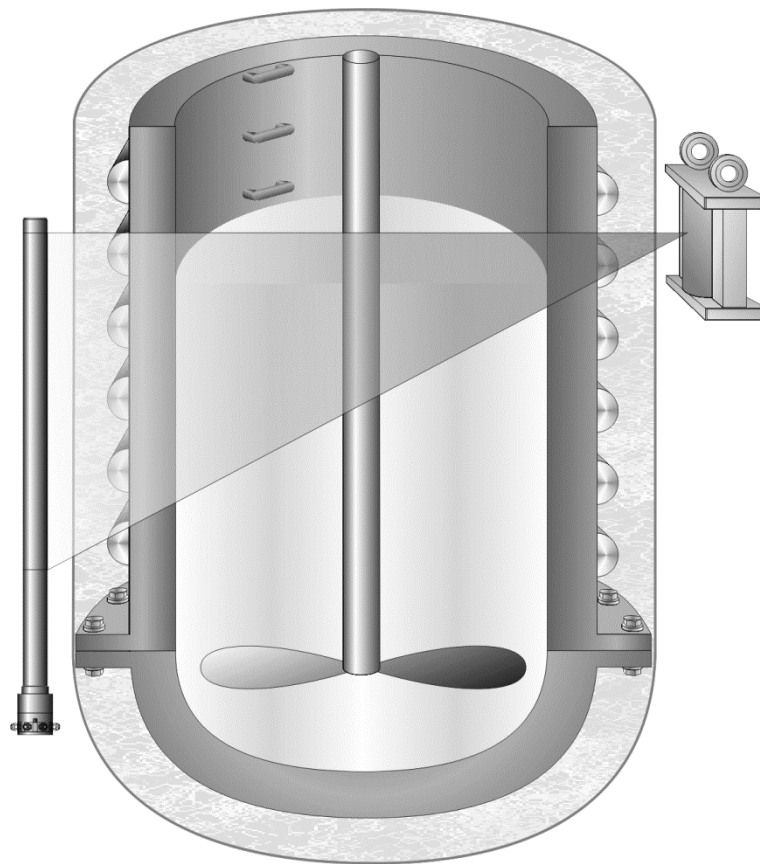


Technical Information

Level LB 480

Level Gauge
Füllstandmessung

Field mounted components
Messstellen-Komponenten

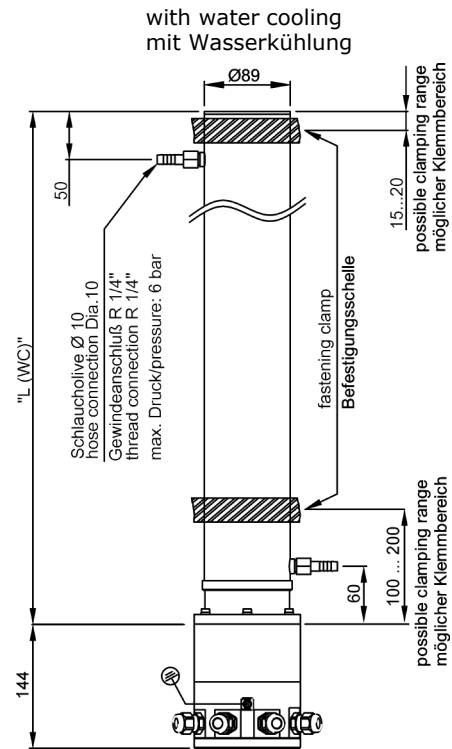
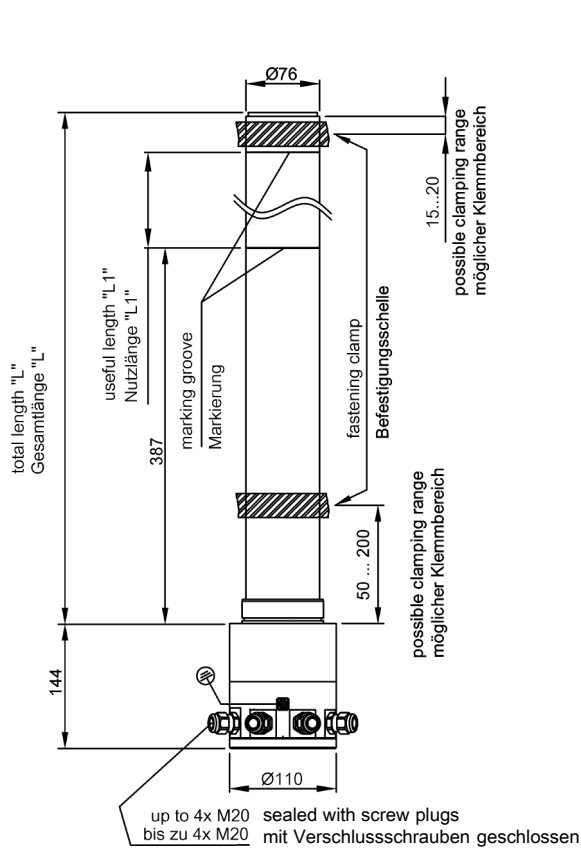


LB 480

Dimensions in mm
Abmessungen in mm

1. UniSENS (Version f. Zone 1/2)

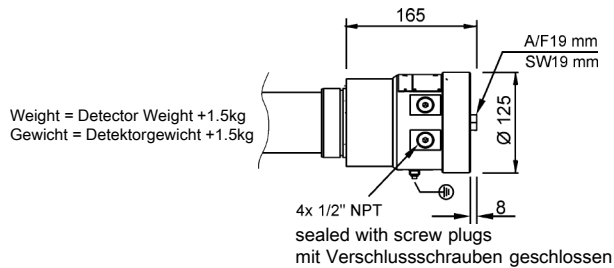
LB 480-2.-0.
LB 480-2.-1.
LB 480-2.-2.
LB 480-2.-Z.



All other dimensions see left hand.
Aller anderen Maße siehe links.

1.1 UniSENS (Version f. Divisions 1/2)

LB 480-2.-F.
LB 480-2.-G.



The detector version for divisions (NEC/CEC) differs only in the terminal housing, compared to the standard version illustrated above. The dimensions of this terminal housing are illustrated in this drawing.

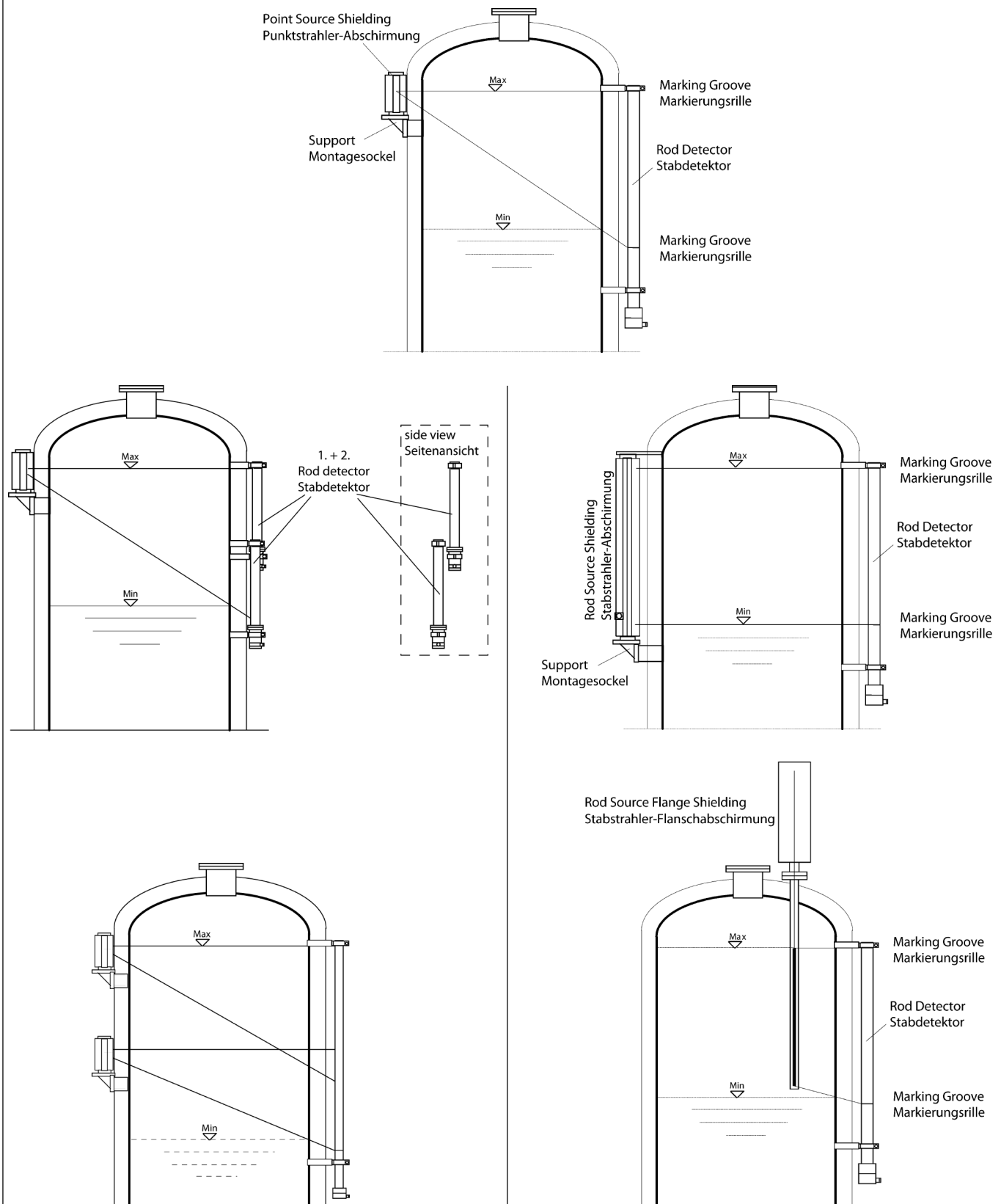
Die Detektor-Version mit Ex-Zulassung für Divisions (NEC/CEC) unterscheidet sich ausschließlich im Anschlusskopf, verglichen mit der oben aufgeführten Standardvariante. Die Abmessungen für den Anschlusskopf sind aus dieser Zeichnung zu entnehmen.

Type Typ	L	L (WC*)	L1	Weight (kg) Gewicht (kg)	Weight with WC* (kg) Gewicht mit WK* (kg)	Weight with WC* + Water (kg) Gewicht mit WK* + Wasser (kg)
LB 480-2A	929		500	16,0		
LB 480-2E	1429		1000	20,5		
LB 480-2I	1929		1500	25,0		
LB 480-2K	2429		2000	29,5		
LB 480-2B		933	500		25.0	25.5
LB 480-2F		1433	1000		32.0	33.0
LB 480-2J		1933	1500		39.5	40.0
LB 480-2L		2433	2000		48.0	50.0

* WC/WK = Water Cooling / Wasserkühlung

LB 480

1.2 Examples for Measurement Arrangements with Rod Detector Beispiele für Messanordnungen mit Stabdetektor

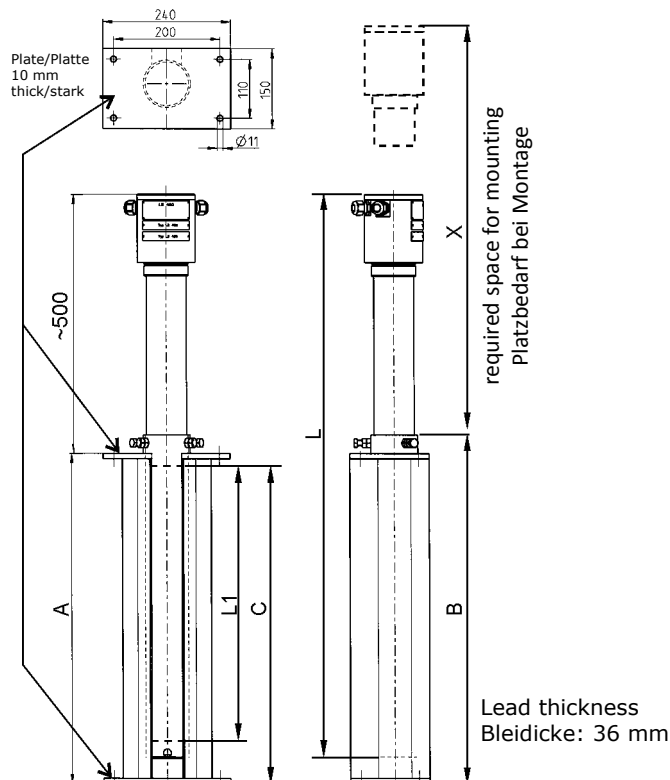


LB 480

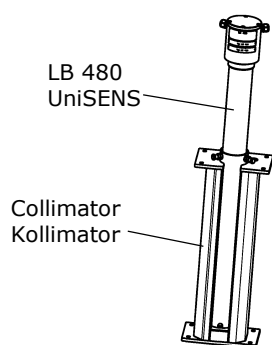
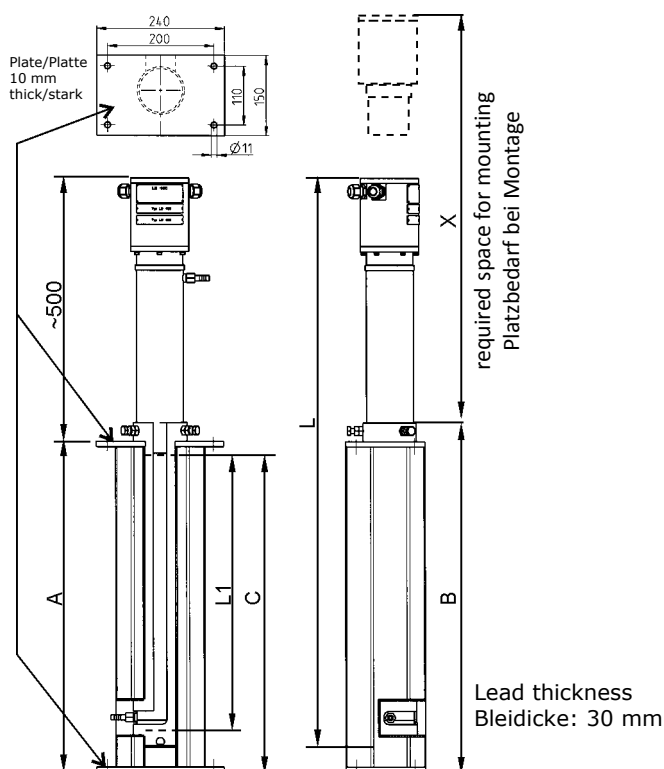
Dimensions in mm
Abmessungen in mm

1.3 Collimator for Rod Detector Kollimator für Stabdetektor

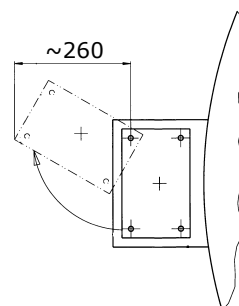
for Detectors without Water Cooling
für Detektoren ohne Wasserkühlung



for Detectors with Water Cooling (WC)
für Detektoren mit Wasserkühlung (WK)

