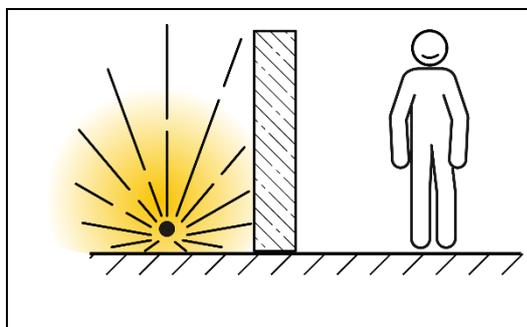
If work close to equipment containing radioactive substances is required, the largest distance possible is to be kept. In particular, this applies to employees not immediately required to work directly with the radioactive devices.

Time



Any work required in the vicinity of radiometric measuring systems is to be prepared for carefully and to be organised in a way that the work can be executed as quickly as possible. Here, providing the correct tools and aids is particularly important.

Shield



When mounting and dismounting the shield, it is to be ensured in advance that the radiation beam outlet is closed.

Exposure of employees to radiation

During installation, maintenance and decommissioning of the shield, employees may be exposed to radiation.

In order to keep such exposure as low as possible, the shield with the source may only be mounted and/or dismantled by authorized employees. Such authorised staff is to be instructed as regards all rules of behaviour when handling radioactive substances in advance.

It is to be ensured that the locking mechanism of the shield is closed and secured in order to prevent the emission of unshielded radiation. Modification or damage to the shield must be avoided at all times.

Work may only be executed according to the instructions and under the supervision of the radiation safety officer, who furthermore has to calculate or estimate the exposure of the employees to radiation in order to ensure that the statutory dose rate limits are not exceeded.

Theft protection

Radioactive substances or equipment containing radioactive substances must be secured in a way that they are protected against access by unauthorized persons. In the case of firmly installed equipment containing radioactive substances, the protection against unauthorized access is generally provided by the firmly attached installation.

Shields with radioactive sources which are decommissioned for a certain period of time must be dismantled and securely stored in a storage room complying with the national regulations for the storage of radioactive substances.

Portable measuring systems must never be left unsupervised. When this equipment is not in use, it is to be protected against access by unauthorised persons.

In the case of fire

The shielding material can melt and leak from the shield if exposed to very high temperatures for an extended period of time. During and after a fire, there is a risk of major long-time consequences for your health due to the incorporation of lead as well as a risk of increased radiation exposure.

When planning the use of radiometric measuring systems, constructional measures ensuring fire prevention are to be provided.

- ▶ In the case of fire, these measures limit the access to this area.
- ▶ Avoid the incorporation, contamination and exposure by keeping sufficient distance.
- ▶ Notify Berthold of the situation; you will promptly receive information on immediate measures.

2.5 Emergency procedure

In case of serious operational trouble, such as fire or explosion, which could adversely affect the radiometric device, it cannot be ruled out that the function of the shielding lock, the shielding efficiency or the stability of the source capsule have been compromised. In this case, it is possible that people in the vicinity of the shield have been exposed to higher levels of radiation.

If you suspect such a severe malfunction, the Radiation Safety Officer has to be notified immediately. He will then investigate the situation immediately and take all necessary actions to prevent further damage and to avoid more exposure of the operating staff to radiation.

The Radiation Safety Officer has to make sure that the measuring system is no longer in operation and then take appropriate steps. He may have to inform the authorities or contact the manufacturer or supplier of the measuring system.

If adequate know-how as well as suitable instruments are available, emergency measures may be taken immediately. In this case, proceed as follows:

1. Locate the shielding.
2. Check the function of the shielding.
3. Check the efficiency of the shielding by measuring the dose rate.
4. Secure and label radiation protection areas.
5. Secure the shielding with source.
6. Document the event and estimate the possible radiation level to which the people involved were exposed.

If you suspect any damage to the source capsule, the following points have to be observed as well:

1. Avoid contamination.
2. Take hold of source using a tool (a pair of pliers or a pair of tweezers) and put both (source and tool) into a plastic bag.
3. Secure them behind an auxiliary shielding (concrete wall, steel or lead plate).
4. Check if the environment is free of contamination.
5. Make sure the radioactive waste is secured and disposed of in compliance with the pertinent regulations.

2.6 Operator's obligations

The operator of the product must regularly train his personnel in the following topics:

- Observation and use of the operating manual and the legal provisions.
- Proper use of the product.
- Observation of the plant security instructions and the operating instructions of the operator

2.7 Source type plate

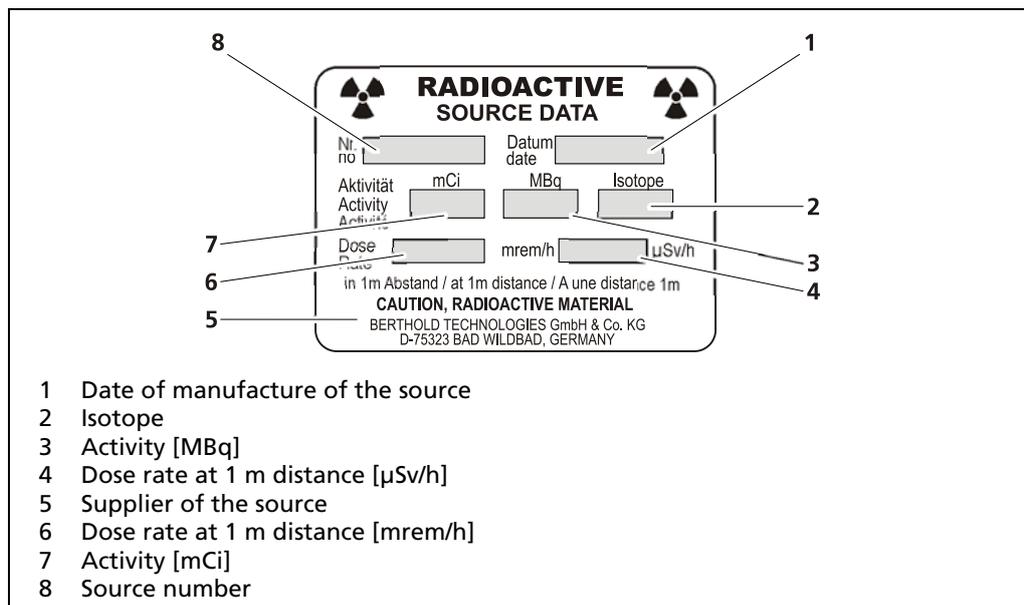


Fig. 1 Type plate of the source

3

System description

The rod source shield is intended as shield/protective container for radioactive rod sources. The radioactive substance is contained in a leak-proof welded source capsule and installed in the rod source shield. Apart from source and shield, additional system components such as detectors and evaluation units are required for a complete measuring system. Usage of these system components is not the subject of this operating manual. Please refer to the individual manuals of the respective system components.

The shield consists of a robust steel housing holding a rotatable cylinder. The rotatable steel cylinder is filled with lead except for the very center which holds the radioactive rod source and a narrow channel along the active length of the source - the radiation beam path. The provided lever rod can be used to rotate the cylinder and its beam path 90 degrees between two positions. The closed position has the beam path pointed towards a vertical lead filled section of the shield frame, blocking the radiation coming through the beam path. The open position has the beam path pointed towards the process unit and detector.

The locking cylinder can be locked in both rotational positions (radiation beam path open, radiation beam path closed) using fixing screws or a padlock.

The locking plate on the upper side of the shield ensures that the source cannot be removed by unauthorized persons. Operators can secure the locking plate by installing a padlock at the right-hand and the left-hand side.

Only the provided bolt holes in the lower mounting plate, upper locking plate (after removing the lifting eye nuts) and vertical side bars can be used to mount the shield!

The vertical side bars and the upper locking plate bolt holes can optionally be used for tilt protection.

The rod source shield is available in different versions. You can find an overview of all variants in the appendix (document "Technical Information" chapter 3).

The Shield has the following functions:

- Shield radiation to a level that is non-hazardous for operating personnel
- Locking the radiation beam path in the CLOSED position for transport and during installation
- Protecting the integrated source capsule from mechanical damage and the effects of the environment