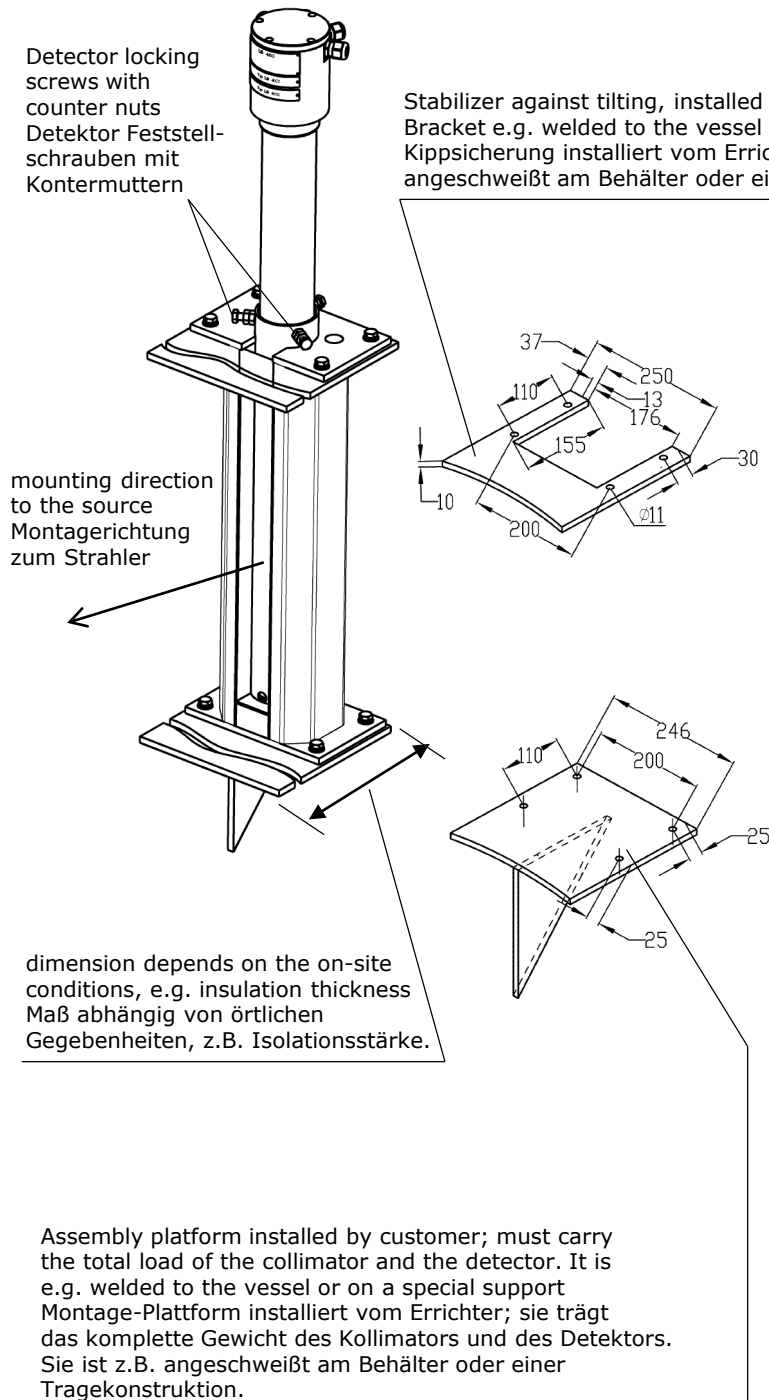


For installation/deinstallation, space for
swiveling the collimator is recommended
Empfohlener Installations-/Deinstallationsraum
zum Ausschwenken des Kollimators

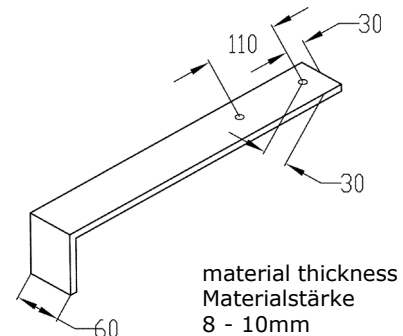


Part No. Id No.	for WC für WK	L1 sensitive length empfindliche Länge	A	B	L detector length Detektorlänge	C	X	Weight Gewicht (kg)
59957-050		500	620	655	1073	590	1100	110
59957-100	-	1000	1120	1155	1573	1090	1600	195
59957-150		1500	1620	1655	2073	1590	2100	280
59957-200		2000	2120	2155	2573	2090	2600	365
60085-050		500	620	655	1077	590	1100	100
60085-100	✓	1000	1120	1155	1577	1090	1600	180
60085-150		1500	1620	1655	2077	1590	2100	255
60085-200		2000	2120	2155	2577	2090	2600	330

1.4 Examples for Mounting Devices and the Installation at the Collimator Beispiele für die Halterung und Montage des Kollimators



Alternative to the stabilizer against tilting in the drawing on the left, you can use a one-sided mounted angle bracket as a stabilizer too.
Alternative zur in der neben stehend dargestellten Kippsicherung, kann auch ein einseitig anmontierter Winkel als Kippsicherung montiert werden.



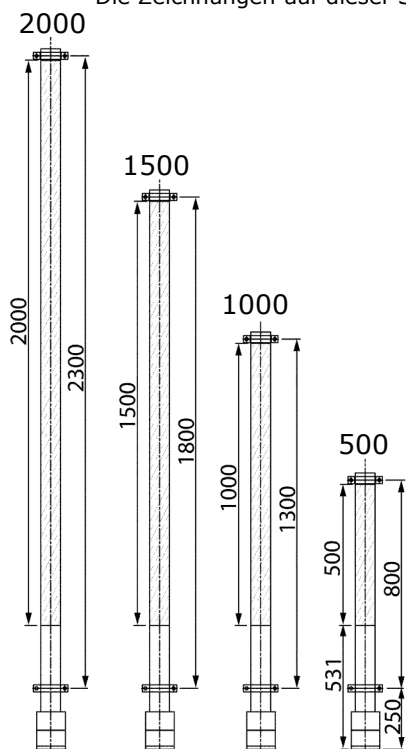
angle bracket as tilt protection, welded
Winkeleisen als Kippsicherung, angeschweißt

Vessel
Behälter

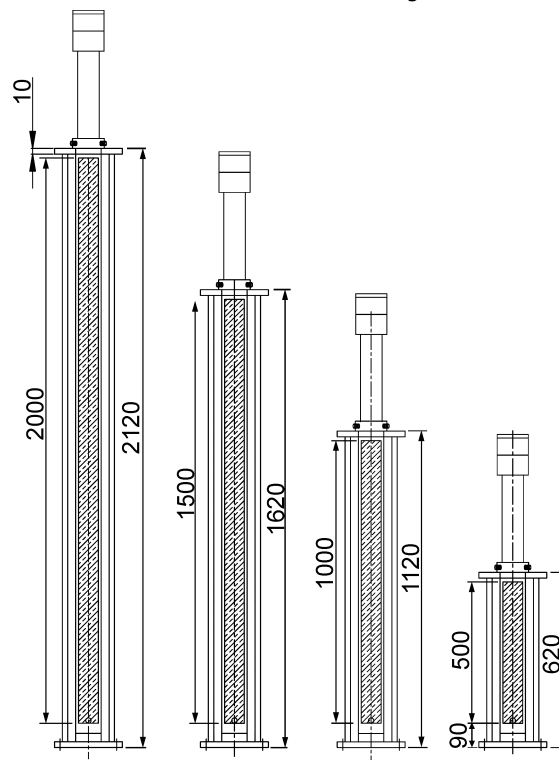
distance to the vessel wall
approx. 100 mm
Abstand zur Behälterwand
ca. 100 mm

1.5 Clamping Position for Rod Detectors Klemmenposition für Stabdetektoren

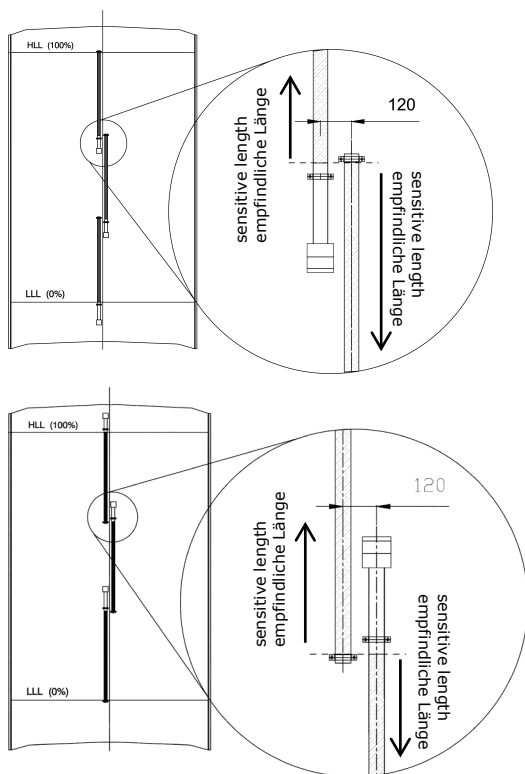
The drawings on this page are valid for Detectors with and without water cooling jacket.
Die Zeichnungen auf dieser Seite sind gültig für Detektoren mit und ohne Wasserkühlung.



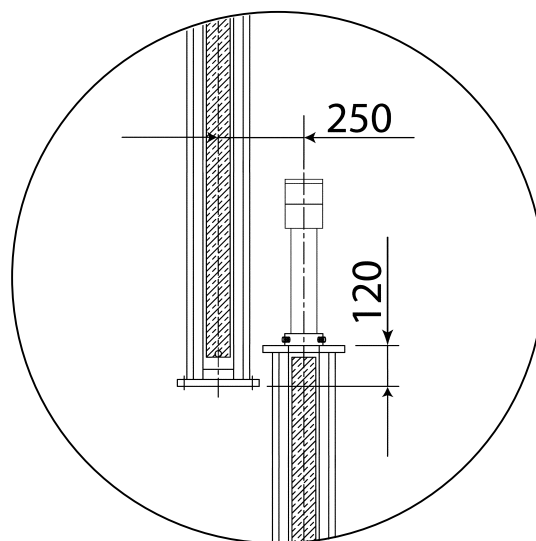
1.6 Mounting Position for Rod Detector Shieldings Klemmenposition für Stabdetektorabschirmungen



1.7 Clamping Position for Multi Detector Arrangement Klemmenposition für Multidetektor-Anordnung



1.8 Mounting Position for Rod Detector Shieldings Klemmenposition für Stabdetektorabschirmungen



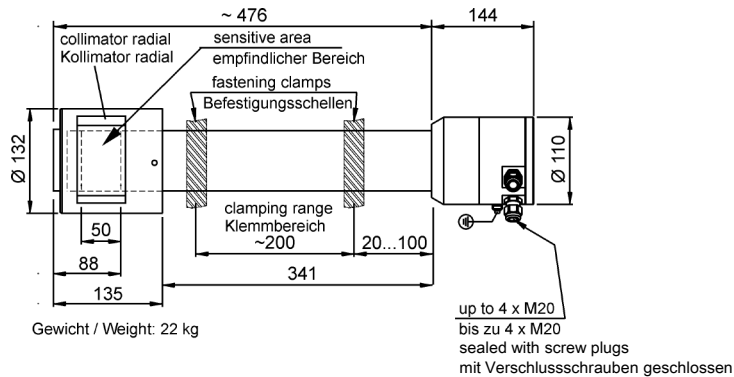
LB 480

Dimensions in mm
Abmessungen in mm

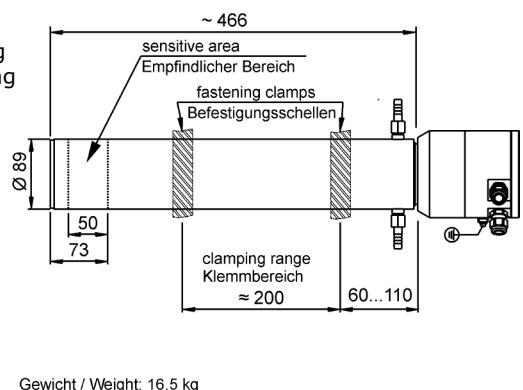
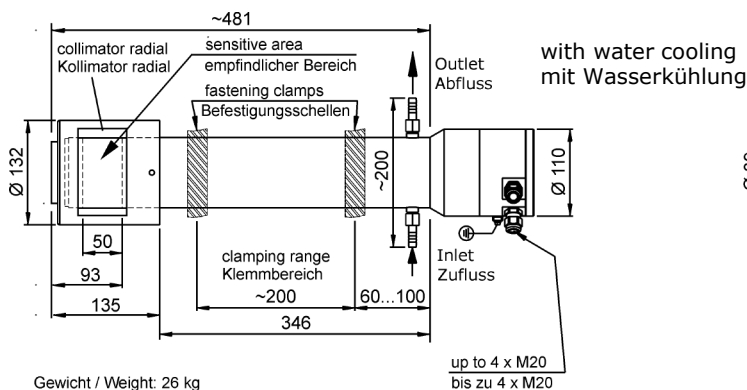
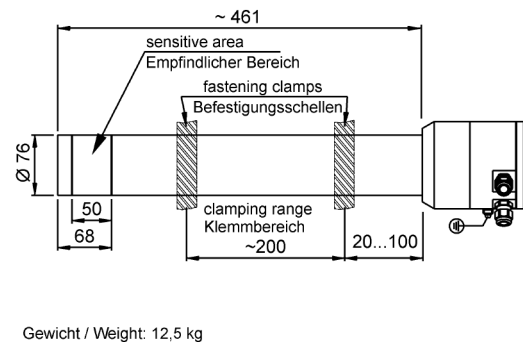
2.0 CrystaSENS (Version f. Zone 1/2)

LB 480-1.-0.
LB 480-1.-1.
LB 480-1.-2.
LB 480-1.-Z.

LB 480-1.-... -.-r.-

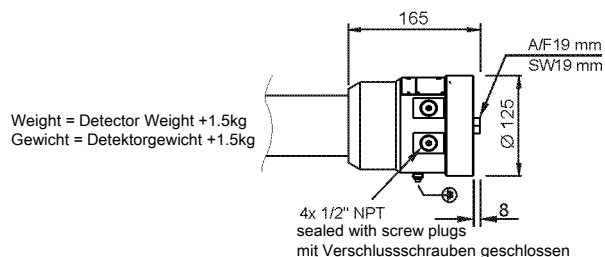


LB 480-1.-... -.-0.-



2.1 CrystaSENS (Version f. Divisions 1/2)

LB 480-1.-F.
LB 480-1.-G.



The detector version for divisions (NEC/CEC) differs only in the terminal housing, compared to the standard version illustrated above. The dimensions of this terminal housing are illustrated in this drawing.

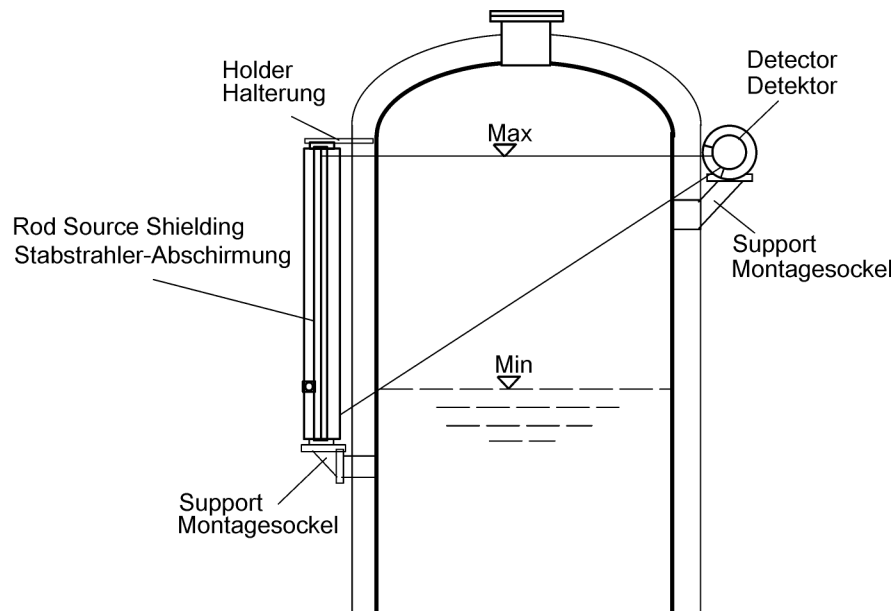
Die Detektor-Version mit Ex-Zulassung für Divisions (NEC/CEC) unterscheidet sich ausschließlich im Anschlusskopf, verglichen mit der oben aufgeführten Standardvariante. Die Abmessungen für den Anschlusskopf sind aus dieser Zeichnung zu entnehmen.

2.2 CrystaSENS Scintillator Size / Szintillatorgröße

Type Typ	Scintillator Size Szintillatorgröße	(Ø/h)	Water cooling Wasserkühlung
LB 480-11	50/50		-
LB 480-12	50/50		✓
LB 480-13	40/35		-
LB 480-14	40/35		✓
LB 480-15	25/25		-
LB 480-16	25/25		✓

LB 480

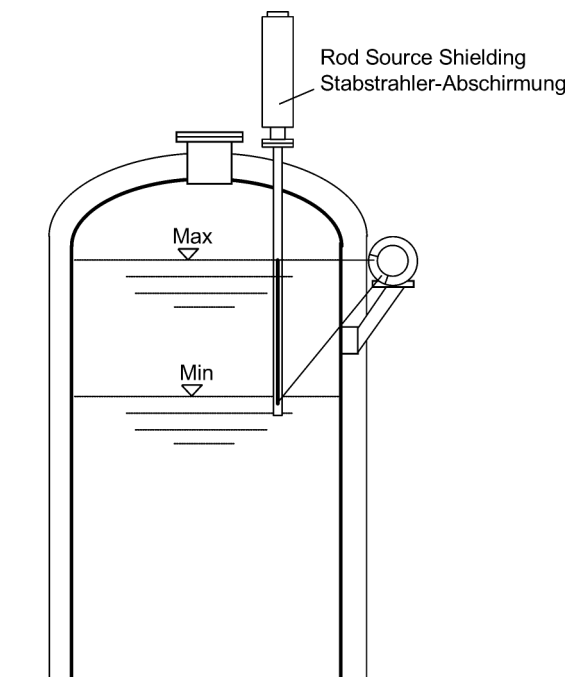
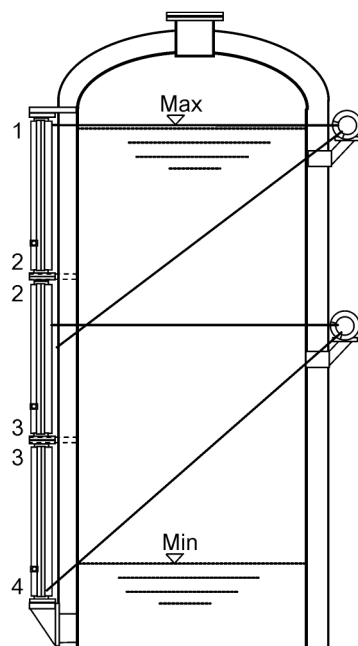
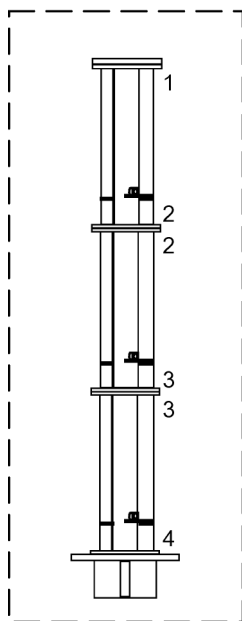
2.3 Point Detector Installation Examples Punktdetektor Installationsbeispiele



Application with multipart rod source shielding Anordnung mit mehrteiliger Stabstrahler-Abschirmung

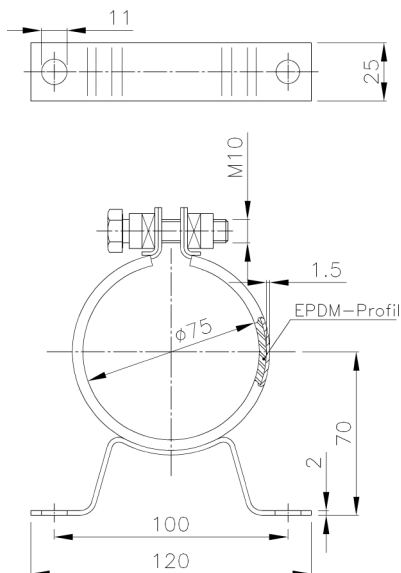
Application with dip pipe Anordnung mit Tauchrohr

Side View
Seitenansicht



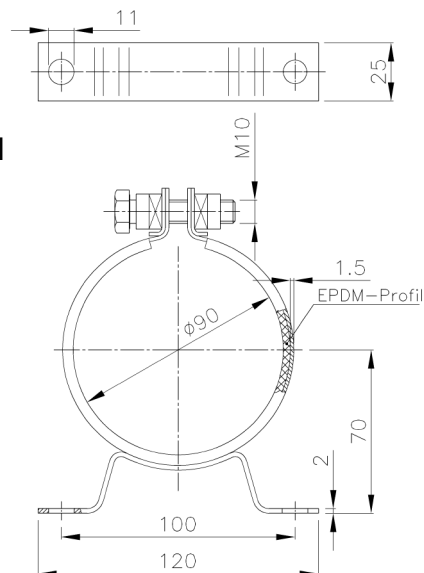
2.4 Mounting Clamps for Detector Befestigungsschellen für Detektor

for Detectors without water cooling
für Detektoren ohne Wasserkühlung



Id. Nr. 31346 (1 set = 2 clamps)
Id. Nr. 31345 (single clamp)

for Detectors with water cooling
für Detektoren mit Wasserkühlung



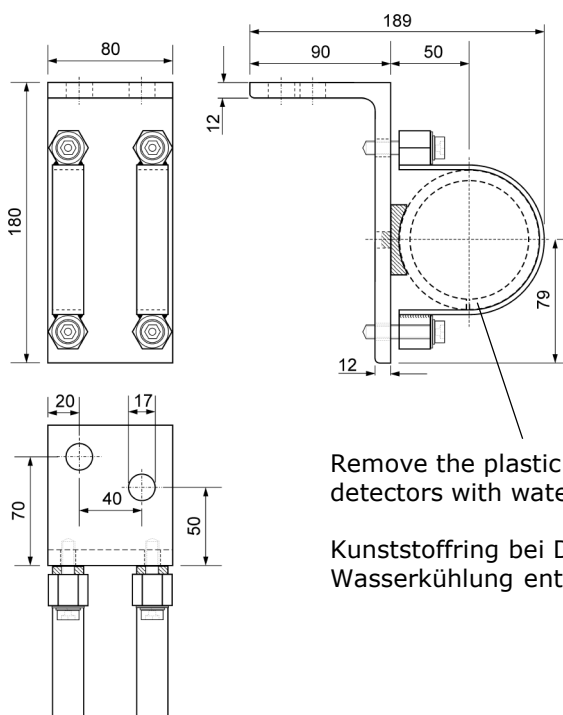
Id. Nr. 31347 (1 set = 2 clamps)
Id. Nr. 31344 (single clamp)

Material
316Ti
1.4571

position for the clamps, see detector drawing
Position für die Schellen-Befestigung siehe Detektor-Zeichnung

Heavy Duty Detector Holder (stainless steel) Robuste Detektor Halterung (Edelstahl)

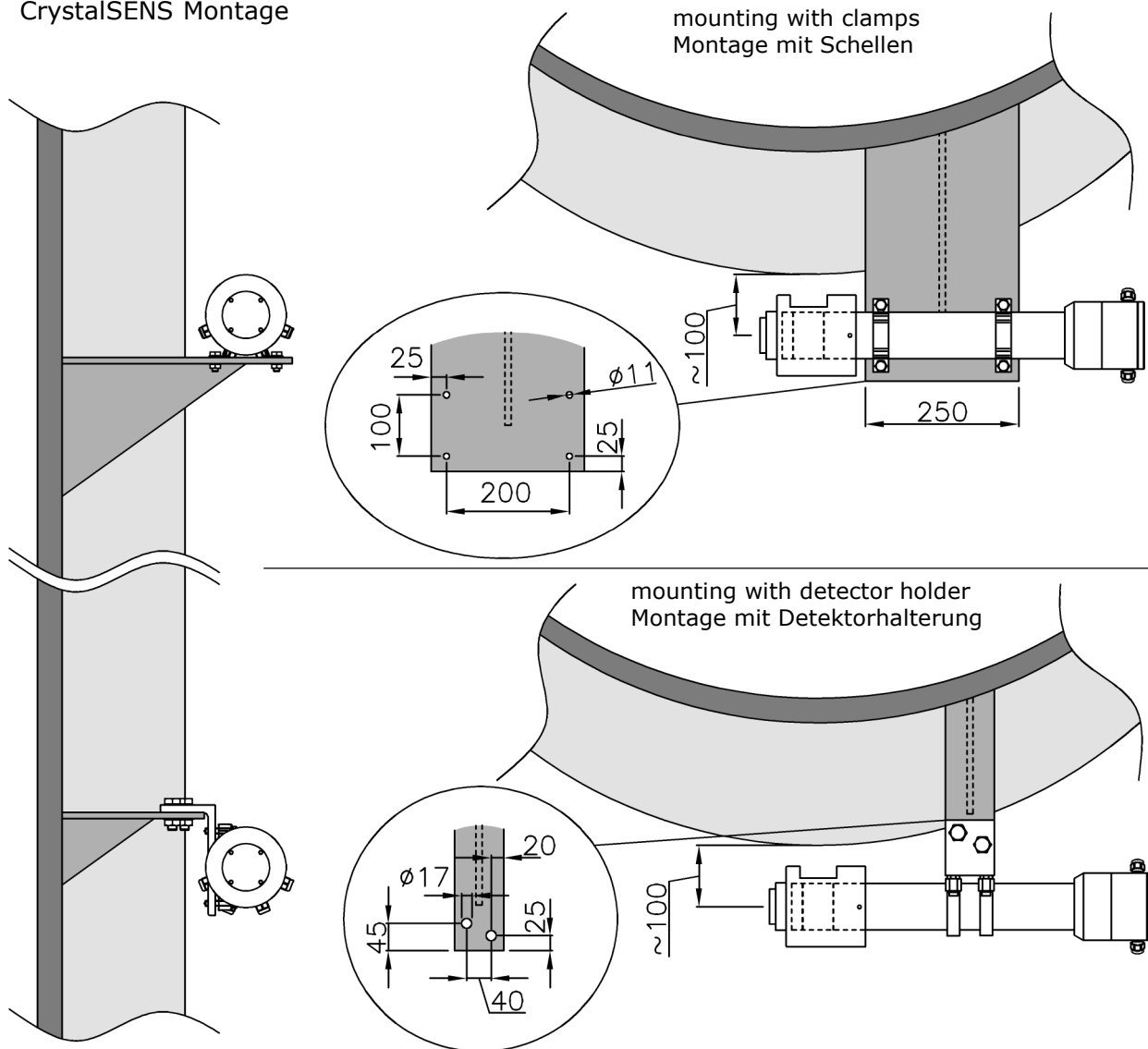
Part No. Id. Nr.	for Detector für Detektor
39246	without water cooling ohne Wasserkühlung
39247	with water cooling mit Wasserkühlung



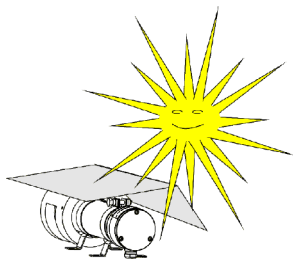
Remove the plastic ring for
detectors with water cooling.

Kunststoffring bei Detektoren mit
Wasserkühlung entfernen.

2.5 CrystalSENS Mounting CrystalSENS Montage



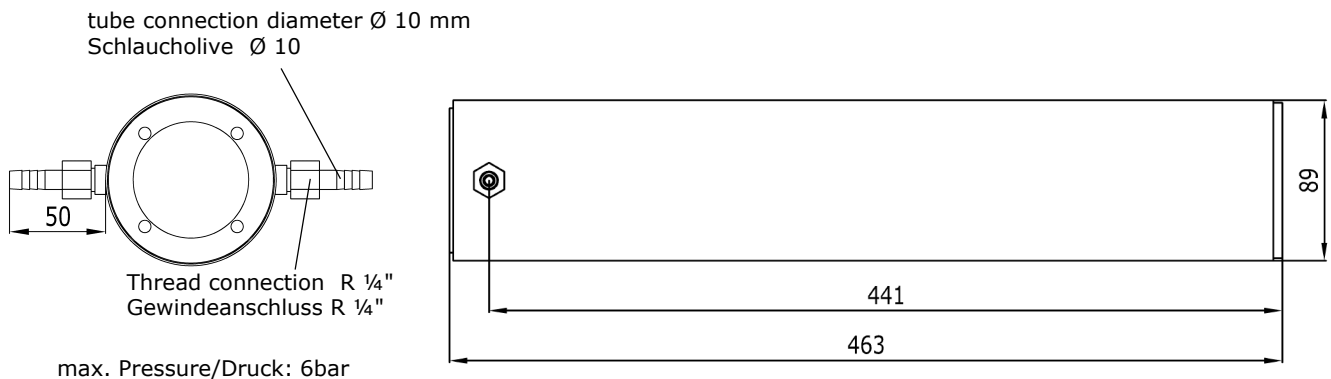
2.6 Sun Roof against Strong Sun Radiation Sonnendach gegen starke Sonneneinstrahlung



Direct sun radiation can overheat the detector. If the detector temperature can reach more than 50°C, a suitable sun roof must be installed. The heating of the detector by thermal radiation from the vessel can also be moderated by a thermal sheet, e.g. by a thin metal plate. For each detector a water cooling (option) is available.