3.1 View

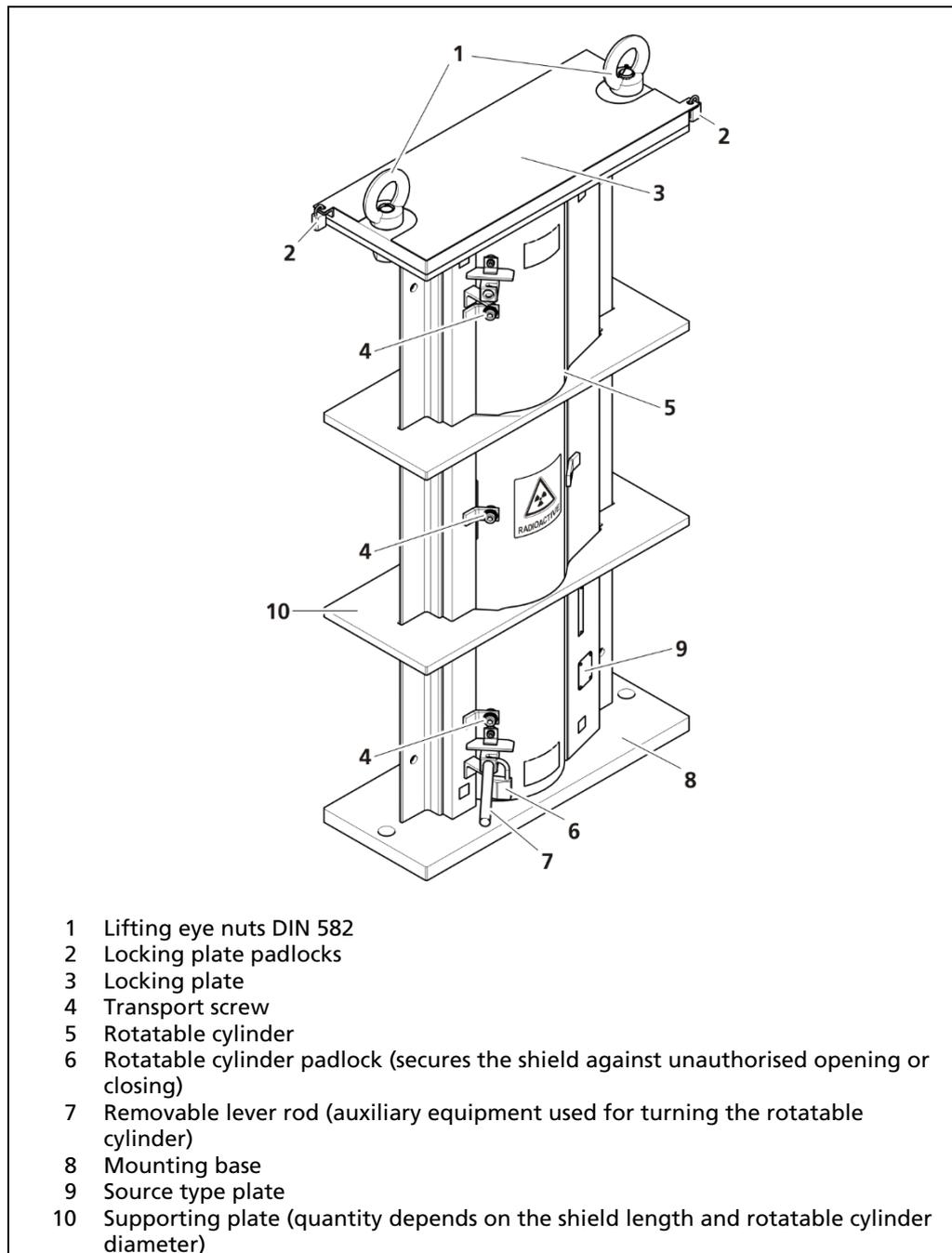


Fig. 2 Basic layout

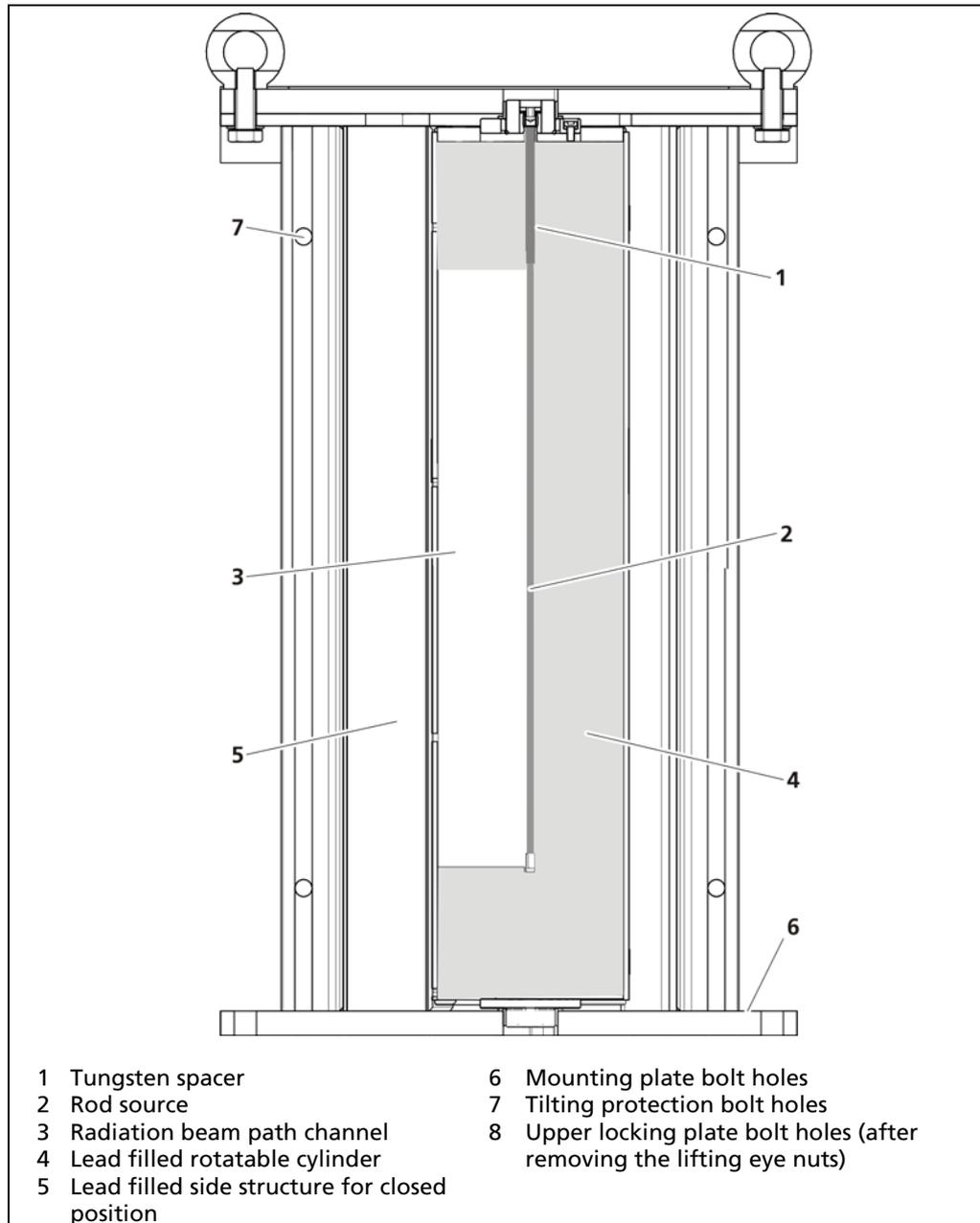


Fig. 3 Sectional view (position CLOSE)

4 Transport

4.1 Safety instructions

WARNING!



Danger of injury by falling loads!

Shields have built-in lead components increasing the weight.



- ▶ Only use the lifting eye nuts provided on the shield to attach the lifting gear.



- ▶ Never stand underneath a lifted or suspended load, keep at a safe distance.



- ▶ Wear head protection and safety shoes.

WARNING!



Danger of injury caused by heavy and bulky system components!

- ▶ Heavy and bulky system components should only be handled using aids and by at least 2 people.
- ▶ Observe the guidelines for safe handling of heavy loads.
- ▶ Ensure stability and use the provided fixing options.

CAUTION!



Danger caused by ionising radiation!

Shields usually contain radioactive sources. An increased exposure to radiation may lead to damage to health.

- ▶ Consult the radiation safety officer responsible for your company.
- ▶ Transport the source exclusively inside the closed and secured shield.
- ▶ The locking mechanism must be in the **CLOSED** position and secured during transport and mounting of the shield.

NOTE



The device may only be transported by competent people (see 2.3 Qualification of the personnel).

The applicable national regulations of the country of use have to be observed. Observe the marking for the center of gravity on the outer packaging, if applicable.

4.2 Packaging

Shields with the radioactive source is delivered in a packaging which complies with the regulations for the transport of radioactive material (Type A packaging).

4.3 Temporary storage of the source

If the source must be temporarily stored at the site of use between delivery and installation, please observe the following notes:

- Store the source exclusively inside a closed and secured shield.
- Store the shield in a lockable and properly marked room. The storage room must comply with the national requirements regarding the storage of radioactive substances.
- Accessible areas of increased radiation exposure must be marked and closed off, if required.

4.4 Scope of delivery

Check the delivery for completeness (material parts) and integrity. In the case of damage, immediately notify the forwarding agent and the manufacturer.

Check delivery

Document any visible transport damage with photos on delivery. In case of damage contact the transport company and the manufacturer immediately.

4.5 Transport to operation site

The transportation may only be executed by construction site personnel experienced in handling heavy components. The construction site personnel is to be instructed by at least by one authorised person. If the shields contains a radioactive source, the radiation protection officer is to be consulted, as well!

For lifting packages or shields weighing more than 25 kg, suitable aids (e.g. forklift truck) are to be used. If shields are lifted without transport packaging (wooden box or pallet), the provided lifting eyes are to be used exclusively for attaching the sling gear.

Please observe the radiation protection regulations for the transportation of radioactive substances.

5 Installation

The shield may only be installed and mounted by competent people (see chapter 2.3 Qualification of the personnel). Possibly the radiation protection officer is to be consulted, as well.

5.1 Safety instructions

WARNING!



Danger of injury by falling loads!

- ▶ Only use the lifting eyes provided on the shield to attach the lifting gear.
- ▶ Never stand underneath a lifted or suspended load, keep at a safe distance.
- ▶ Only use tested sling gear components appropriate for the transport weight.
- ▶ Observe the marking for the centre of gravity on the outer packaging, if applicable.
- ▶ Wear head protection and safety shoes.



WARNING!



Danger of injury caused by heavy and bulky system components!

- ▶ Heavy and bulky system components should only be handled using aids and by at least 2 persons.
- ▶ Observe the guidelines for safe handling of heavy loads.
- ▶ Ensure stability and use the provided fixing options.

CAUTION!



Danger caused by ionising radiation!

Shields usually contain radioactive sources. An increased exposure to radiation may lead to damage to health.

- ▶ Consult the radiation safety officer responsible for your company.
- ▶ Transport the source exclusively inside the closed and secured shield.
- ▶ The locking mechanism must be in the **CLOSED** position and secured during transport and installation of the shield.

IMPORTANT



The applicable national regulations of the country of use have to be observed.

5.2 Ambient conditions during installation

Mount shield using the lifting eyes (Fig. 2, item 1). The minimum permissible temperature during installation is **-20 °C**.

5.3 Temporary storage of the source

If the source must be temporarily stored at the site of use between delivery and installation, please observe the following notes:

- ▶ Store the source exclusively inside a closed and secured shield.
- ▶ Store the shield in a lockable and properly marked room. The storage room must comply with the national requirements regarding the storage of radioactive substances.
- ▶ Accessible areas of increased radiation exposure must be marked and closed off, if required.

5.4 Preparing the installation

NOTE



The shutter must be turned and secured in the **CLOSED** position during transport and installation of the shield.

IMPORTANT



Size and position of the applicable measuring range are specified in the project planning phase and determined by means of drawings, sketches or written notes. During installation, these specifications must be strictly observed since deviations may lead to malfunction of the measuring system.

To avoid an unnecessary exposure to radiation, the handling period of the source (even if it is shielded) must be kept as short as possible during installation. It is therefore recommended to perform the following steps prior to delivery of the source:

- Installation planning (including estimation of radiation exposure)
- Instruction of the employees
- Organising and cleaning the installation site
- Preparing the required tools and hoisting devices
- Removing all obstacles which might hinder the installation of the shield or the source

5.5 Checking the delivery

1. Check the delivery for completeness using the packing list.
2. Clean the parts, if required
3. In the case of damage, immediately notify the forwarding agent and the manufacturer.

5.6 Installation of the shield

WARNING!



Risk of injury due to tilting shield!

There is a risk of injury to people when working on non-fixed shields.

- ▶ Always secure upright shields against falling by fasten securing straps to the lifting eyes.

NOTE



Malfunctions and / or damage to the shield due to impermissible ambient conditions!

The mounting location and mounting position are determined during the project planning and defined by drawings, sketches or written instructions. These instructions must be observed strictly during assembly.

- ▶ Also note the information in the document "Technical Information" (see Technical Information - chapter 3).

To avoid unnecessary exposure to radiation, install the shield as final system component. The shield is placed on a mounting base at the measuring point and fixed. The mounted shield is secured against tilting by means of tipping protection.

Tip



If the shield contains a radioactive source, the measurement of the background count rate of the detector should take place before the shield is installed.

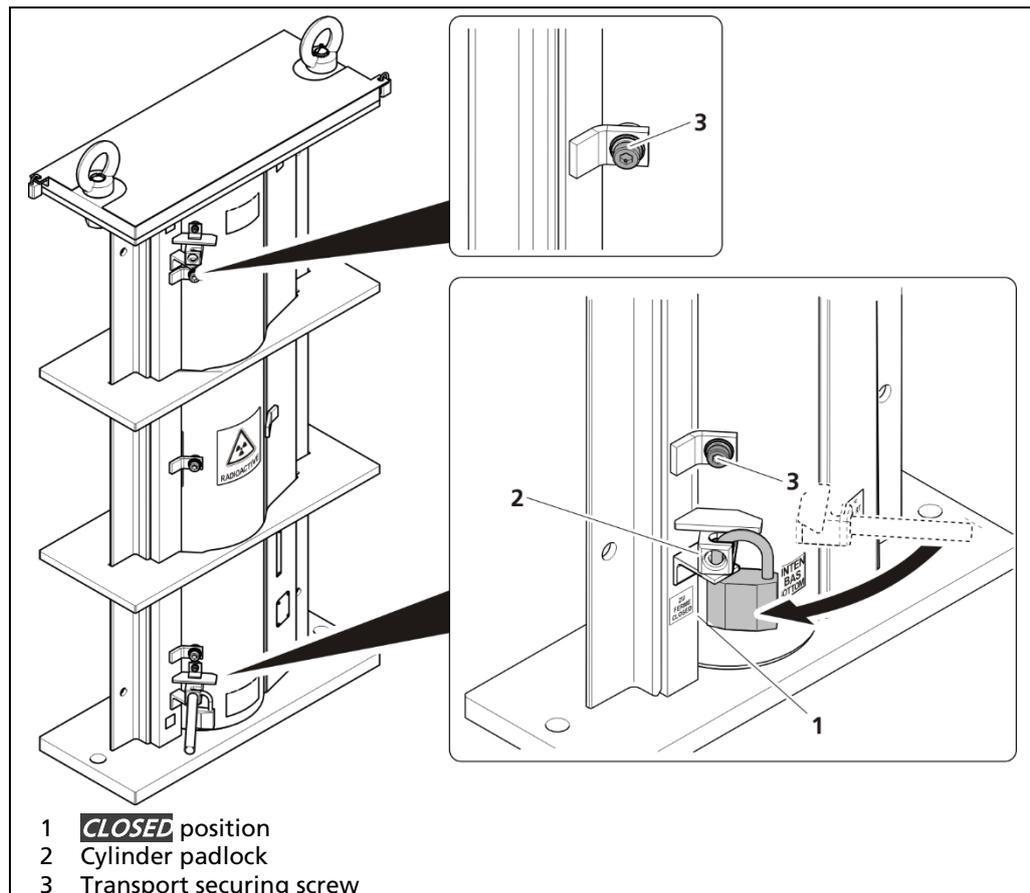


Fig. 4 Shield closed (**CLOSED** position)

1. Ensure that the shield has been closed and secured
 - ▶ the rotatable cylinder is set to the **CLOSED** position
 - ▶ the transport screw and, if applicable, the cylinder padlock are in place.

⚠ WARNING!



Danger of injury by falling loads!

Falling loads can endanger people.

- ▶ Never stand underneath a lifted or suspended load, keep at a safe distance.
- ▶ Cordon off the danger zone with barrier tape.

⚠ WARNING!



Danger of crushing by moving the shields!

Due to the high weight of the shield, there is a risk of injury.

- ▶ Heavy and bulky system components should only be handled using aids and by at least 2 people.
- ▶ Do not reach between the shield and the container or mounting base.

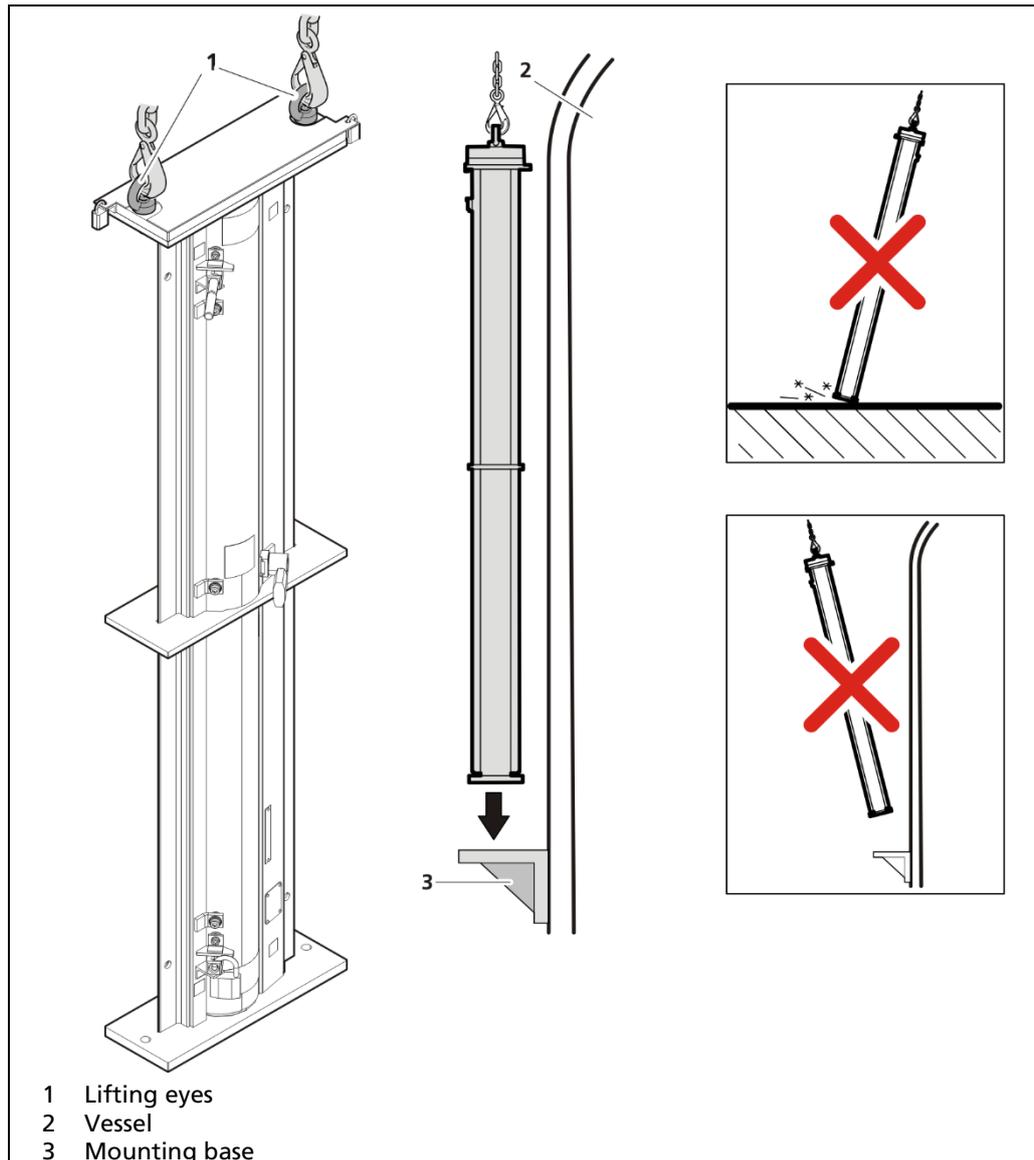


Fig. 5 lift and lower the shield

2. Fasten the shield with sufficiently dimensioned lifting straps at the lifting eyes (Fig. 5, item 1).
3. Raise the shield and place it vertically in the mounting position.
4. Carefully place the shield on the mounting base (Fig. 5, item 3).
5. Make sure the cylinder padlock side faces out and the beam path side faces the vessel.

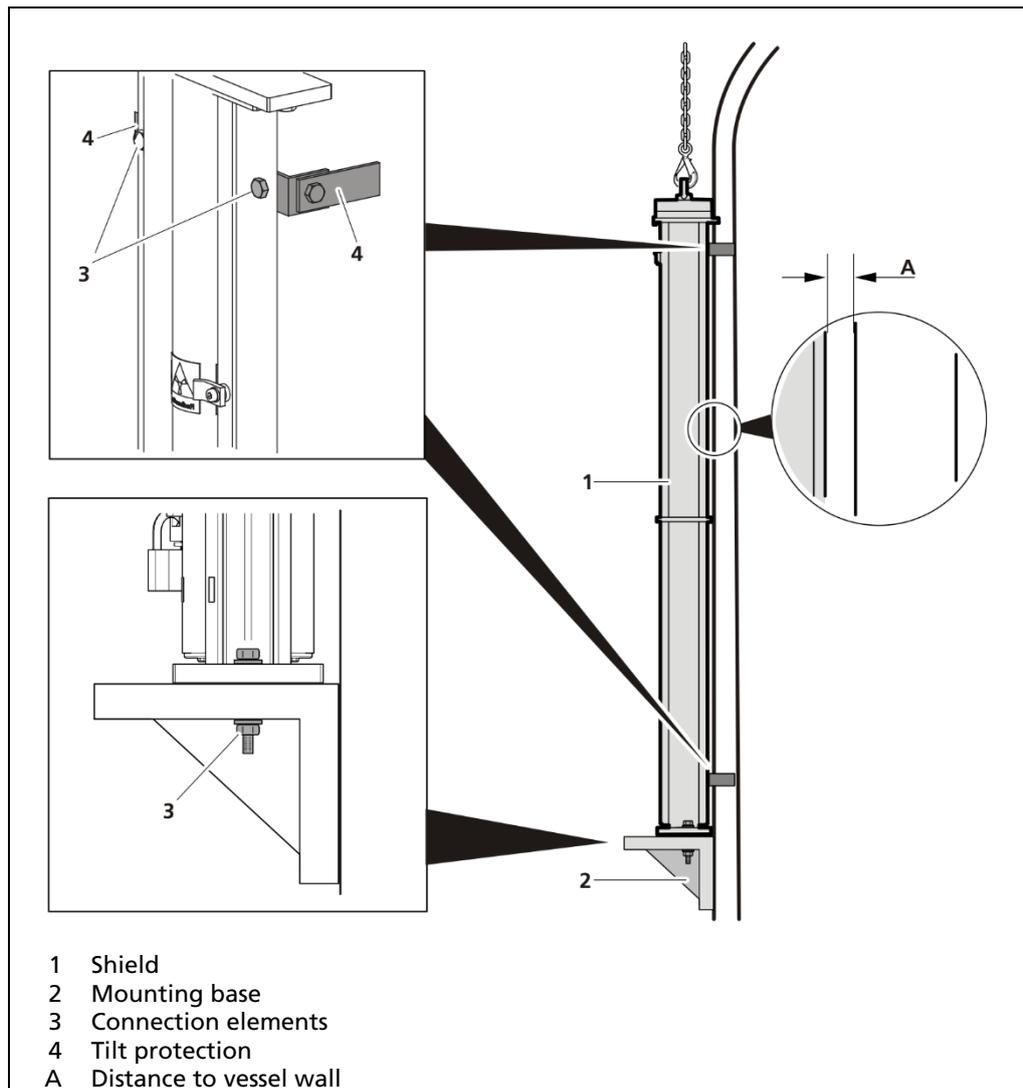


Fig. 6 Installation on vessels

⚠ WARNING!**Incorrect assembly may result in accidents!**

Incorrect selection of the fastening elements endangers the stability of the shield.

- ▶ Use correctly sized and approved mounting elements (mounting base, anti-tilt, nuts and bolts, screw lock).

6. Fasten the shield (Fig. 6, item 1) to the mounting base (Fig. 6, item 2). The plant operator is responsible for selecting the connection elements (Fig. 6, item 3), the screw locking and the tightening torque.
7. Secure the shield to the container wall with the anti-tilt device on both sides (Fig. 6, Pos.4). The plant operator is responsible for selecting the connection elements (Fig. 6, item 3), the screw locking and the tightening torque.

NOTICE

In the design of the shield it was assumed that the shield is mounted with a distance A of 20 mm, see Fig. 6.

NOTICE

Shields can be stacked or staggered when a larger measuring range is needed than one shield can provide. Please refer to the chapters 4.1 and 4.2 "Mounting Variants" in the document "Technical Information".

- ▶ The installation is complete.

5.7 Marking

The radiation warning sign (Fig. 7, item 1) indicates the beginning of the control area as long as the control area is outside the shield.

If the control area is within the shield, then the radiation warning sign (Fig. 7, item 2) already applied to the shield is sufficient. Fig. 7 shows an exemplary identification of the control area according to German law. Please observe the local legislation of the country of install.

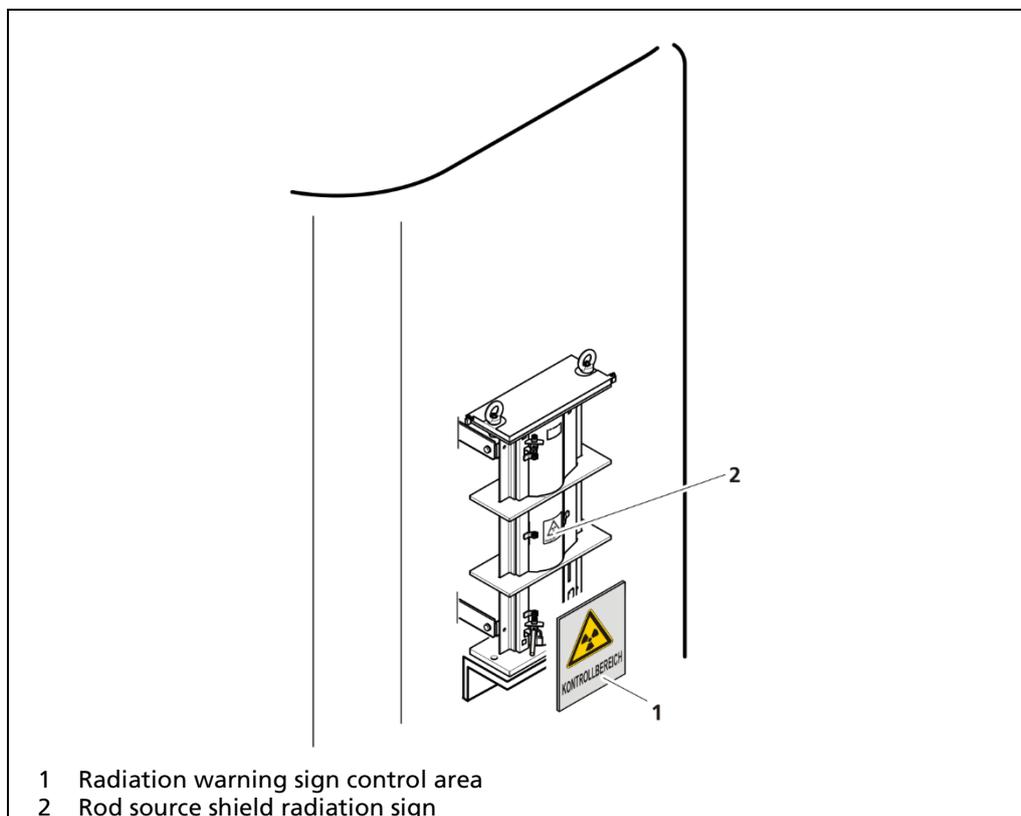


Fig. 7 Example marking

6 Usage

The shield may only operate by competent people (see 2.3 Qualification of the personnel). When operating the shield, unexpected changes at the process control system can occur. Before opening the locking mechanism, the guidelines of the system operator must be observed.

6.1 Safety instructions

CAUTION!



Danger caused by nuclear radiation!

Shields usually contain radioactive sources. An increased exposure to radiation may lead to damage to health.

- ▶ Ensure that the shield does not show any signs of damage or functional limitations.

6.2 Opening and closing the locking mechanism

The locking mechanism is used for opening and closing the beam path.

NOTICE



During transport and installation of the shield, the rotary cylinder must be set to **CLOSED** and be secured.