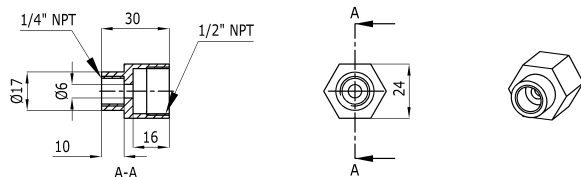
Wird durch Sonneneinstrahlung eine Detektortemperatur von über 50°C erreicht, so ist ein geeigneter Sonnenschutz zu montieren. Auch die Aufheizung des Detektors durch Wärmeabstrahlung vom Behälter kann durch ein dünnes Wärmeableitblech gemildert werden. Für jeden Detektor steht auch eine geeignete Wasserkühlung (Option) zur Verfügung.

2.7 CrystalSENS Water Cooling Jacket and Adaptor Fittings CrystalSENS Wasserkühlung und Adapter Anschlussstücke



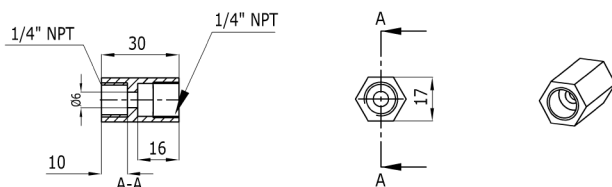
Fitting adaptor for standard water cooling Rp 1/4" → 1/2" NPT
stainless steel 304, part no: 47189

Adapter für Standard Wasserkühlung Rp 1/4" → 1/2" NPT
Edelstahl 1.4301, Id. Nr.: 47189



Fitting adaptor for standard water cooling Rp 1/4" > 1/4" NPT
stainless steel 304, part no: 46743

Adapter für Standard Wasserkühlung Rp 1/4" > 1/4" NPT
Edelstahl 1.4301, Id. Nr.: 46743



Further fitting adaptors for standard water cooling jacket:
Rp 1/4" > 1/2" NPT male, stainless steel 304, part no: 06352
Rp 1/4" > 1/4" NPT male, stainless steel 304, part no: 06349

Weitere Adapter für die Standard-Wasserkühlung:
Rp 1/4" > 1/2" NPT Außengewinde, 1.4301, Id. Nr.: 06352
Rp 1/4" > 1/4" NPT Außengewinde, 1.4301, Id. Nr.: 06349

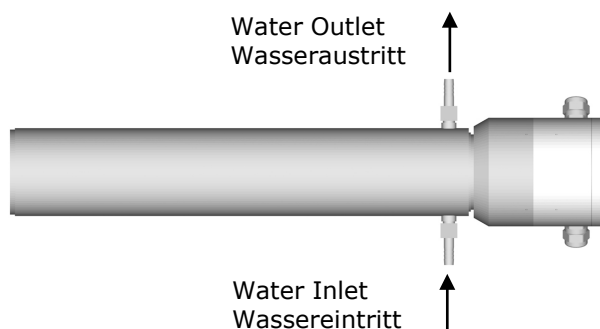
The above mentioned water cooling jackets and adaptor fittings offers following connection versions:
Die oben aufgeführten Wasserkühlungen und Adapter ermöglichen folgende Anschlussvarianten:

Fitting Connection	Anschluss-Stutzen	part no. (material) Id.Nr. (Werkstoff)
R 1/4" pipe connection, male European standard Whitworth pipe thread	R 1/4" Außengewinde für Rohrverschraubung europäisches Standard Whitworth-Rohrgewinde	21326 (304/1.4301) 38055 (Carbon Steel St37)
10 mm hose connection for water hose connection ID 10 mm	Schlauchstutzen für Schlauch- Innendurchmesser 10 mm	21326 (304/1.4301) 38055 (Carbon Steel St37)
fitting adaptor 1/2" NPT female	Adapter mit 1/2" NPT Innengewinde	47189 (304/1.4301)
fitting adaptor 1/4" NPT female	Adapter mit 1/4" NPT Innengewinde	46743 (304/1.4301)
fitting adaptor 1/2" NPT male	Adapter mit 1/2" NPT Außengewinde	06352 (304/1.4301)
fitting adaptor 1/4" NPT male	Adapter mit 1/4" NPT Außengewinde	06349 (304/1.4301)

LB 480

2.8 Water Cooling Installation Instruction Anweisung zur Installation der Wasserkühlung

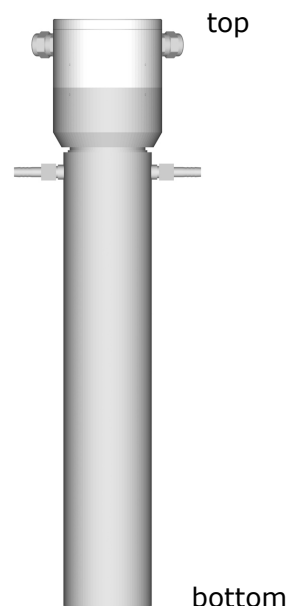
Horizontal Detector Installation Horizontale Detektor Installation



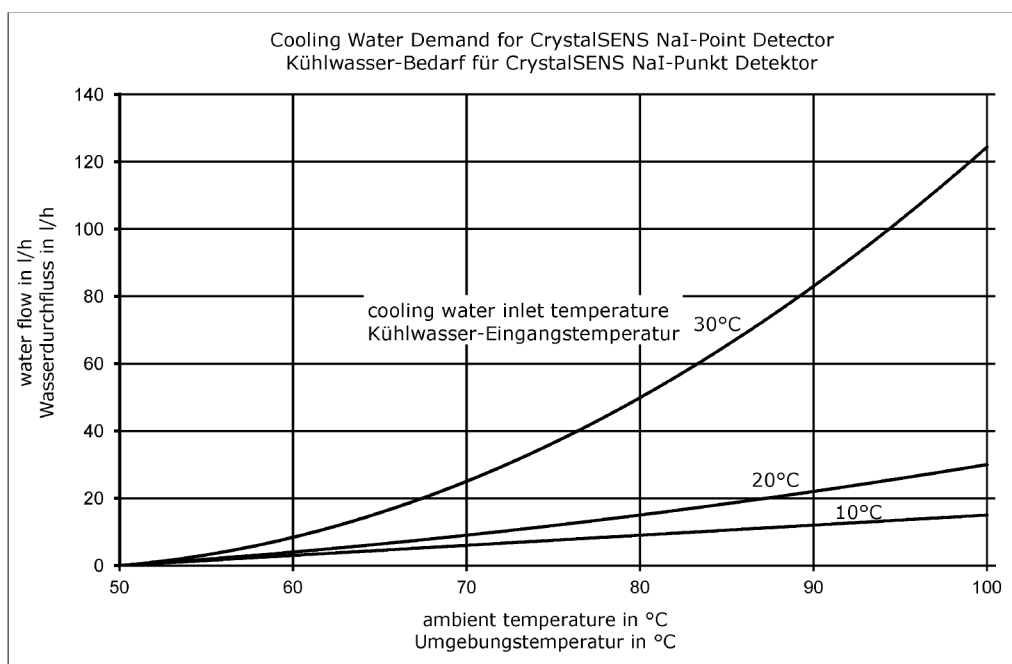
In order to fill the entire water cooling jacket, incoming water must enter from the bottom. Damit sich die Wasserkühlung vollständig mit Wasser füllt, muss der Wasserzufluss von unten erfolgen.

Vertical Detector Installation Vertikale Detektor Installation

Install the Detector with the connection at the top. Installieren Sie den Detektor mit dem Anschlussgehäuse oben.

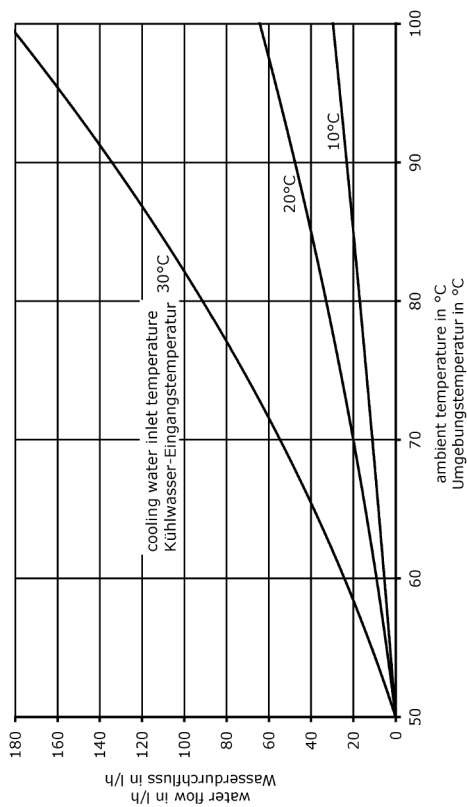


2.9 Detector Cooling Water Demand Detektor Kühlwasserbedarf

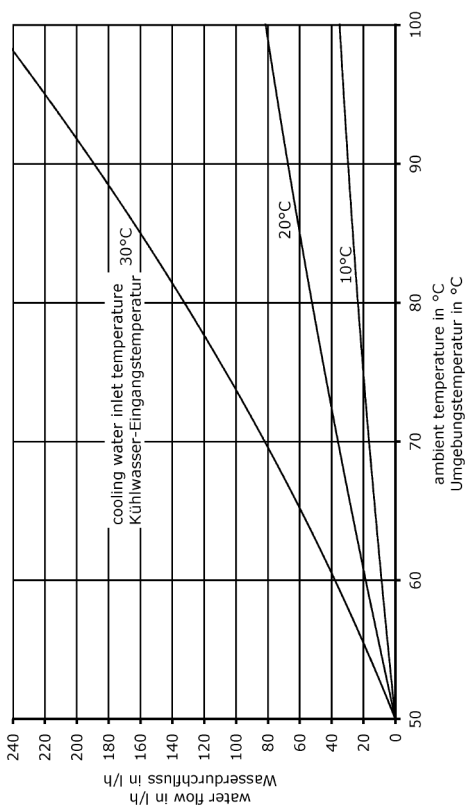


LB 480

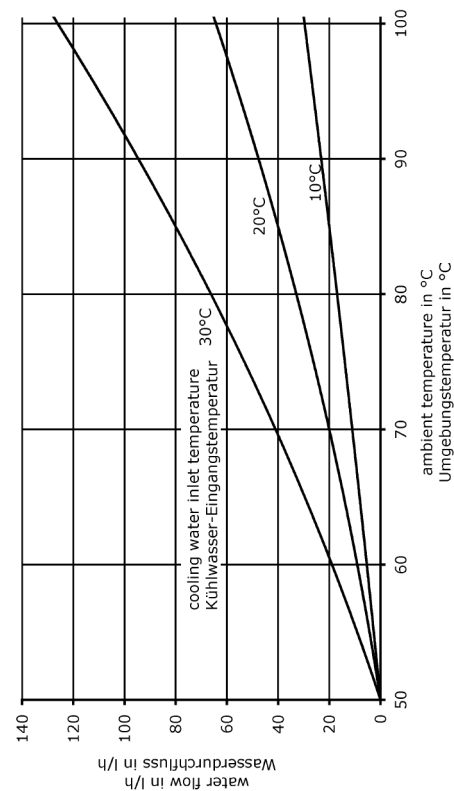
Cooling Water Demand for UniSENS 1000mm Rod Detector
Kühlwasser-Bedarf für UniSENS 1000mm Stabdetektor



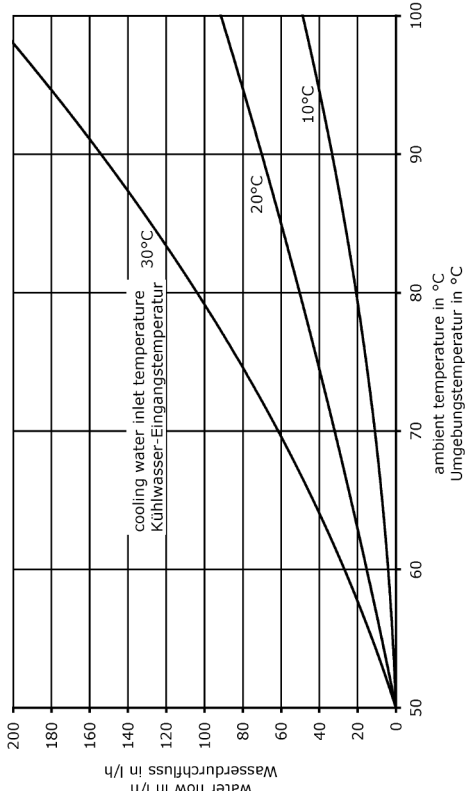
Cooling Water Demand for UniSENS 2000mm Rod Detector
Kühlwasser-Bedarf für UniSENS 2000mm Stabdetektor



Cooling Water Demand for UniSENS 500mm Rod Detector
Kühlwasser-Bedarf für UniSENS 500mm Stabdetektor



Cooling Water Demand for UniSENS 1500mm Rod Detector
Kühlwasser-Bedarf für UniSENS 1500mm Stabdetektor