Opening the locking mechanism

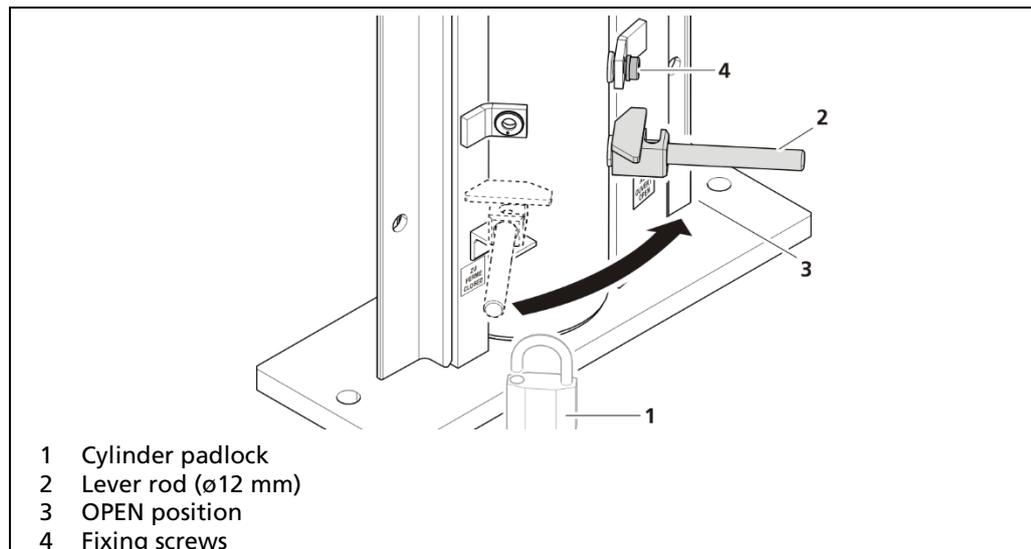


Fig. 8 Shield open (**OPEN** position)

1. If applicable, remove the cylinder padlock.
 2. Remove the fixing screws.
 3. Turn the rotary cylinder to the **OPEN** position with the help of the lever rod.
 4. Secure the rotatable cylinder in the new position using fixing screws or the locking tool set (Fig. 9) with padlock.
- The radiation beam outlet is now open.

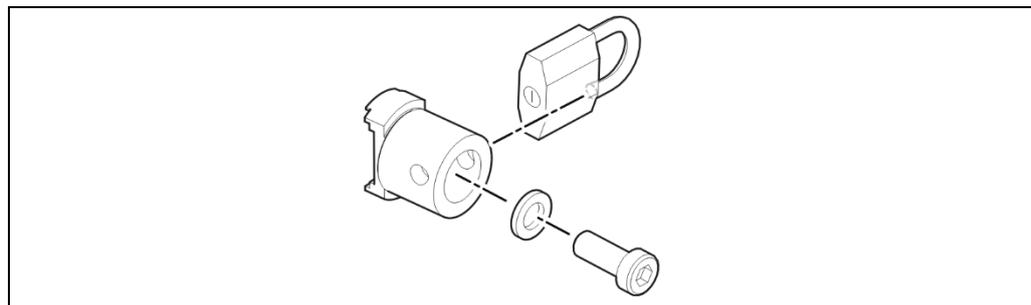


Fig. 9 Locking tool set (Mat. No. 73883)

Closing the locking mechanism

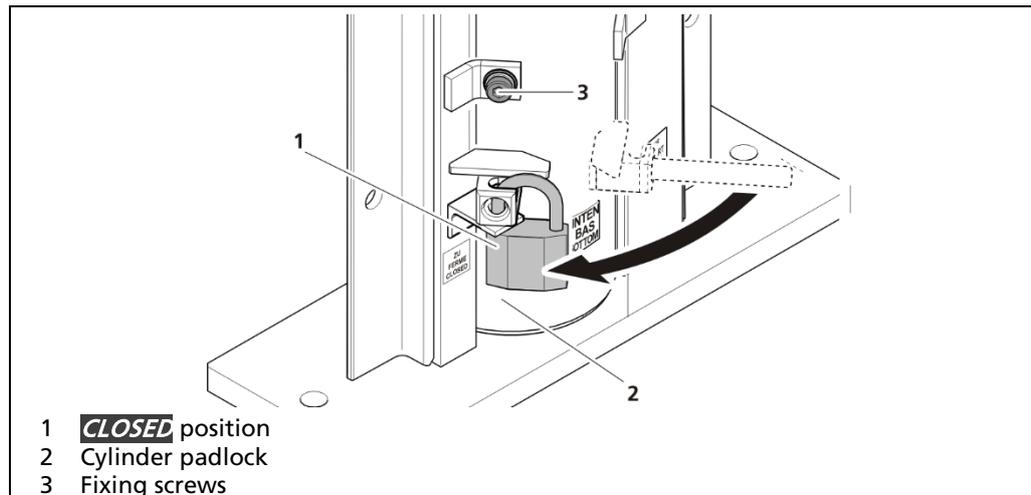


Fig. 10 Shield closed (**CLOSED** position)

1. Remove the fixing screws or the locking tool (Fig. 9).
 2. Turn the rotatable cylinder to the **CLOSED** position using the lever rod.
 3. Secure the rotatable cylinder in the new position using fixing screws and the cylinder padlock.
- ▶ The radiation beam outlet channel is now closed.

7 Maintenance and repair

The maintenance may only be carried out by competent people (see 2.3 Qualification of the personnel), which is supervised by experts or authorized persons. For repair work the radiation protection officer is to be consulted.

7.1 Safety instructions

⚠ CAUTION!



Danger caused by nuclear radiation!

Shields usually contain radioactive sources. An increased exposure to radiation may lead to damage to health.

- ▶ Consult the radiation safety officer responsible for your company.
- ▶ Ensure that no one is in the radiation beam during operation.
- ▶ Longer repair and maintenance work on shields must not be carried out with the source installed. If a removal of the source is not possible, please contact the manufacturer.

IMPORTANT



The applicable national regulations of the country of use have to be observed. Document the results of the tests you conducted using the check list in the appendix. This documentation as well as the seal test certificates must be kept during the entire life cycle of the shield.

NOTICE



Adjust the intervals of the visual inspection and the functional check to the ambient conditions. If the ambient conditions are especially rough, the atmosphere is corrosive and / or there is a serious threat of contamination, the intervals should be shortened accordingly.

NOTICE



If damage affecting the secure function of the shield is detected during handling, maintenance or regular checks, the damage must be repaired after consultation with Berthold before the system is used again. Only spare parts specified by Berthold may be used and required repair work may be carried out exclusively by authorized persons.

7.2 Visual inspection

The visual inspection must be carried out by people who are at least employees with general knowledge at the following intervals:

- before initial commissioning
- with every repair that may be required
- regularly at least every six months
- before shipping

When determining the intervals for the visual inspection, the following conditions are to be considered:

- ambient conditions (outdoors, rain, sunlight, wind)
- operating conditions (degree of utilisation of the plant, misuse)

Performing the Visual Inspection:

1. Check the shield for obvious damage (dents, cracks, holes etc.) and corrosion.
2. Only before shipping: Check if the shield is in closed position and if the cylinder padlock as well as the fixing screws are installed.

Tip



If there is any doubt regarding the actual position of the beam path: Check the dose rate at the shield using a dose rate measuring device. If the shield is closed, the dose rate in the direction of the beam may not be significantly higher than at other parts of the shield.

3. If any defects are identified during visual inspection, inform the radiation safety officer who will initiate the measures required for repairing the defects.

For further information, please contact Berthold.

7.3 Leak test

Depending on the supervisory authority responsible for the area where the source is used, regular leak tests must be carried out according to ISO 9978. These tests are to be carried out at the discretion of the competent supervisory authority either by an authorized expert or by the manufacturer. For this test, the corresponding source documentation must be made available.

Required documents

- Inventory list of the sources to be tested including indication of previous leak tests.
- **Source certificate containing the following information:** Nuclide, activity, procurement date, physical/chemical form, description of the enclosure and type of sealing, resistance against mechanical and thermal influences or classification of the source design. The source certificate is provided together with the source.
- Information on the location, application as well as on the maximum customary mechanical and thermal strains.
- If the source is installed in a device, a drawing is to be provided. In the drawing, the position of the source and of all parts intended to protect the source against external influences must be clearly indicated. Recommendations regarding the most appropriate test method should be given, e.g. by indicating alternative test areas. If required, recommendations should be given on how the necessary test can be conducted without affecting the functionality of the system or device.

Nearest accessible surface

The nearest accessible surface is defined as the area, which is most likely to be contaminated if a source is leaking. This area should be used primarily for performing the leak test.

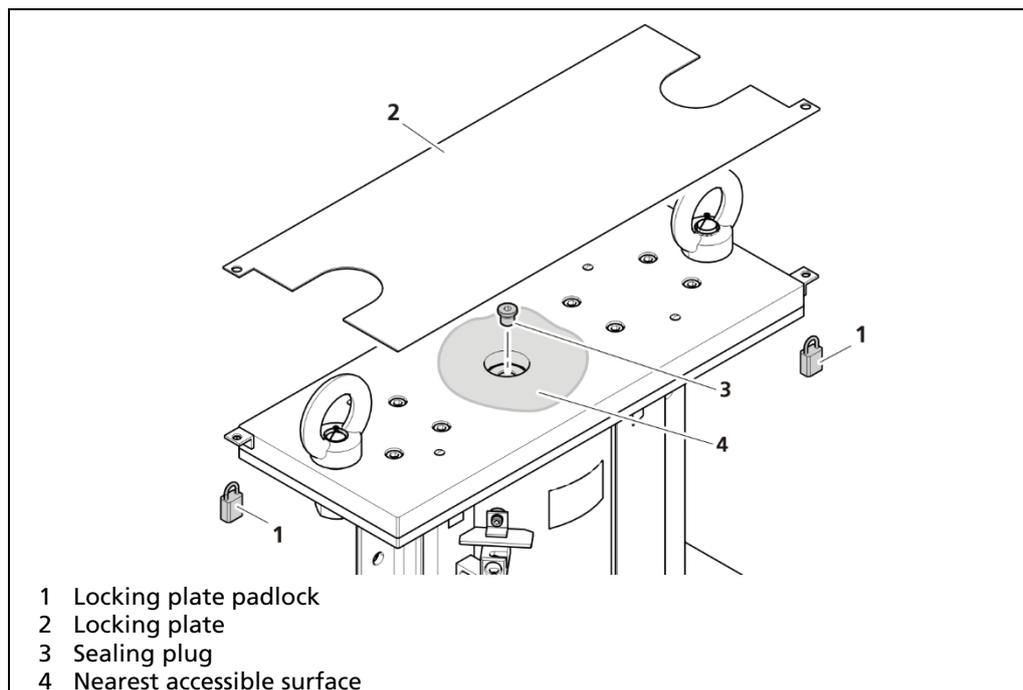


Fig. 11 Nearest accessible surface for leak test

The following instruction should be used:

- the area marked in grey in Fig. 11.
1. Remove the locking plate padlock and the locking plate
 2. Unscrew the sealing plug.
 3. Swab the area underneath the sealing plug and the surrounding area highlighted in Fig. 11.

7.4 Shutter check procedure

The functional test of the rotating mechanism must be carried out by a person who is at least an employee with general knowledge. We recommend conducting the test of the rotating mechanism at the same intervals as the visual inspection if the operating condition of the system allows.

1. Ensure that no employees are present in the direction of the radiation beam outlet. Otherwise, these persons will be exposed to the radiation beam when the locking mechanism is opened.
2. Open and close the shield several times (see chapter 6) while observing the measuring signal of the corresponding measurement system. The shield works properly if the measured value approximately corresponds to the maximum value when the locking mechanism is closed (**CLOSED** position).

Tip



Alternatively, you can measure the dose rate in the beam path. In this case, the dose rate in the **CLOSED** position must be clearly lower than in the **OPEN** position. At the same time, the dose rate in the **CLOSED** position may not be significantly higher than at other parts of the shield.

3. Return the rotatable cylinder to its starting position (**OPEN** position or **CLOSED** position).
 4. If applicable, reinstall the fixing screws and/or the cylinder padlock.
 5. Immediately report any malfunction or stiffness of the rotating mechanism to the competent radiation safety officer.
- ▶ The shutter check is completed.

7.5 Removing and installing the source

Before conducting extended maintenance work at the shield, you have to remove the source. After the maintenance work is finished, you can reinstall the source. This does require the direct handling of the unshielded source.

You must clarify with the responsible supervisory authority in advance whether your license for handling radioactive substances includes handling unshielded sources.

In any case, the source may only be installed/removed by authorised people. The radiation safety officer is responsible for planning and monitoring the work.

NOTICE



In order to keep the exposure to radiation during the removal / installation of the source as low as possible, all people involved should familiarise themselves with the exact procedure in advance.

NOTICE



Only remove / install the source in areas with closed floors. Any apertures which are larger than the external diameter of the source (7 mm) are to be covered before commencing any work.

7.5.1 Removing the source

To install the source, the following tools are required:

- pliers to securely grip the source
- 12 mm hexagon wrench to unscrew the sealing plug of the tungsten spacer
- M3 threaded rod (approx. 50 mm) for screwing into tungsten spacer and pulling out the attached rod source

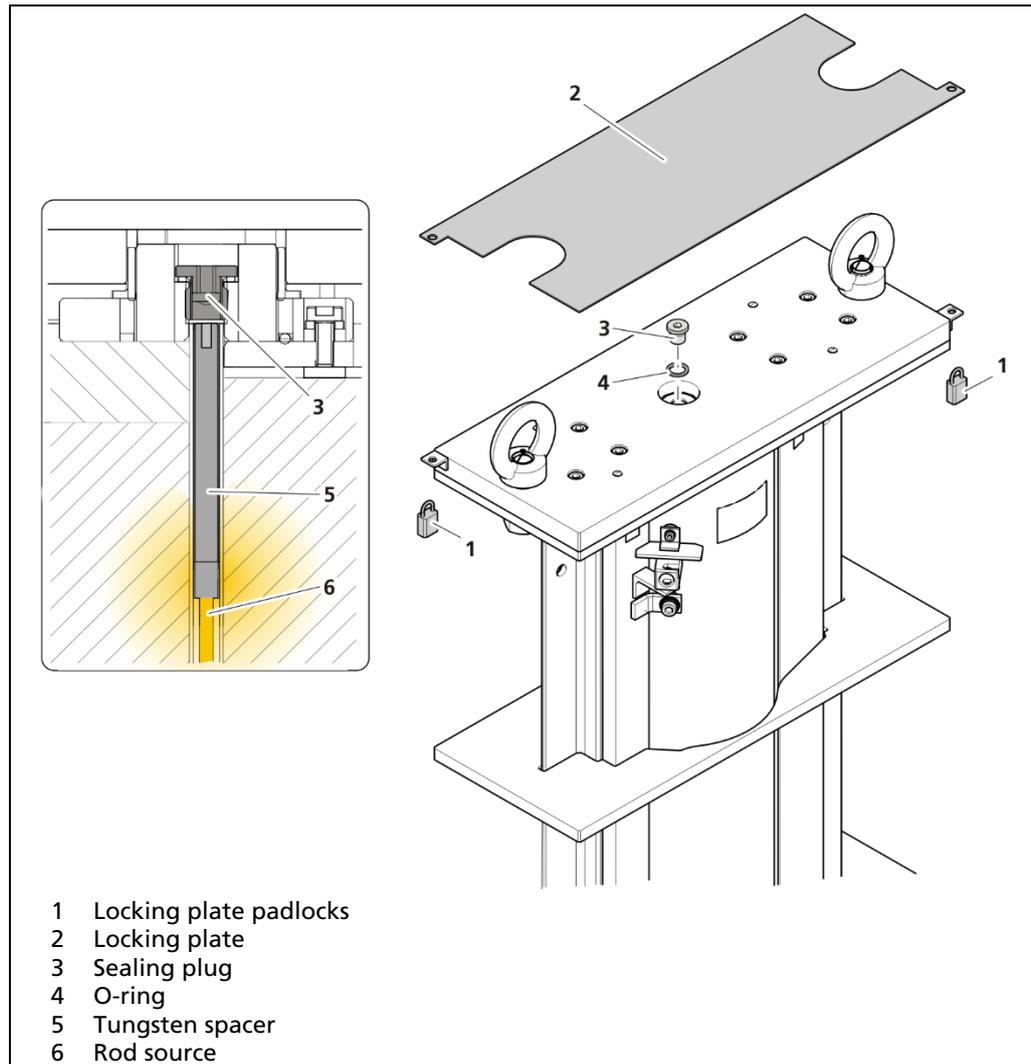


Fig. 12 Make the source accessible

To remove the source, proceed as follows:

1. Prepare the transport shield (Fig. 13, item 4) for the source, remove the cover (Fig. 13, item 6) and the locking plug (Fig. 13, item 5).
2. Remove the locking plate padlocks (Fig. 12, item 1) and remove the locking plate (Fig. 12, item 2).
 - ▶ You can now see the sealing plug (Fig. 12, item 3).
3. Unscrew the screwed sealing plug (Fig. 12, item 3).
4. Remove the O-ring (Fig. 12, item 4).

⚠ CAUTION!



Danger caused by ionising radiation!

When the source is removed, the unshielded source must be handled as quickly as possible. An increased exposure to radiation may lead to damage to health.

- ▶ Hold the source only with pliers.
- ▶ Keep the source far from the body.
- ▶ Ensure that no people are in the field of radiation during the work.

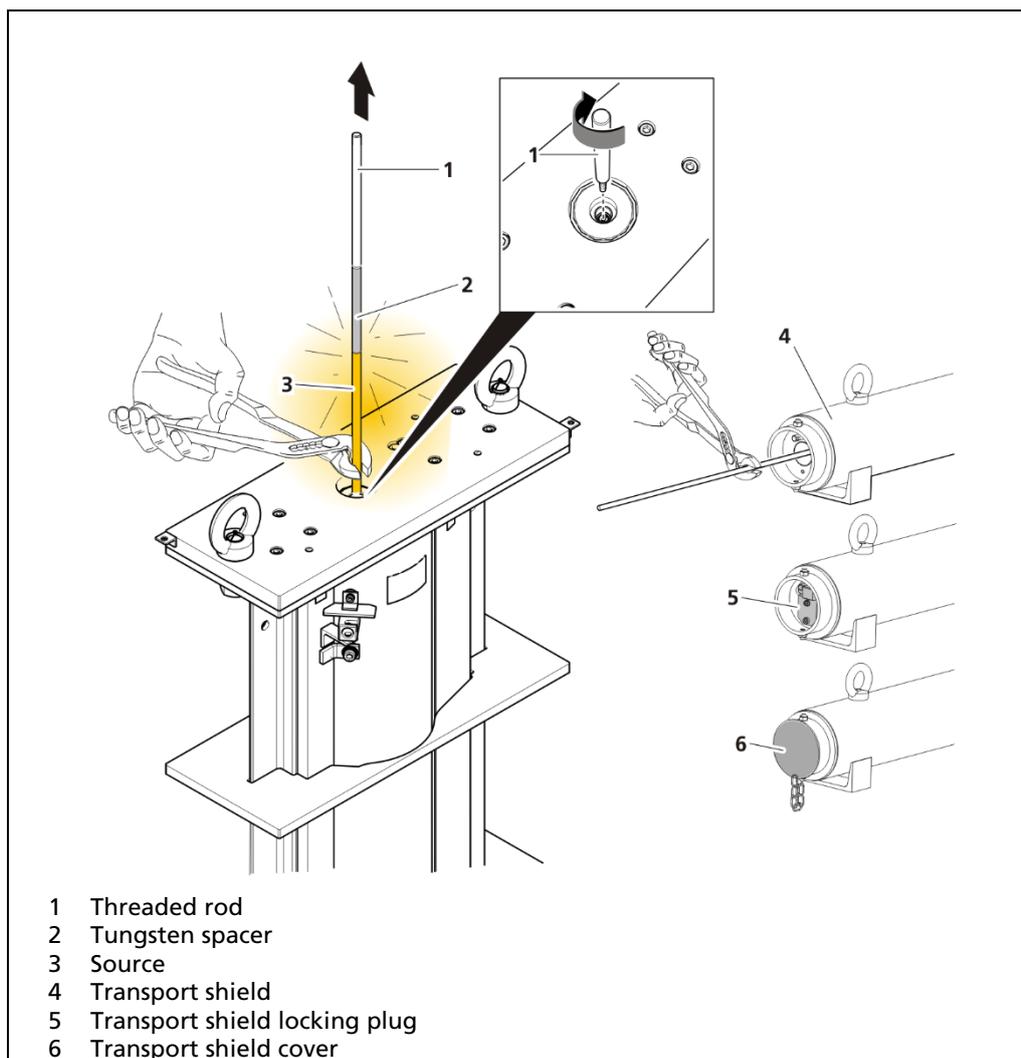


Fig. 13 Remove source out of shield

5. Screw the threaded rod (Fig. 13, item 1) into the thread of the tungsten spacer (Fig. 13, item 2).
6. Pull the source out of the shield with the threaded rod. Hold the source (Fig. 13, item 3) with pliers.
7. Insert the source as quickly as possible into the provided and opened transport shield (Fig. 13, item 4) and close the locking plug (Fig. 13, item 5) and the cover (Fig. 13, item 6).

7.5.2 Installing the source

To install the source, the following tools are required:

- pliers to securely grip the source
- 12 mm hexagon wrench to unscrew the sealing plug of the distance piece

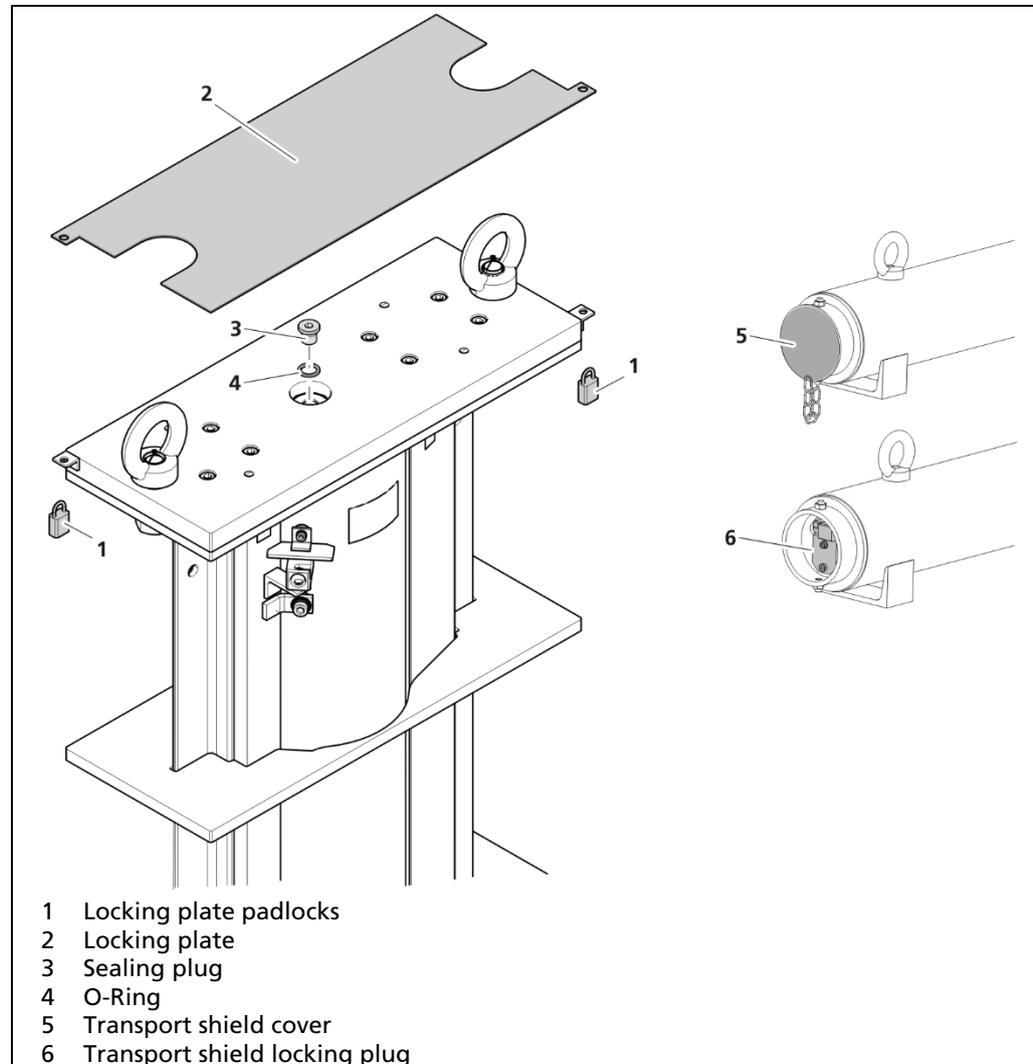


Fig. 14 Prepare the installation

To install the source, proceed as follows:

1. If necessary, remove the locking plate padlocks (Fig. 14, item 1) the locking plate (Fig. 14, item 2).
2. Unscrew the sealing plug (Fig. 14, item 3).
3. Remove the O-ring (Fig. 14, item 4).
4. Prepare the transport shield (Fig. 13, item.4) with the source to be installed and remove the cover of the transport shield (Fig. 14, item 6).
5. Remove the locking plug (Fig. 14, item 7) to open the transport shield.

⚠ CAUTION**Danger caused by ionising radiation!**

When the source is removed, the unshielded source must be handled as quickly and little as possible. An increased exposure to radiation may lead to damage to health.

- ▶ Hold the source only with pliers.
- ▶ Keep the source far from the body.
- ▶ Ensure that no people are in the field of radiation during the work.

6. Grasp the tungsten spacer (Fig. 15, item 3) and pull the rod source out of the transport shield with the pliers (Fig. 15, item 5).