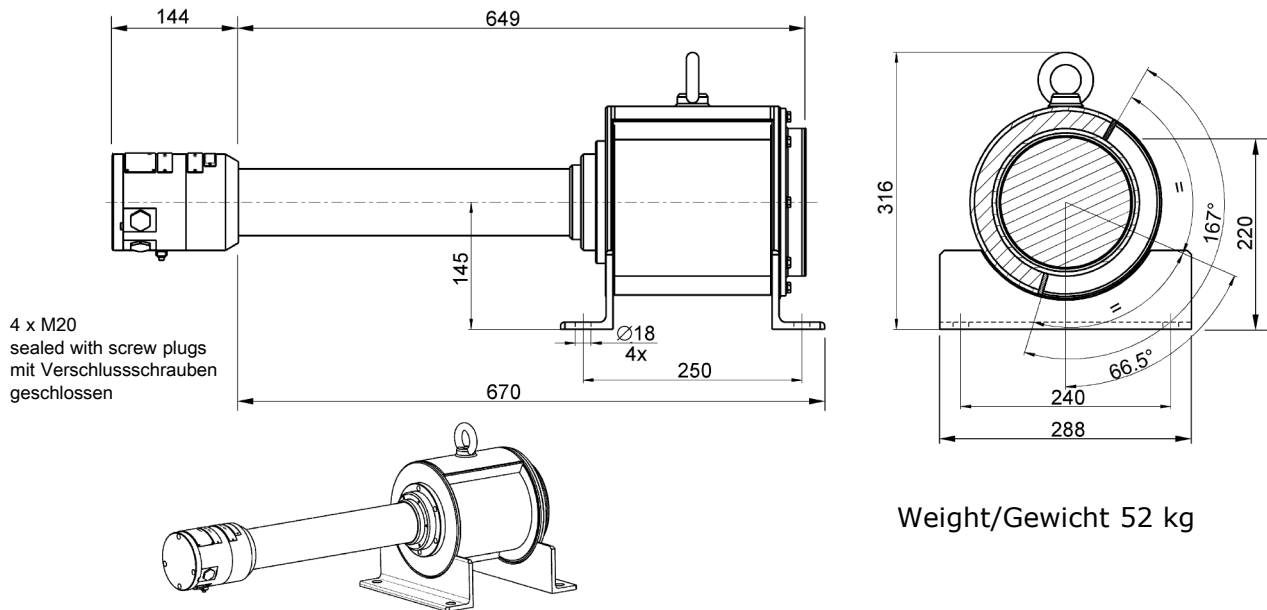
LB 480

Dimensions in mm
Abmessungen in mm

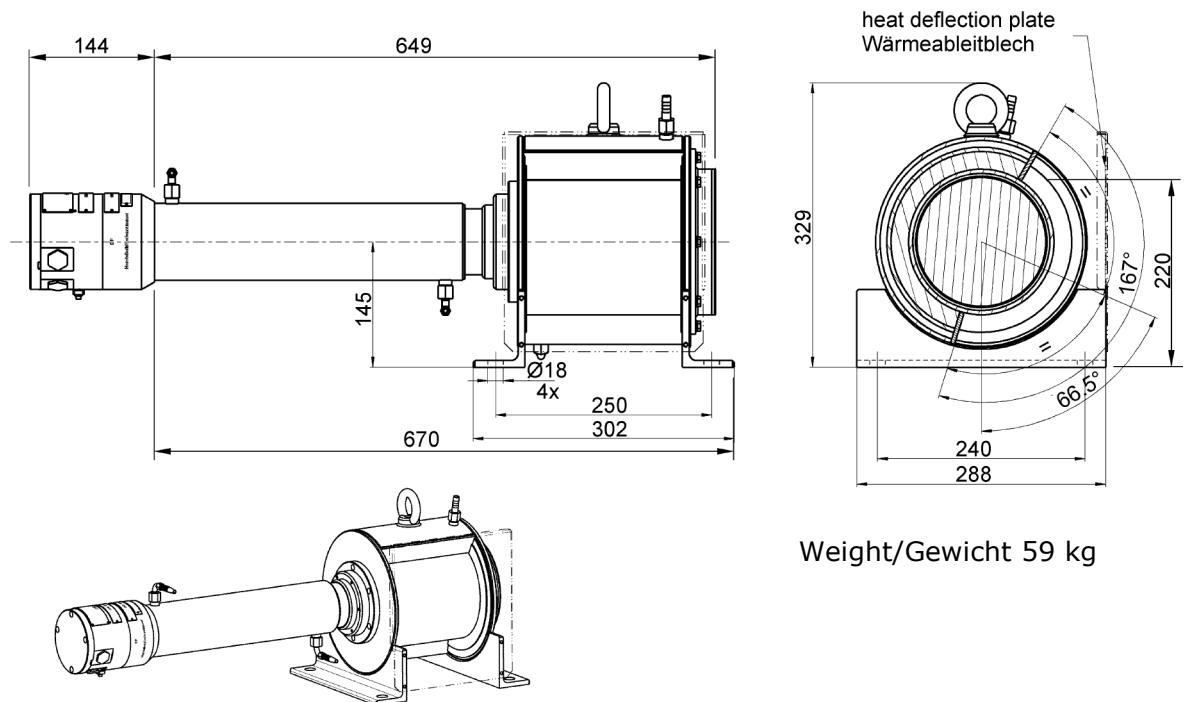
3. SuperSENS (Version f. Zone 1/2)

LB 480-3.-0.
LB 480-3.-1.
LB 480-3.-2.
LB 480-3.-Z.

3.1 with Side Irradiation mit seitlicher Einstrahlung LB 480-31-...-r.



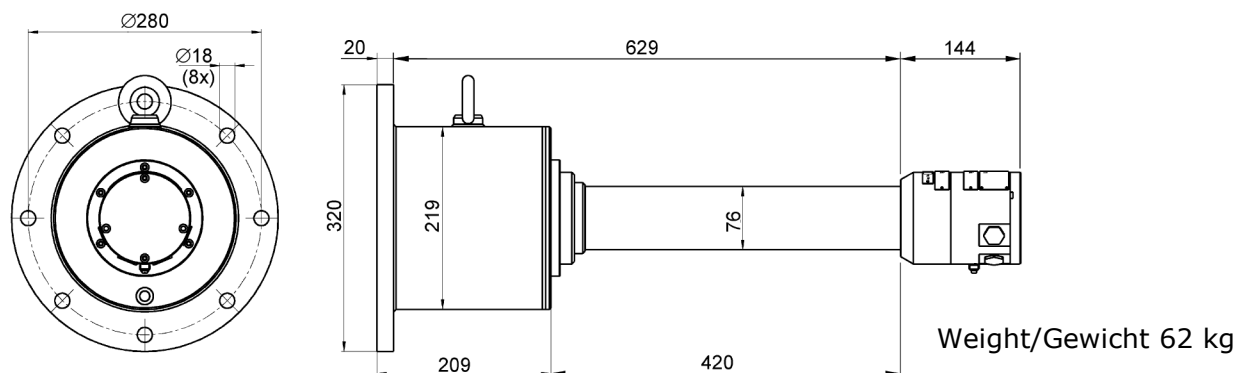
3.2 with Side Irradiation and Water Cooling mit seitlicher Einstrahlung und Wasserkühlung LB 480-32-...-r.



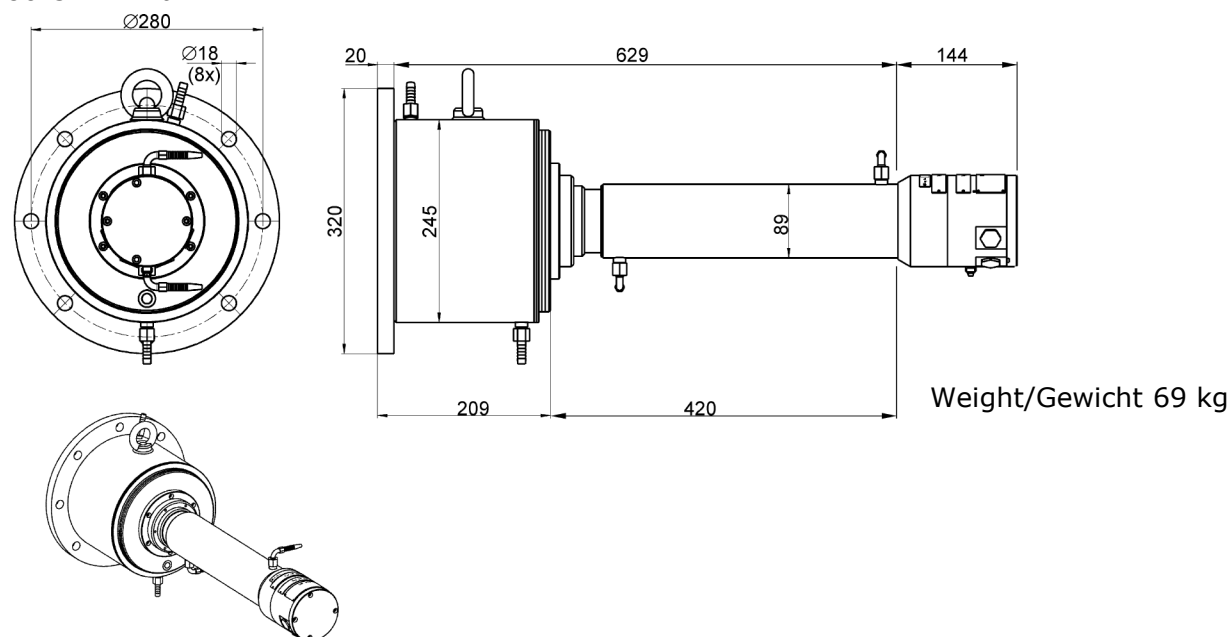
LB 480

Dimensions in mm
Abmessungen in mm

3.3 with Axial Irradiation mit frontaler Einstrahlung LB 480-31-...-a.

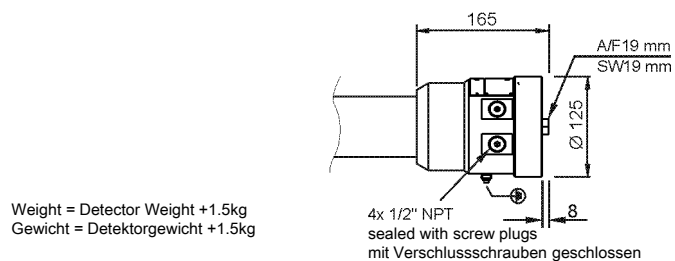


3.4 with Axial Irradiation and Water Cooling mit frontaler Einstrahlung und Wasserkühlung LB 480-32-...-a.



3.5 SuperSENS (Version f. Divisions 1 + 2)

LB 480-3.-F.
LB 480-3.-G.



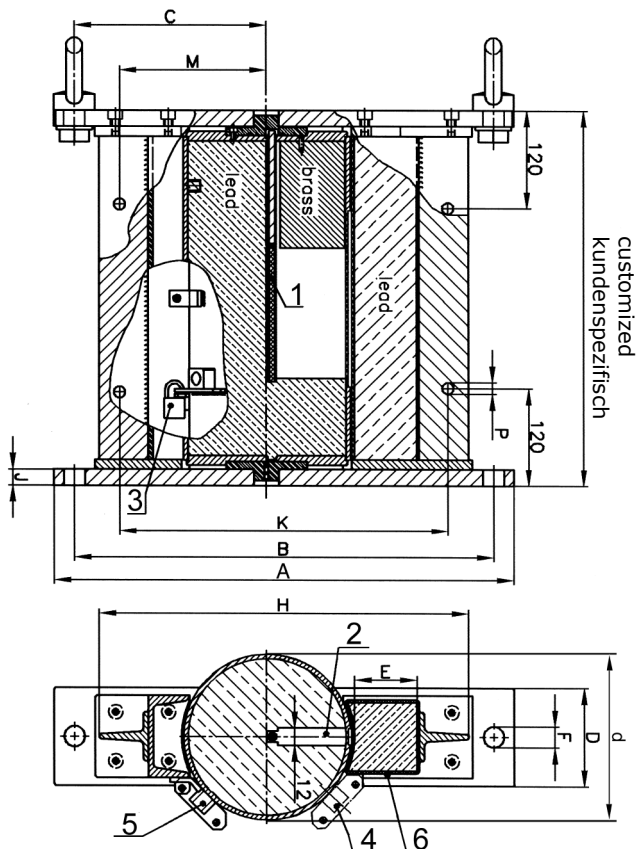
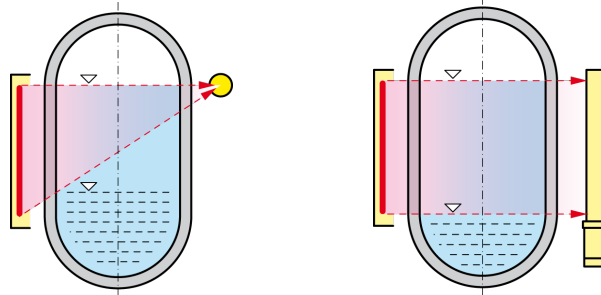
The detector version for divisions (NEC/CEC) differs only in the terminal housing, compared to the standard version illustrated above. The dimensions of this terminal housing are illustrated in this drawing.

Die Detektor-Version mit Ex-Zulassung für Divisions (NEC/CEC) unterscheidet sich ausschließlich im Anschlusskopf, verglichen mit der oben aufgeführten Standardvariante. Die Abmessungen für den Anschlusskopf sind aus dieser Zeichnung zu entnehmen.

LB 480

Dimensions in mm
Abmessungen in mm

4. Shielding for Rod Source Abschirmung für Stabstrahler



1	rod source	Stabstrahler
2	radiation channel	Strahlenaustritt
3	pad lock	Vorhängeschloss
4	position open	Stellung auf
5	position closed	Stellung zu
6	name plate	Typenschild

Multiple Arrangement of Rod Sources Mehnteilige Anordnung von Stabstrahlern

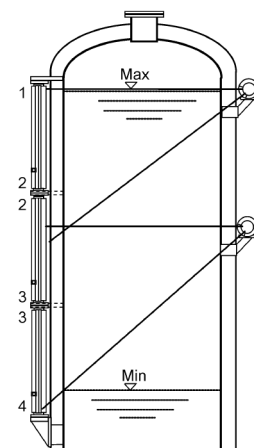
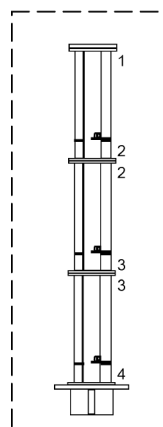
Holes for tilt protection
Bohrungen für Kippsicherung



The shields are marked with TOP and BOTTOM to rule out any side-inverted installation. Multi-part shieldings are additionally marked with the numbers 1, 2, 3, ... from top to bottom.

Die Abschirmungen sind markiert mit OBEN und UNTEN um eine seitenverkehrte Installation auszuschließen. Mehrteilige Abschirmungen sind zusätzlich von oben nach unten mit den Zahlen 1, 2, 3, ... markiert.

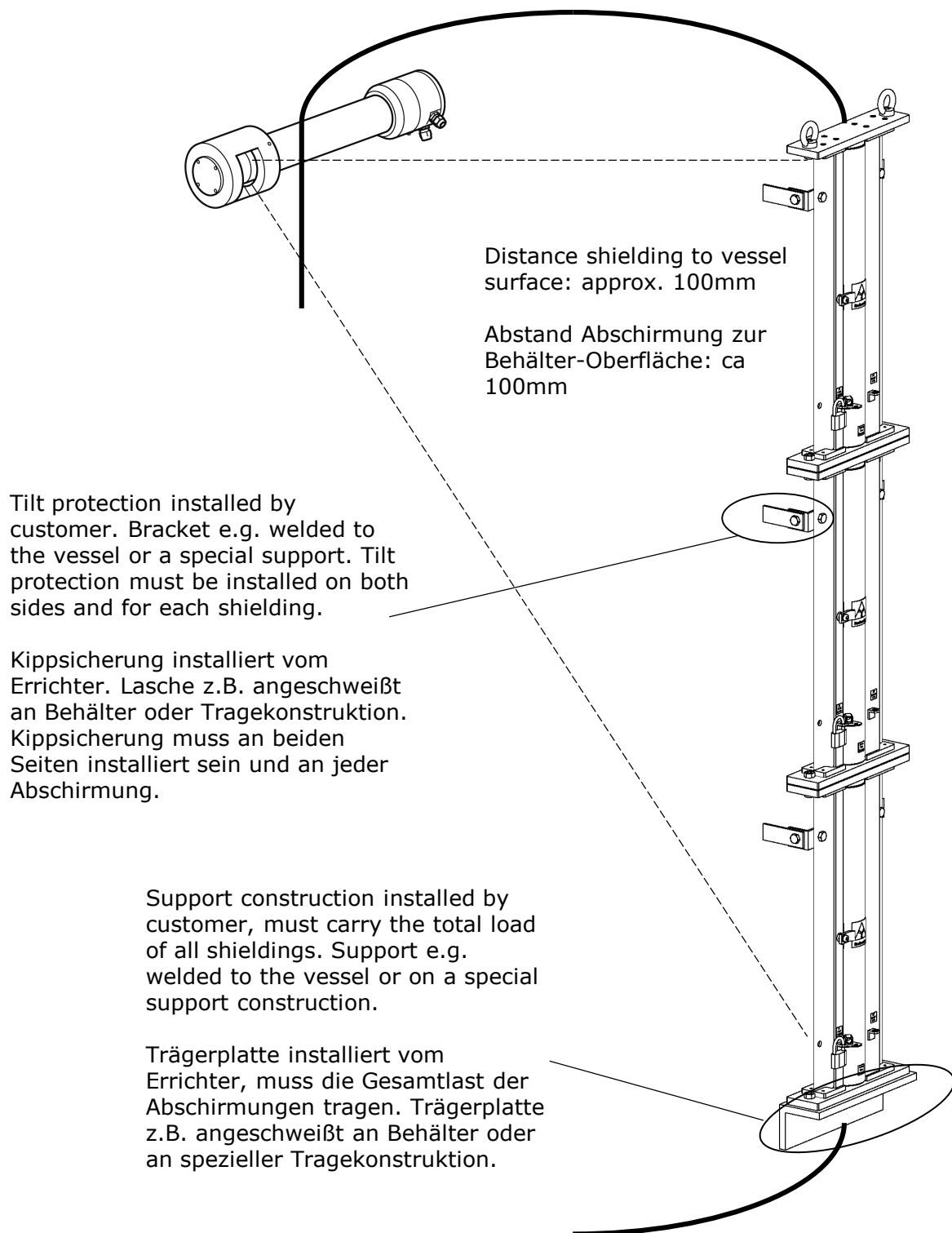
Side View
Seitenansicht



Type Typ	d	kg/m approx	A	B	C	D	E	F	H	J	K	M	P
80	85	100	320	285	130	80	41	14	235	15	204.5	90	11.5
100	105	130	340	305	140	80	41	14	255	15	224.5	100	11.5
120	127	190	400	360	165	80	47	18	310	20	270	120	11.5
150	159	300	465	425	195	100	65	18	375	20	334	150	14
200	203	450	560	510	235	120	75	25	450	25	400	180	14
270	267	780	580	510	230	125	100	26	545	25	503	222.5	14

LB 480

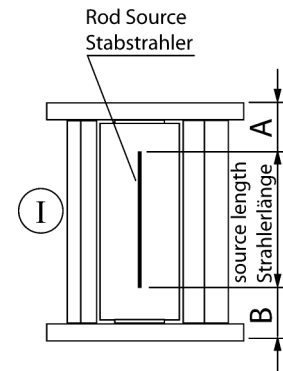
4.1 Multipart Installation of Rod Source Shieldings Mehrteilige Installation von Stabstrahler-Abschirmungen



4.2 Single Part Arrangement for Rod Source Shieldings Einteilige Anordnung von Stabstrahler-Abschirmungen

For single part rod source arrangement only the shielding with the "arrangement I" is necessary

Eine einteilige Stabstrahler-Anordnung benötigt lediglich die Abschirmung mit der Montageart I.



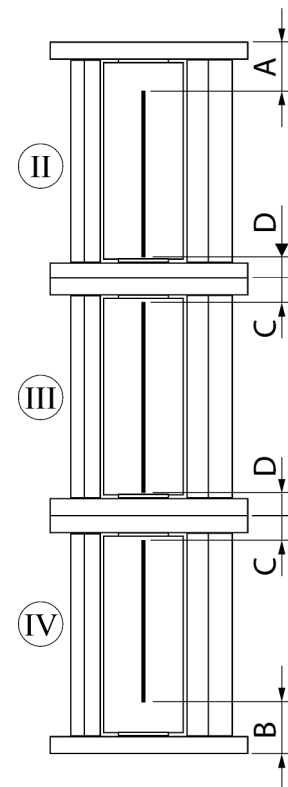
4.3 Multipart Arrangement for Rod Source Shieldings Mehrteilige Anordnung von Stabstrahler-Abschirmungen

In order to establish a continuous source length with a multiple source arrangement, different types of shieldings must be used for interconnection. The different shielding types are mentioned with roman figures in the drawing.

Damit eine mehrteilige Stabstrahler-Anordnung einen möglichst durchgängigen Strahler bildet, sind unterschiedliche Abschirmungen zum Zusammenbau zu verwenden. Die dazu notwendigen einzelnen Abschirmungstypen sind im Bild oben mit römischen Ziffern beschriftet.

2-part arrangement: II + IV
3-part arrangement: II + III + IV
4-part arrangement: II + III + III + IV
etc.

Anordnung mit 2 Abschirmungen: II + VI
Anordnung mit 3 Abschirmungen: II + III + IV
Anordnung mit 4 Abschirmungen: II + III + III + IV
etc.

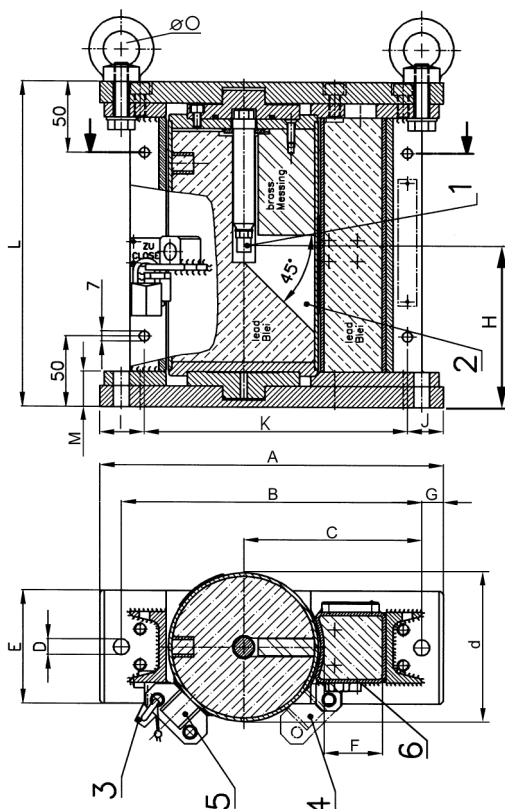
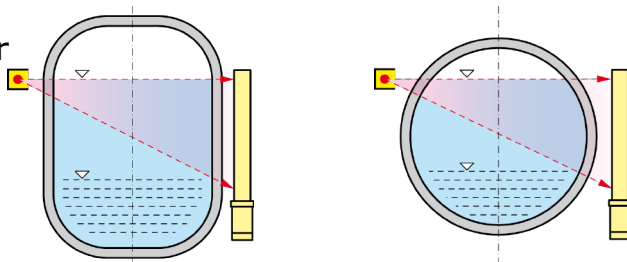


Shielding / Abschirmung		A	B	C	D
Ø	drawing no. Zeichnungs-Nr.				
80	21156.000-000	78	78	18	24
100	21157.000-000	92	92	18	24
120	21158.000-000	113	89	23	30
150	21159.000-000	132	106	21	32
200	21160.000-000	169	132	26	37
270	21161.000-000	202	140	30	40

LB 480

Dimensions in mm
Abmessungen in mm

5. Shielding for Point Source Abschirmungen für Punkt Strahler

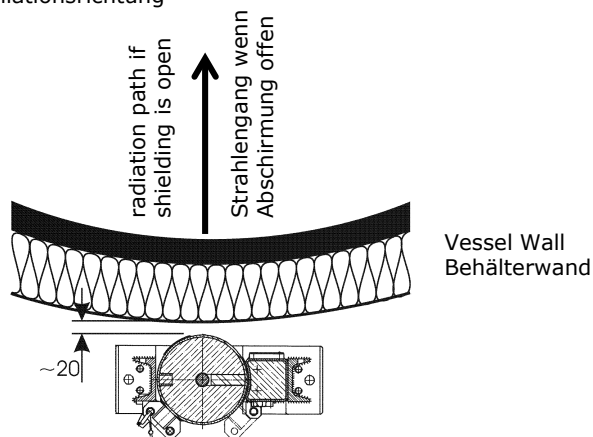


The shielding are marked with BOTTOM to rule out any side-inverted installation.

Die Abschirmungen sind markiert mit UNTEN um eine seitenverkehrte Installation auszuschließen.

1	point source	Punktstrahler
2	radiation channel	Strahlenaustritt
3	pad lock	Vorhängeschloss
4	position open	Stellung auf
5	position closed	Stellung zu
6	name plate	Typenschild

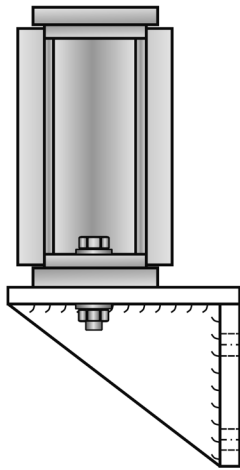
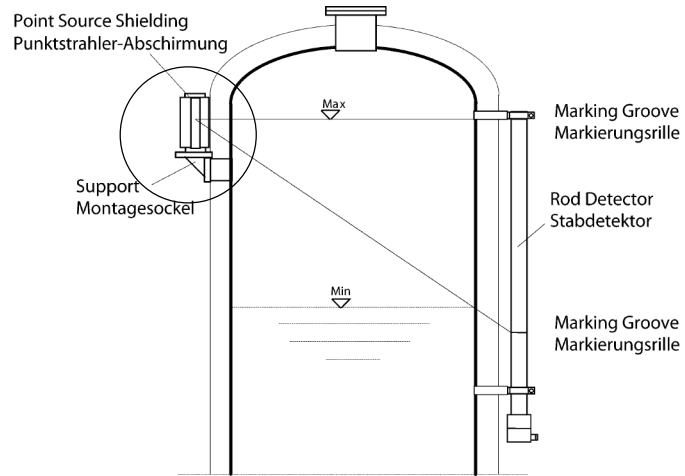
Direction of Installation
Installationsrichtung



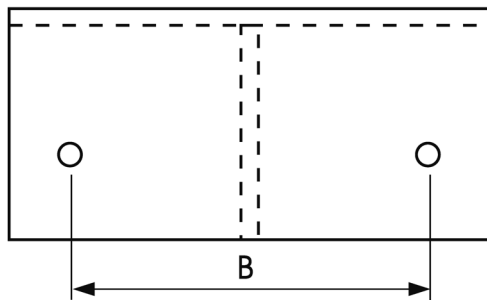
Type Typ	d	A	B	C	D	E	F	G	H	I	J	K	L	M	øO	Weight Gewicht
100	105	240	210	125	11	80	40	15	110	30	25	185	225	25	25	32 kg
150	159	360	320	195	18	100	65	20	150	29	29	302	265	25	35	70 kg
200	203	410	360	220	23	120	75	25	185	22	22	366	340	30	40	150 kg
270	267	580	510	315	26	300	101	35	280	45	55	436	480	45	50	370 kg

LB 480

5.1 Proposal for Support Vorschlag für Montagesockel

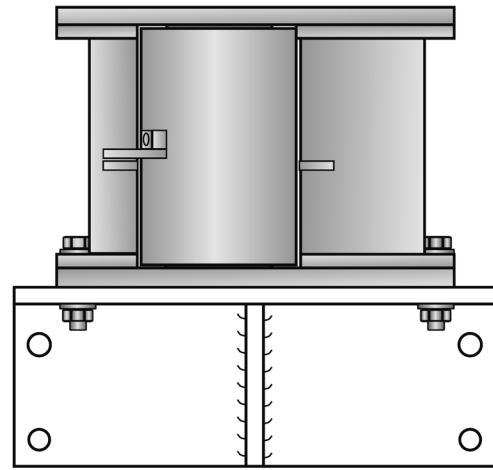


1

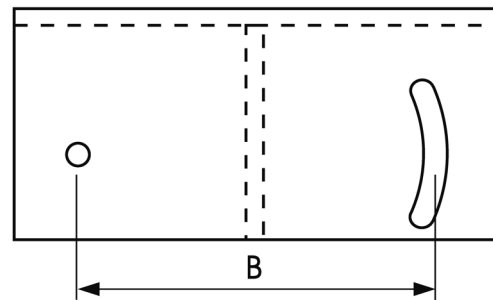


The design 1 of the support plate is easier to manufacture.

Der Montagesockel 1 ist leichter herzustellen.



2

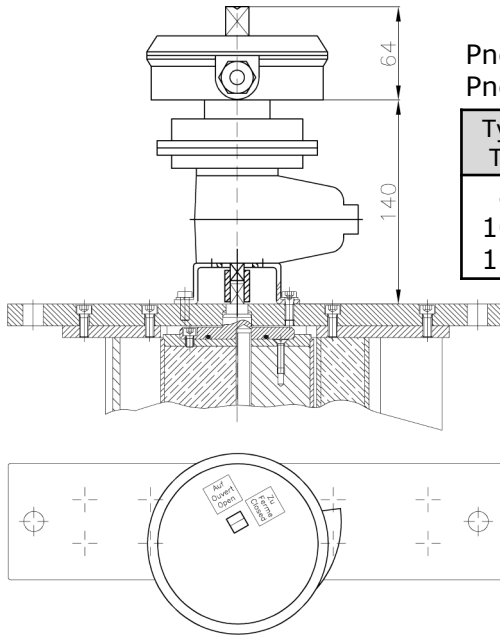


The design 2 of the support plate with the oblong slot, allows an adjustment of the radiation beam to the detector.

Die Ausführung 2 des Montagesockels mit dem Langloch, ermöglicht die Ausrichtung des Strahlers auf den Detektor.

LB 480

6. Pneumatic Actuator for Rod and Point Source Shielding (Type 80 ... 270) Pneumatik für Stab- und Punktstrahler-Abschirmungen (Typ 80 ... 270)



Pneumatic actuator for **rod source** shielding
Pneumatischer Antrieb für **Stabstrahler**-Abschirmung

Type Typ	without Limit-Switch ohne Endschalter	with Limit-Switch mit Endschalter	with Limit-Switch Exe mit Endschalter Exe
80	49943	49942	50013
100	38169	38171	51444
150	51446	51445	51316

Pneumatic actuator for **point source** shielding
Pneumatischer Antrieb für **Punktstrahler**-Abschirmung

Type Typ	without Limit-Switch ohne Endschalter	with Limit-Switch Exe mit Endschalter Exe
100	40085	40084
150	41689	41690
200	41681	41691
270	-	50531

Other shieldings with pneumatic on request.
Andere Abschirmungen mit Pneumatik auf Anfrage.