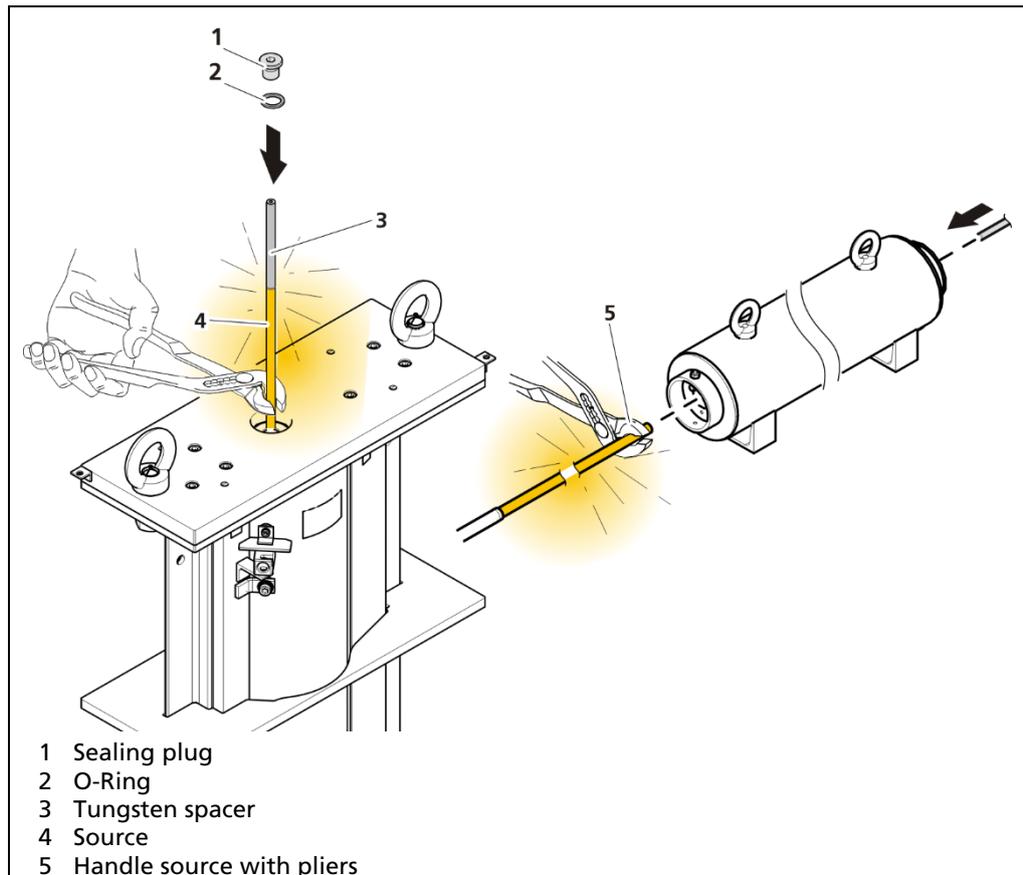


Fig. 15 Installing the source in the shield

7. Insert the source as quickly as possible into the shield.
8. Attach the O-ring (Fig. 15, item 2) and screw the sealing plug (Fig. 15, item 1).
9. Place the locking plate (Fig. 14, Pos.2) and secure locking plate with locking plate padlocks (Fig. 14, item 1).
10. Turn the rotatable cylinder to the starting position (position **OPEN** or position **CLOSED**).
11. Replace the fixing screws and the cylinder padlock.
 - ▶ The installation is complete.

7.6 Source replacement

When to replace the source

Generally, the radioactive source used enables a service life between 5 and 10 years. A replacement of the source is only required if the statistical fluctuations of the output signal are inadmissibly large and compensation by increasing the time constant is no longer possible, e.g. for control-technical reasons.

Manufacturer's number for reordering

When reordering, always indicate the manufacturer's number of the source used since the new source must correspond to the original version.

The manufacturer's number can be found on the type plate of the shield as well as on the individual seal test certificate of each source.

7.6.1 Replacing the source

Replacing the source does require the direct handling of the unshielded source.

You must clarify with the responsible supervisory authority in advance whether your license for handling radioactive substances covers handling unshielded sources.

In any case, the source may only be installed/removed by authorised people. The radiation safety officer is responsible for planning and monitoring the work.

NOTICE



In order to keep the exposure to radiation during the removal / installation of the source as low as possible, all persons involved should familiarise themselves with the exact procedure in advance.

NOTICE



Only remove / install the source in areas with closed floors. Any apertures which are larger than the external diameter of the source (7 mm) are to be covered before commencing any work.

To replace the source, the following tools are required:

- pliers to securely grip the source
- 12 mm hexagon wrench to unscrew the screwed sealing plug of the tungsten spacer
- M3 threaded rod (approx. 50 mm)

To replace the source, proceed as follows:

1. Prepare the transport shield and open it.
2. Remove the old source from the shield and insert the old source in the transport shield, see chapter 7.5.1.
3. Check the shield for damage and wear and repair it if required (see chapter 7.2).
4. Remove the type plate of the old source from the shield and rivet on the new type plate.
5. Install the new source in the shield, see chapter 7.5.2.

8 Uninstallation

The shield may only be dismantled and put out of operation by competent people (see 2.3 Qualification of the personnel). Possibly the radiation protection officer is to be consulted, as well.

8.1 Safety instructions

WARNING!



Danger of injury by falling loads!

- ▶ Never stand underneath a lifted or suspended load, keep at a safe distance.
- ▶ Only use tested sling gear components appropriate for the transport weight.
- ▶ Observe the marking for the centre of gravity on the outer packaging, if applicable.
- ▶ Wear head protection and safety shoes.



WARNING!



Danger of injury caused by heavy and bulky system components!

- ▶ Heavy and bulky system components should only be handled using aids and by a at least 2 people.
- ▶ Observe the guidelines for safe handling of heavy loads.
- ▶ Ensure stability and use the provided fixing options.

CAUTION!



Danger caused by ionising radiation!

Shields usually contain radioactive sources. An increased exposure to radiation may lead to damage to health.

- ▶ Consult the radiation safety officer responsible for your company.
- ▶ Transport the source exclusively inside a closed and secured shield.

IMPORTANT



The applicable national regulations of the country of use have to be obeyed.

8.2 Dismounting the shield

⚠ WARNING!



Danger of injury by falling loads!

Falling loads can endanger people.

- ▶ Never stand underneath a lifted or suspended load, keep at a safe distance.
- ▶ Cordon off the danger zone with barrier tape.

⚠ WARNING!



Danger of crushing by moving the shields!

Due to the high weight of the shield, there is a risk of injury.

- ▶ Heavy and bulky system components should only be handled using aids and by a at least 2 people.
- ▶ Do not reach between the shield and the container or mounting base.

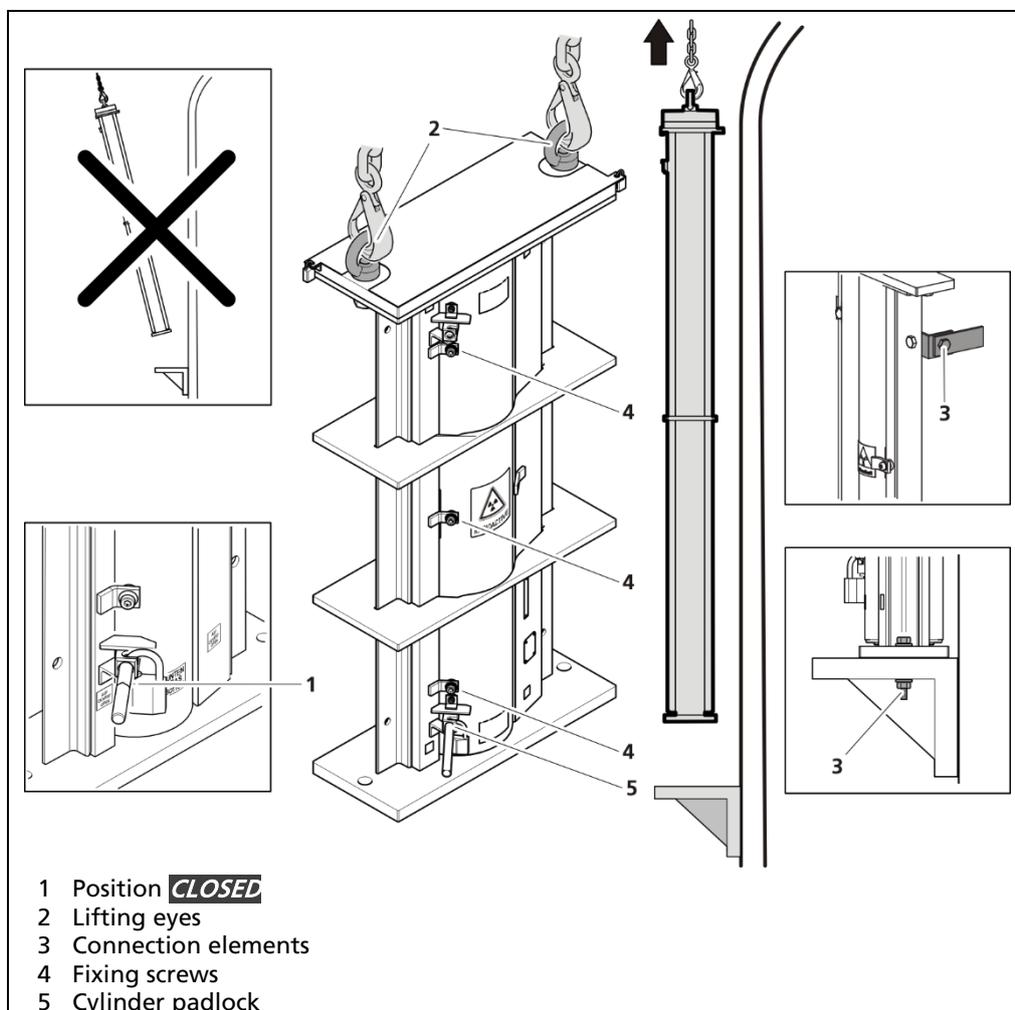


Fig. 16 Dismantle the shield

1. Ensure that the shield has been closed and secured.
 - ▶ The rotatable cylinder is in to the **CLOSED** position
 - ▶ The fixing screws (Fig. 16, item 4) and, if applicable, the cylinder padlock (Fig. 16, item 5) are installed.

2. Fasten the shield with sufficiently sized lifting straps at the lifting eyes (Fig. 16, item 2).
3. Remove all connection elements (Fig. 16, item 3).
4. Raise and set down the shield vertically.

8.3 Disposal of radioactive substances

NOTICE



Should you have any questions regarding the transportation or disposal of radioactive substances, please contact the manufacturer.

Generally, each country has a depot where radioactive material is accepted and can be disposed of.

8.4 Shipment of radioactive substances

If you wish to ship radioactive material, you must observe the international regulations for the transport of dangerous goods and for radiation protection as well as any applicable national regulations. It is the full responsibility of the sender to comply with these regulations.

Please also note the following:

- Dose rate at the surface of the packaging must be lower than 2000 $\mu\text{Sv/h}$.
- Dose rate at a distance of 1m from the surface of the packaging must be lower than 100 $\mu\text{Sv/h}$.
- The packaging must be marked with a reference for dangerous goods in accordance with the applicable regulations with the UN number.
- Furthermore, transport documents with the correct description of the content as well as an accident procedures sheet according to the ADR regulations are required.
 - With regard to transport by air, the IATA DGR.
 - IAEA as the basis for all regulations.
- Packaging must comply with the current ADR regulations (for example European Agreement concerning the International Carriage of Dangerous Goods by Road).
- The cargo must be secured in the transport vehicle according to the applicable national and international regulations.
- Before dispatch, each shield containing a source must be subjected to a visual inspection by the user (chapter 7.2). The shields may only be dispatched if the test requirements are met.
- At the time of shipment, a valid leak test lab result for the source must be available.

Preconditions for returning shields to BERTHOLD TECHNOLOGIES GmbH & Co. KG

- Radioactive substances and their shields must not be damaged in any way and a respective, valid leak test certificate must be provided. The leak test certificate issued before the arrival of the radioactive substances in Germany may not be older than 6 months.
- If radioactive sources with isotope Am-241 or Cm-244 are returned, the special form certificate must be attached.
- At all times, any radioactive material sent to us must be sufficiently labelled with your name and address. If we have sent you a quotation in advance, our quotation number must be indicated, as well.
- Radioactive substances may only be returned after the respective approval by Berthold. We are pleased to send you a quotation regarding the costs.
- Radioactive substances must be sent to Bad Wildbad, Germany, carriage paid. Berthold will not assume any costs for customs clearance or transportation.
- Berthold is to be informed about any return transport in advance. Berthold will reject any radioactive substances sent to Berthold without prior notice. Any storage costs accrued in such case shall be borne by the sender.
- A Before shipment of a radioactive source, a notification form with information about the source needs to be filled out and sent to Berthold. We will be pleased to send you the current version of the notification form.
- It is mandatory to attach a copy of the enclosed notification form and the valid leak test certificate to the shield. The original documents must be enclosed with the shipping documents. Before shipment, the documents must be sent to Berthold via e-mail.