For remote control a pneumatic shutter system is available as an option. The pneumatic is available with a limit switch open/close indication. The limit switch box is available in IP 65 or as an Ex-proofed system.

Attention: Multipart rod source shieldings can not be mounted on top of one another, but must be arranged sidewise out-of-line. For each shielding a separate pneumatic actuator is necessary. For the rod source shielding with pneumatic, use the shielding type I only.

Data for Pneumatic Actuator

Compressed Air: min. 4×10^5 Pa (4 bar) max. 7×10^5 Pa (7 bar) connection: G 1/8
Air Quality: clean as usual for compressed air tools, oil free
Temperature Range: -20°C ... + 80°C

Indication Open/Closed via Limit Switch

Option I: IP65, 2 contacts (Open/Closed) 48 VDC, 1 A
Option II: 2 contacts (Open/Closed) max. 250V AC, 1 A Ex-Protection of the limit switches: EEx d IIC T6 Ex-Protection of the limit switch housing: EEx e IIC T6

Sofern Sie den Verschlussmechanismus fernsteuern wollen ist eine pneumatische Verschlusseinrichtung optional erhältlich. Die Pneumatik ist zusätzlich mit Endschalter für die Positionsrückmeldung erhältlich.

Achtung: mehrteilige Stabstrahler-Abschirmungen mit Pneumatik können nicht direkt aufeinander montiert werden, sondern müssen seitlich zueinander versetzt angeordnet werden. Für jede Abschirmung wird eine kpl. Pneumatik benötigt. Die Pneumatik darf nur auf Abschirmungen vom Typ I montiert werden.

Daten für pneumatischen Verschlussantrieb

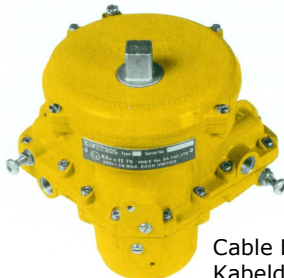
Druckluft: min. 4×10^5 Pa (4 bar) max. 7×10^5 Pa (7 bar) Anschluss: G 1/8
Luftqualität: Saubere wie bei Druckluft-Werkzeugen üblich, ölfrei
Temperaturbereich: -20°C ... + 80°C

Signalisierung AUF/ZU durch Endschalter

Option I: IP65, 2 Kontakte (AUF/ZU) 48 VDC, 1 A
Option II: 2 Kontakte (AUF/ZU) max. 250V AC, 1 A Schutzart der Microeinbautaster: EEx d IIC T6 Gehäuseschutzart: EEx e IIC T6

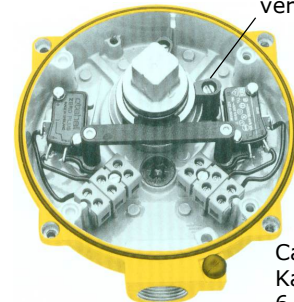
6.1 Components for Pneumatic on Rod and Point Source Shielding Einzelteile für Pneumatik an Stabstrahler- und Punktstrahlerabschirmung

Limit Switch
Endschaltereinheit
Ex de Zone 1



Cable Diameter
Kabeldurchmesser
6 ... 12 mm

Limit Switch
Endschaltereinheit
IP 65 (SPDT)



adjustable cams
verstellbare Nocken

Cable Diameter
Kabeldurchmesser
6 ... 12 mm

Contact Rating
Kontaktbelastbarkeit

Volt		Load / Last (A)	
AC	DC	R	L
250		7	5
125		7	5
	12	7	5
	24	1	1
	48	0,5	0,06
	250	0,25	0,03

Position marked with a red coloured cut in the top.
Position mit eingefrästem roten Strich markiert.

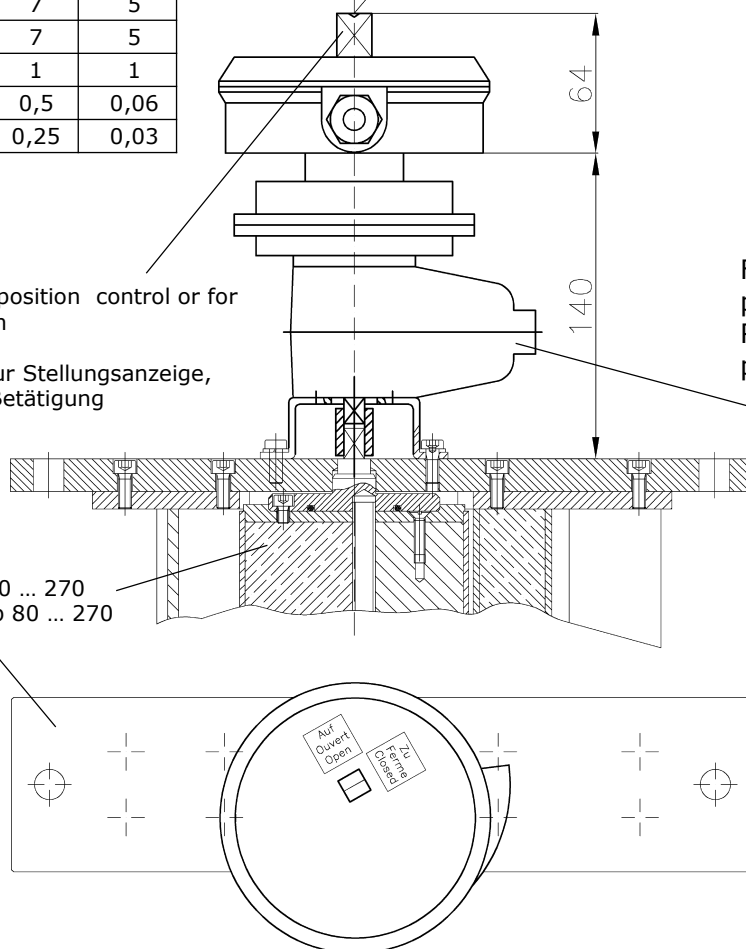
Contact Rating
Kontaktbelastbarkeit

Volt		Load / Last (A)		
AC	DC	R	L	Lamp
250		15	3	1,5
125		15	3	1,5
	12	15	3	1,5
	24	10	2	1
	48	3	0,6	0,3
	250	0,25	0,05	0,025

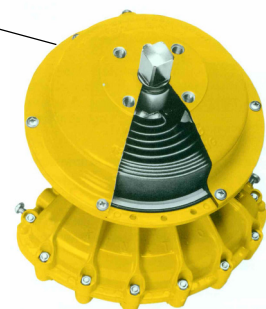
Square neck for position control or for manual operation

Außenvierkant zur Stellungsanzeige, oder manueller Betätigung

Shielding Type 80 ... 270
Abschirmung Typ 80 ... 270



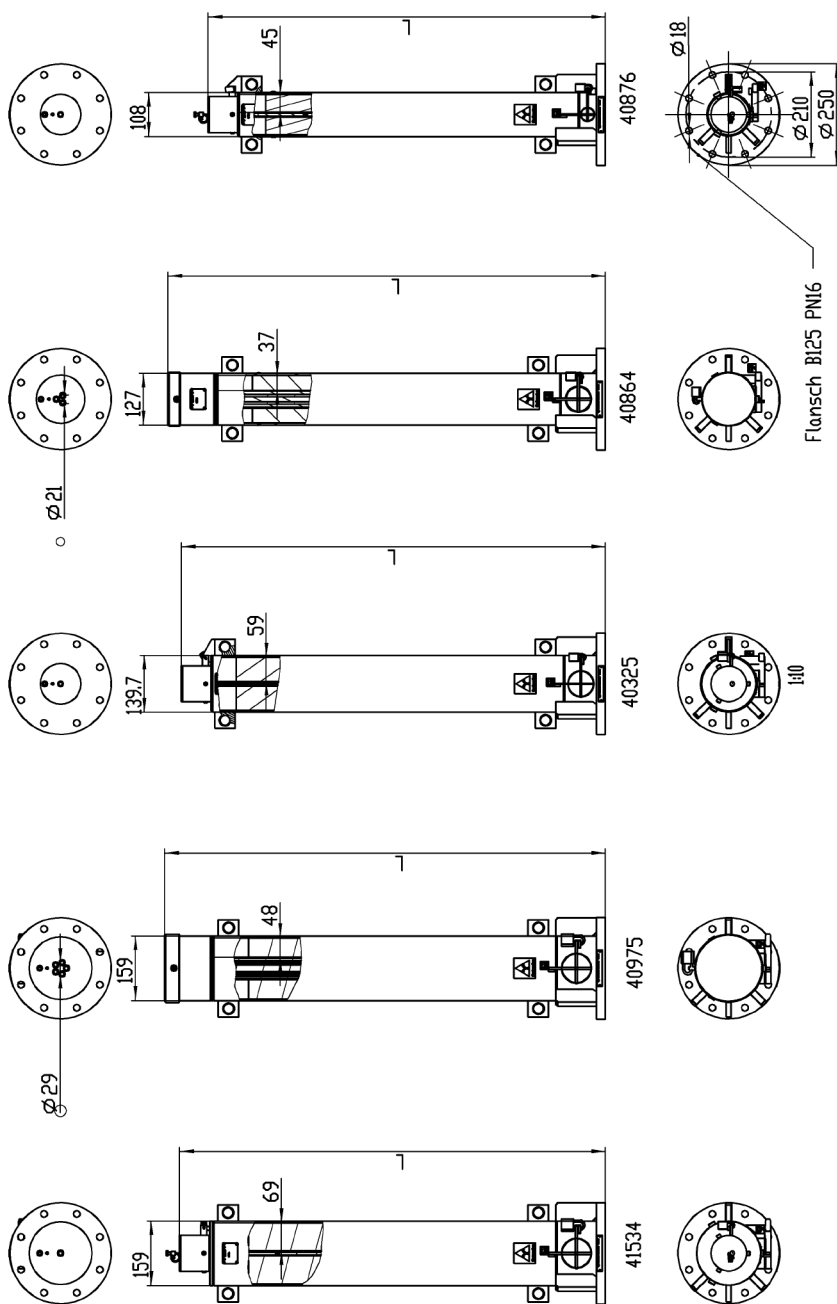
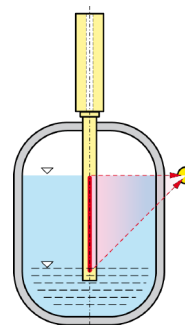
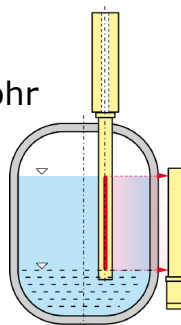
Fail safe spring return with pneumatic actuator.
Rückstellfeder (Fail-Save) mit pneumatischem Antrieb



LB 480

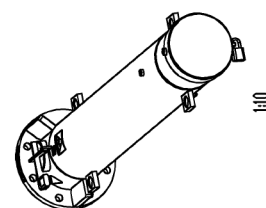
Dimensions in mm
Abmessungen in mm

7. Flange Shielding for Rod Source on Dip Pipe Abschirmungen für Stabstrahler am Tauchrohr



CAUTION
Do not close the shutter, if the source is in measuring position.
Do not remove the steel cord until the source is in home position and the shutter is closed.

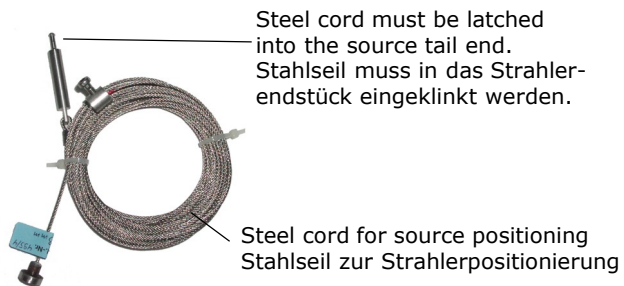
ACHTUNG
Bei Strahler in Messposition, den Verschluss nicht betätigen!
Vor dem Entfernen des Seiles die Strahler sichern und Verschluss schließen.



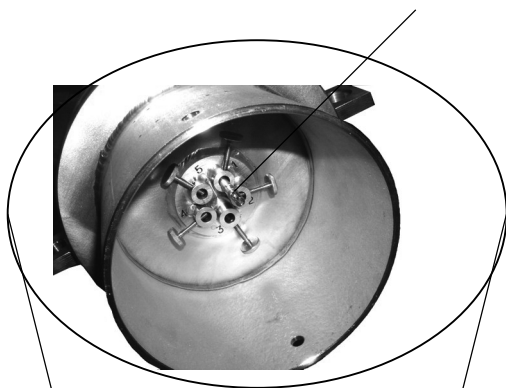
part no. Id. Nr.	41534	40975	40325	40864	40876
effective lead thickness wirksame Bleidicke	69	47.5	58.85	36.5	45
max. Anzahl der Strahler max. number of sources	1	5	1	3	1
Source length [mm]	1046	1085	1042	1075	976
Strahlerlänge [mm]	1246	1285	1242	1275	1176
	1446	1485	1442	1475	1376
	1646	1685	1642	1675	1576
	1846	1885	1842	1875	1776
	2046	2085	2042	2075	1976
	2246	2285	2242	2275	2176

LB 480

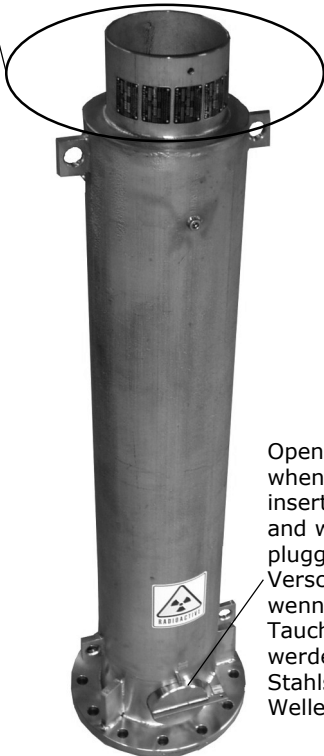
7.1 Operation of Flange Shielding Bedienung der Flanschabschirmung



source tail end (inactive)
Strahlerendstück (inaktiv)



shielding head
Abschirmungskopf

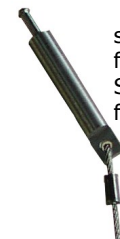


Open the shutter only when the source shall be inserted into the dip pipe and when the steel cord is plugged in.
Verschluss erst öffnen wenn der Strahler in das Tauchrohr gefahren werden soll und das Stahlseil, bzw. die Wellenseele eingeklinkt ist.