NOTICE



Observe the applicable regulations when taking them out of operation and disposal.

9 Appendices

9.1 Check lists for tests

Visual inspection

Requirement	met	not met
The shield must be free from damage.		
The type plates must be legible.		
If there is corrosion, it must not affect the function and stability of the shield.		
Only before shipment: The beam path must be closed.		
Only before shipment: The locking plate must be mounted and secured.		
Only before shipment: The transport securing device(s) must be installed.		
Name of inspector:	Date:	Signature:

Checking the functions of the locking mechanism

Requirement	met	not met
It must be possible to move the locking mechanism completely back and forth between the two positions (OPEN / CLOSED).		
The measuring signal of the measurement system must approximately correspond to the maximum value in the CLOSED position.		
If alternatively a dose rate measurement device is used, the dose rate in the CLOSED position must be clearly lower than in the OPEN position.		
Only with manual actuation: It must be possible to turn the rotatable cylinder using one hand and the lever rod.		
Only with pneumatic actuation: The air pressure required for moving the locking mechanism may not exceed the maximum permissible air pressure for the pneumatic actuator (see technical data of the pneumatic actuators)		
Name of inspector:	Date:	Signature:

Modifications due to technical advancement reserved.

BERTHOLD TECHNOLOGIES GmbH & Co. KG

ROD SOURCE SHIELDS LB 8300

Technical Information



Table of contents

1.	Maximum permissible activities during transport	3
2.	Ambient conditions during mounting	3
3.	Technical Data	4
3.1.	Variant 100.....	4
3.2.	Variant 150.....	5
3.3.	Variant 200.....	6
3.4.	Variant 270.....	7
3.5.	Dimensions Variant 100 / 150	8
3.6.	Dimensions variant 200 / 270	9
3.7.	Mounting base for LB 8300	10
3.8.	Mounting Kit Adapter Plate	11
4.	Rod Source Cascade	12
4.1.	Mounting Variant Stacked Rod Source Cascaded	13
4.2.	Mounting variants staggered rod source cascade	14

1. Maximum permissible activities during transport

The activities indicated here are the maximum activities with which the individual shields can be loaded according to the international dose rate limit for the transport of radioactive substances. National dose rate limits for the operation of measurement systems are usually significantly lower and only allow loading with significantly lower activities. The national provisions must be complied with.

Shield model	Maximum activity (Co-60)		Maximum activity (Cs-137)	
	MBq/mm	mCi/mm	MBq/mm	mCi/mm
Variant 100	1.38	max. 0.037	49.71	max. 1.34
Variant 150	11.94	max. 0.32	1657.29	44.79
Variant 200	67.44	max. 1.82	43166	1166
Variant 270	475.60	max. 12.85	43166	1166
	calculated values			

2. Ambient conditions during mounting

The shield is lifted and positioned using the two lifting eyes. The minimum permissible temperature during assembly is -20 °C. Follow the instructions in the operating manual.

3. Technical data

3.1. Variant 100

Variant	Housing	Shielding	Number of support plates	Source length	Weight	Operating temperature	Coating	Fire proof
ID: 69540-01	C-Stahl S235	43 mm lead (rotatable cylinder), 40 mm (lead shielding profile)	-	500 mm	97 kg	-40°...+200°C	RAL 1004 2K polyurethane varnish	NO
ID: 69540-02			-	600 mm	110 kg			
ID: 69540-03			-	700 mm	123 kg			
ID: 69540-04			-	800 mm	136 kg			
ID: 69540-05			1	900 mm	150 kg			
ID: 69540-06			1	1000 mm	163 kg			
ID: 69540-07			1	1100 mm	176 kg			
ID: 69540-08			1	1200 mm	189 kg			
ID: 69540-09			1	1240 mm	193 kg			

3.2. Variant 150

Variant	Housing	Shielding	Number of support plates	Source length	Weight	Operating temperature	Coating	Fire proof
ID: 69550-01	Carbon steel S235	70 mm lead (rotatable cylinder), 63 mm lead (shielding profile)	-	500 mm	249 kg	-40° ... +200°C	RAL 1004 2K polyurethane varnish	NO
ID: 69550-02			-	600 mm	274 kg			
ID: 69550-03			-	700 mm	303 kg			
ID: 69550-04			-	800 mm	338 kg			
ID: 69550-05			1	900 mm	368 kg			
ID: 69550-06			1	1000 mm	397 kg			
ID: 69550-07			1	1100 mm	426 kg			
ID: 69550-08			1	1200 mm	456 kg			
ID: 69550-09			1	1240 mm	466 kg			

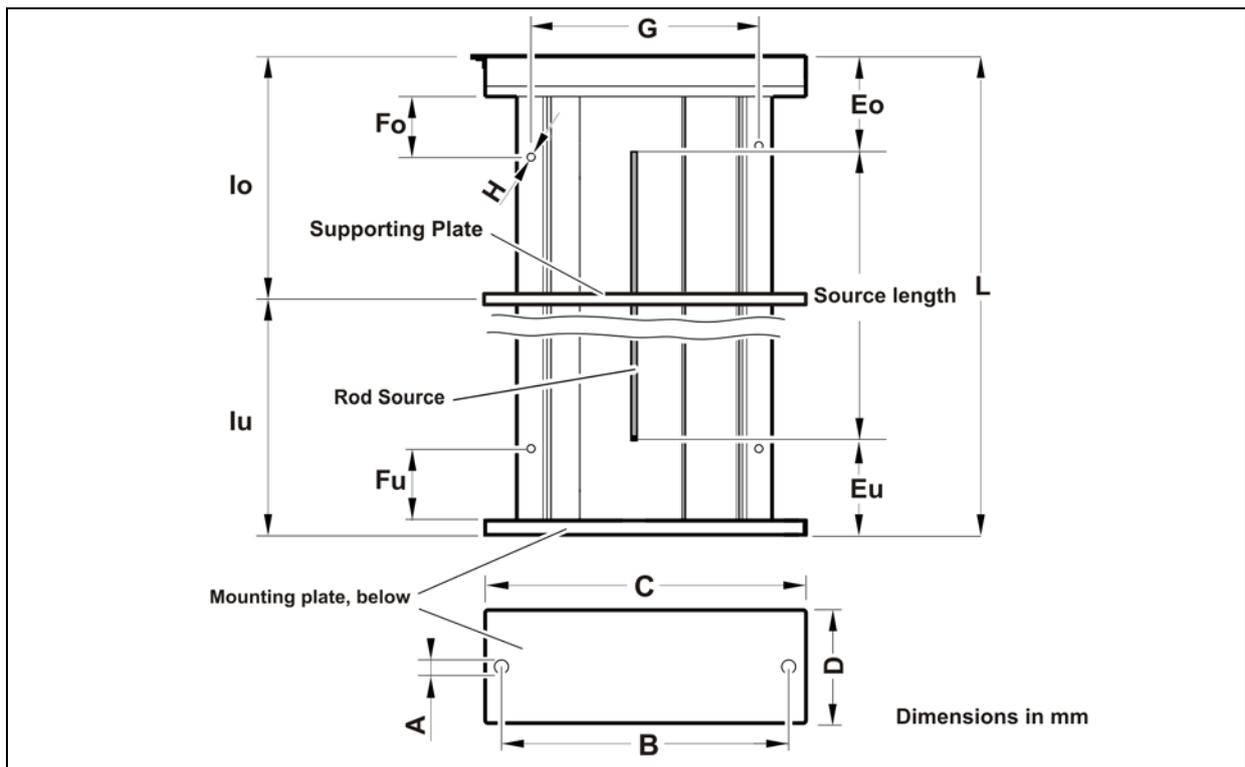
3.3. Variant 200

Variant	Housing	Shielding	Number of support plates	Source length	Weight	Operating temperature	Coating	Fire proof
ID: 69560-01	Carbon steel S235	88 mm lead (rotatable cylinder), 73 mm (lead shielding profile)	-	500 mm	412 kg	-40° ... +200°C	RAL 1004 2K polyurethane varnish	NO
ID: 69560-02			-	600 mm	456 kg			
ID: 69560-03			-	700 mm	500 kg			
ID: 69560-04			-	800 mm	563 kg			
ID: 69560-05			2	900 mm	607 kg			
ID: 69560-06			2	1000 mm	652 kg			
ID: 69560-07			2	1100 mm	697 kg			
ID: 69560-08			2	1200 mm	741 kg			
ID: 69560-09			2	1240 mm	757 kg			

3.4. Variant 270

Variant	Housing	Shielding	Number of support plates	Source length	Weight	Operating temperature	Coating	Fire proof
ID: 69570-01	Carbon steel S235	121 mm lead (rotatable cylinder), 100 mm lead (shielding profile)	-	500 mm	761 kg	-40° ...+200°C	RAL 1004 2K polyurethane varnish	NO
ID: 69570-02			-	600 mm	838 kg			
ID: 69570-03			2	700 mm	938 kg			
ID: 69570-04			2	800 mm	1016 kg			
ID: 69570-05			2	900 mm	1094 kg			
ID: 69570-06			2	1000 mm	1171 kg			
ID: 69570-07			2	1100 mm	1249 kg			
ID: 69570-08			2	1200 mm	1327 kg			
ID: 69570-09			2	1240 mm	1355 kg			

3.5. Dimensions variant 100 / 150

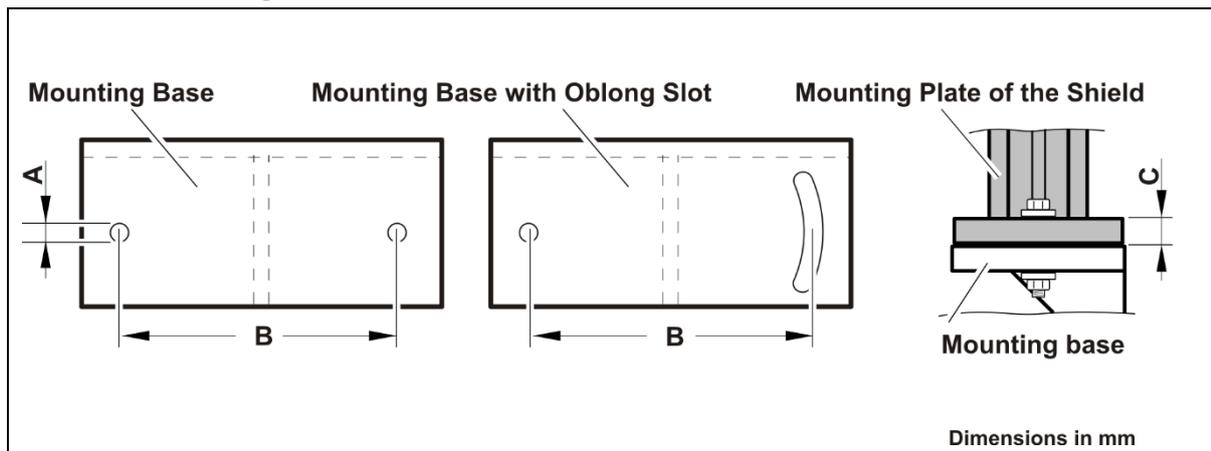


TI-Abb. 1 Dimension Variant 100 / 150

Variant	ID	Source length	L	A	B	C	D	Eu	Eo	Fu	Fo	G	H	io*	lu*
100	69540-01	500	650											-	-
	69540-02	600	750											-	-
	69540-03	700	850											-	-
	69540-04	800	950											-	-
	69540-05	900	1050	Ø14	305	340	120	77	73	120	132	225	Ø12	531	519
	69540-06	1000	1150											581	569
	69540-07	1100	1250											631	619
	69540-08	1200	1350											681	669
	69540-09	1240	1390											701	689
150	69550-01	500	700											-	-
	69550-02	600	800											-	-
	69550-03	700	900											-	-
	69550-04	800	1000											511	489
	69550-05	900	1100	Ø18	425	465	180	102	98	125	137	332	Ø14	561	539
	69550-06	1000	1200											611	589
	69550-07	1100	1300											661	639
	69550-08	1200	1400											711	689
	69550-09	1240	1490											731	709

* Dimensional deviation due to welded construction

3.7. Mounting base for LB 8300



TI-Abb. 3 Mounting base dimensions for LB 8300 Rod Source Shields

Variant	Mounting base	A	B	C
100	> 340 x 120	> Ø 14	305	15
150	> 465 x 180	> Ø 18	425	20
200	> 560 x 220	> Ø 26	510	25
270	> 600 x 300	> Ø 26	510	30

NOTICE



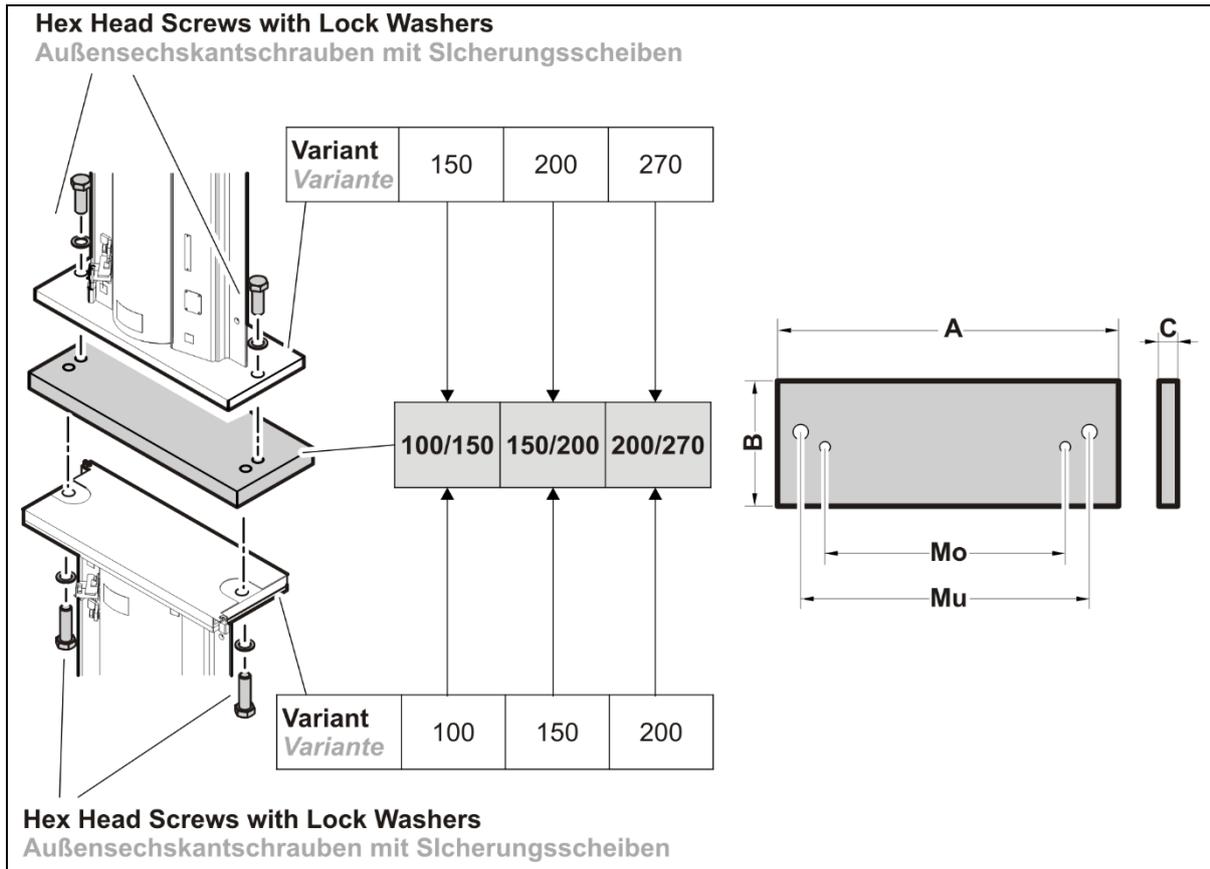
Mounting base installed by the installation contractor or end user. The base must be designed for the total load of the shield(s) and is designed for the total load of the shield(s) and be secured to the vessel or the supporting structure, e.g. welded.

IMPORTANT



The mounting base with the oblong slot allows the shield to be rotated to better aim the radiation beam at the detector.

3.8. Mounting kit adapter plate



TI-Abb. 4 Mounting Kit Adapter Plate

Ident. No.	Description	A	B	C	Mo	Mu	Connection Elements
73291	Mounting Kit Adapter Plate 100/150	465	120	20	305	425	M12x35, M16x45 Lock washer NL12, NL16
73295	Mounting Kit Adapter Plate 150/200	560	180	25	425	510	M24x60, M16x45 Lock washer NL24, NL16
73301	Mounting Kit Adapter Plate 200/270	600	220	40	510	510	M24x60 Lock washer NL24

NOTICE

- i** Only use the supplied screws and lock washers. Tightening torque recommendation:
- ▶ M12: 56 Nm
 - ▶ M16: 136 Nm
 - ▶ M24: 460 Nm

4. Rod source cascade

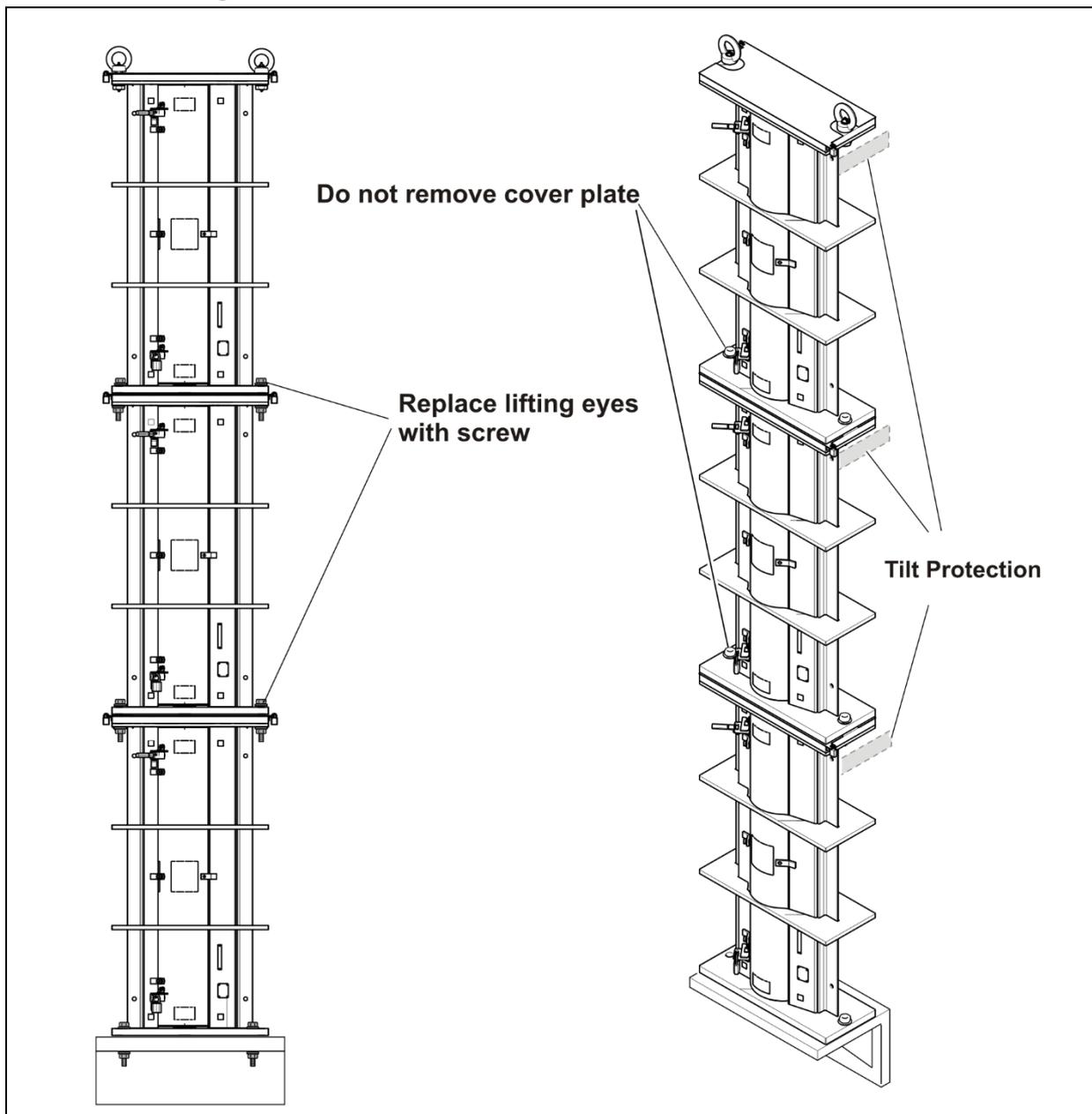
If several rod sources are needed, the individual shields are arranged directly on top of each other or staggered. Multi-part shields are marked by the letters A, B, C etc. from top to bottom. Each part must be secured adequately, e.g. by using a flat iron as anti-tipping device.

NOTICE



Following installation of the shields, the function of the positioning mechanism has to be tested. The shutter check has to be carried out every six months and recorded.

4.1. Mounting variant stacked rod source cascaded



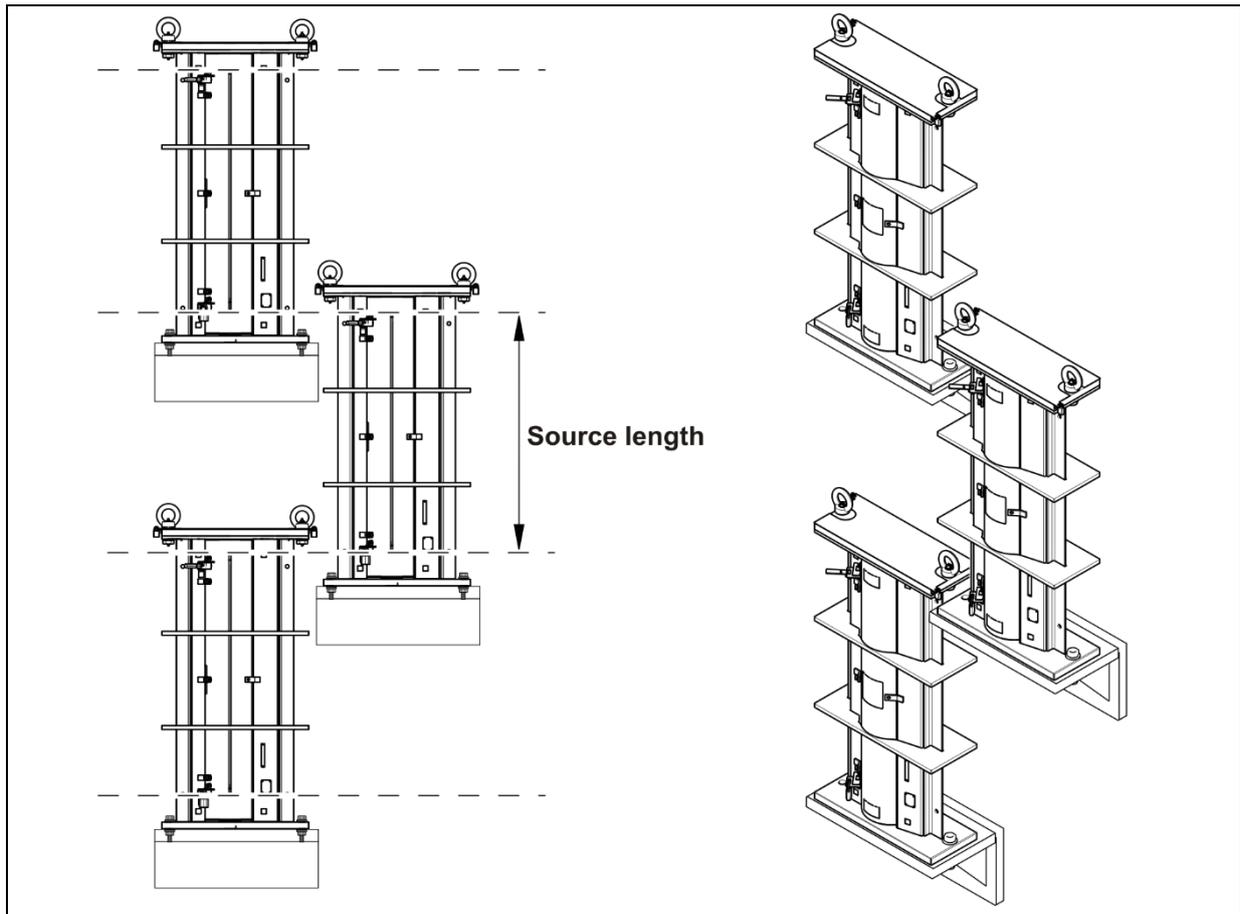
TI-Abb. 5 Mounting variant stacked (example illustration))

NOTICE



Tilt-protections should be installed by the installation contractor or end user, and affixed to the vessel or the supporting structure, e.g. welded. The anti-tipping measures must be installed on both sides and on each shield.

4.2. Mounting variants staggered rod source cascaded



TI-Abb. 6 Mounting variants staggered (example illustration)