Operation of the coupling device Bedienung der Kupplung

A)

source tail end with catch bolt
Strahlerendstück mit Kupplung



steel cord with latch for catch bolt
Stahlseil mit Klinke für Kupplung

B)

plug the latch into the catch bolt
Klinke in Kupplung einklinken



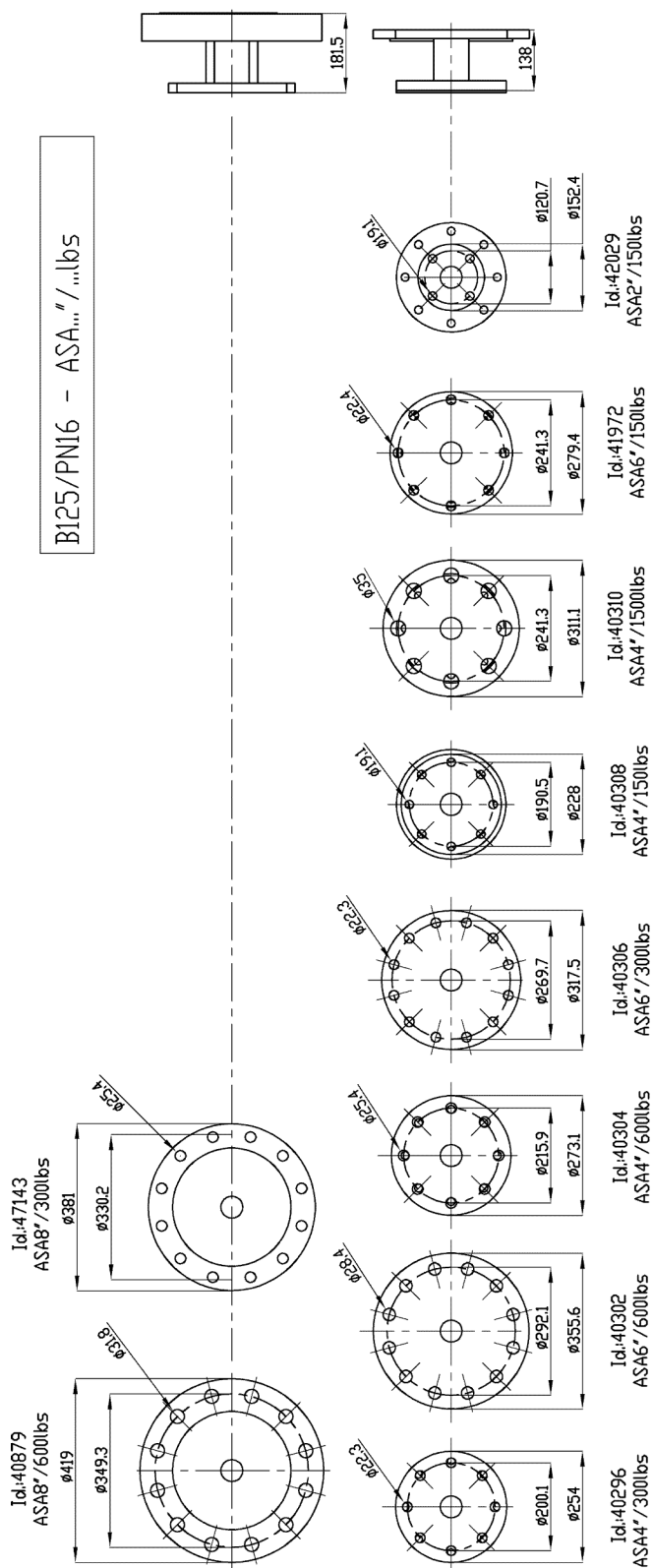
C)

latch plugged into the catch bolt
Klinke in Kupplung eingeklinkt

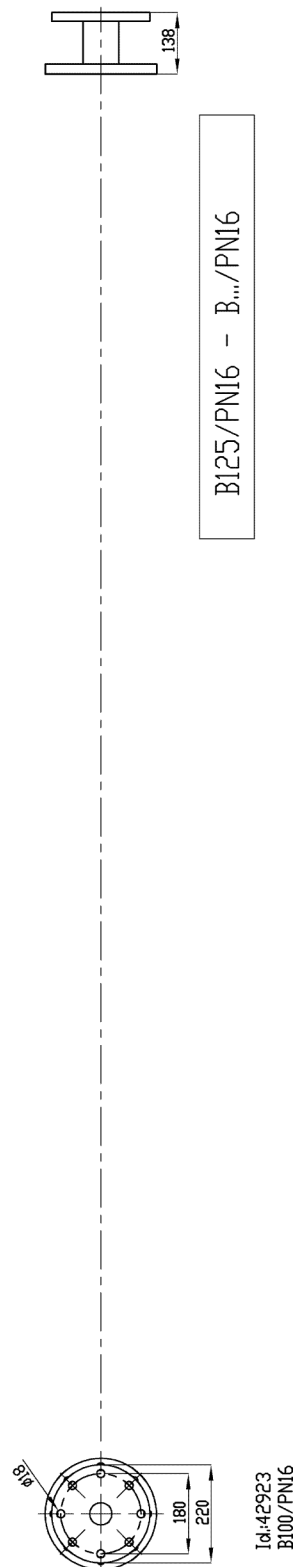


LB 480

7.2 Flange Adaptors for Rod Source Shieldings Flanschadapter für Stabstrahler-Abschirmungen



B125/PN16 - ASA.../...lbs

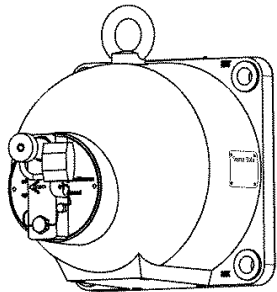
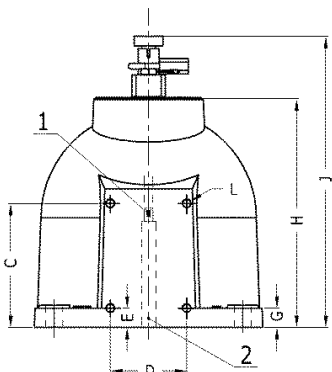
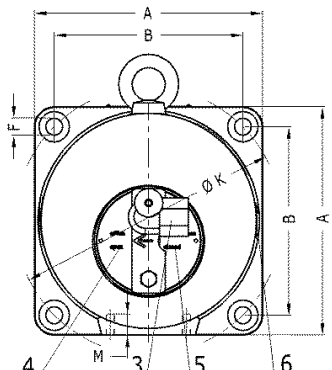
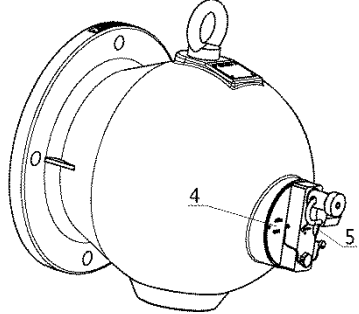
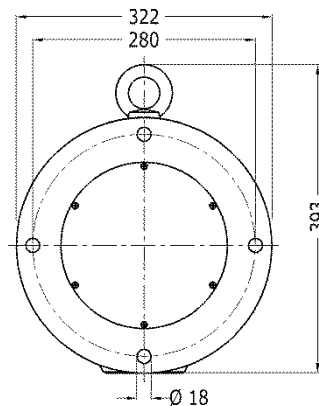
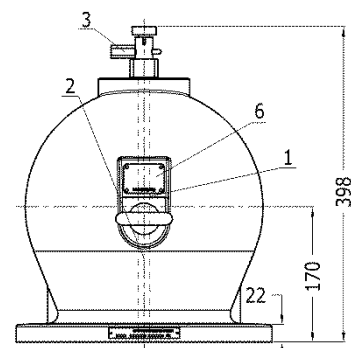


B125/PN16 - B.../PN16

LB 480

Dimensions in mm
Abmessungen in mm

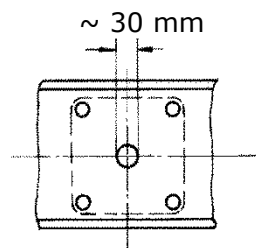
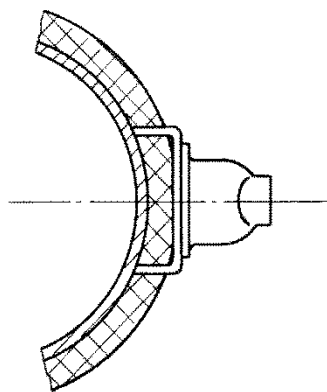
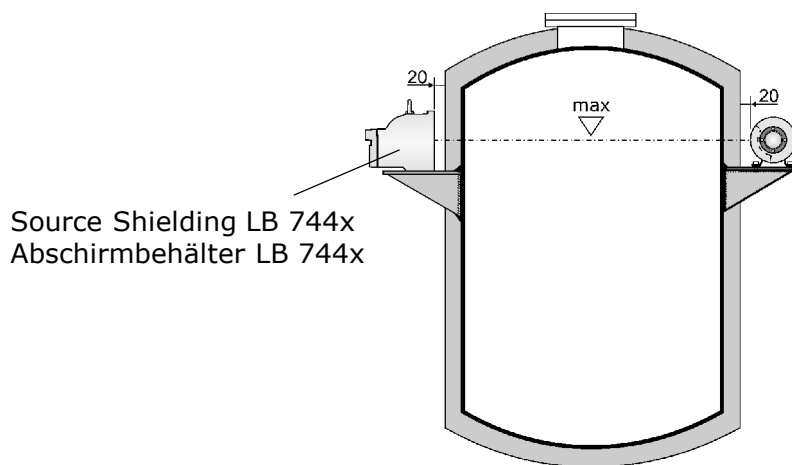
8. Point Source Shielding LB 744x Punktstrahler-Abschirmbehälter LB 744x

LB 7440 / 7442 / 7445 / 7446		LB 7444	
  		  	
1 Point Source 2 Radiation Exit 3 Lock 4 Position Open 5 Position Closed 6 Type Label 1 Strahler 2 Strahlenausgang 3 Schloss 4 Position Offen 5 Position Geschlossen 6 Typenschild			

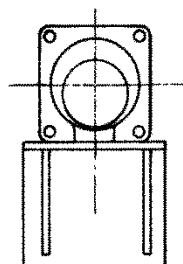
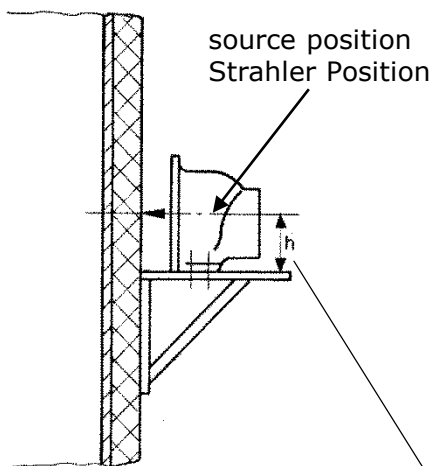
Housing Gehäuse	Type Typ	A	B	C	D	E	FØ	G	H	J	KØ	L	M		kg
Carbon Steel St 37	LB 7440-F-CR LB 7445-F-CR	180	142	75	60	15	18	20	173	238	200	M8	12	16°	31
Super Duplex UNS 32750 SAF 2507 1.4410	LB 7440-FE-CR LB 7445-FE-CR														
	LB 7442-FE-CR	240	198	130	80	20	18	20	242	306	280	M10	14	9°	81
Carbon Steel St 37	LB 7442-F-CR LB 7446-F-CR														
	LB 7444-CR	Dimensions in drawing / Abmessungen in Zeichnung												6°	170

 Radiation Angle of the Shielding / Abstrahlwinkel der Abschirmung

8.1 Mounting Proposal for Source Shielding LB 744x Montagevorschlag für Abschirmbehälter LB 744x



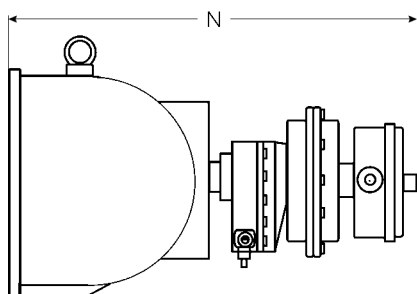
Flange Installation
Flanschinstallation



Pedestal Installation
Sockelinstallation

	LB 7440 LB 7445	LB 7442 LB 7446	LB 7444
h	90	120	161

8.2 Pneumatic Actuator for Source Shielding LB 744x Pneumatischer Verschlussantrieb für Abschirmbehälter LB 744x



N approx / ca.	Type / Typ
390	LB 7440-F-CR LB 7440-D-CR LB 7440-FE-CR LB 7440-DE-CR
460	LB 7442-F-CR LB 7442-D-CR LB 7442-FE-CR LB 7442-DE-CR
570	LB 7444-CR

Protection Schutz	Part No. Ident.	Description Beschreibung
IP 65	36119	Pneumatic Actuator with Limit Switch Option I Pneumatischer Verschlussantrieb mit Endschalter Option 1
Ex de IIC T6	80919	Pneumatic Actuator with Limit Switch Option II Pneumatischer Verschlussantrieb mit Endschalter Option II
Ex ia IIC T6	43698	Pneumatic Actuator with Limit Switch Option III Pneumatischer Verschlussantrieb mit Endschalter Option III

Technical Data for Pneumatic Actuator Technische Daten für pneumatischen Verschlussantrieb

Compressed Air Druckluft	min. 4×10^5 Pa (4 bar) max. 7×10^5 Pa (7 bar) Connection / Anschluss: G 1/8
Air Quality Luftqualität:	clean as usual for air compressed tools, oil free Saubere wie für Druckluft-Werkzeuge üblich, ölfrei
Temperature Range Temperaturbereich:	-20°C ... +80°C

Limit Switch, Option for Signaling OFF / CLOSED Endschaltereinheit, Optionen für Signalisierung AUF / ZU

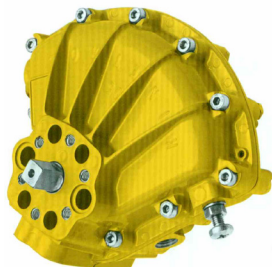
Option I:	IP 65, 2 contacts (OFF/CLOSED) IP 65, 2 Kontakte (AUF/ZU)
Option II:	2 contacts (OFF/CLOSED) Protection for internal micro switches: EEx d IIC T6 Housing protection: EEx e II T6 2 Kontakte (AUF/ZU) Schutzart der Microeinbautaster: EEx d IIC T6 Gehäuseschutzart: EEx e II T6
Option III:	2 Proximity switches EEx ia IIC T6, for intrinsically safe power supply 2 Näherungsinitiatoren EEx ia IIC T6, für eigensichere Speisung

LB 480

Dimensions in mm
Abmessungen in mm

8.3 Components for Pneumatic Actuator Einzelteile für Pneumatischen Verschlussantrieb

Pneumatic Actuator
Pneumatischer Antrieb

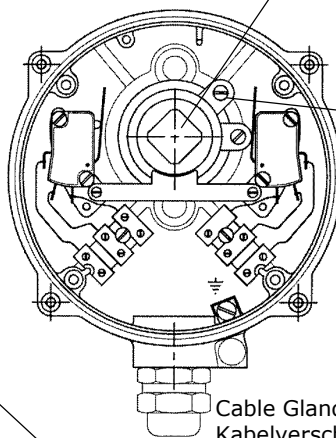
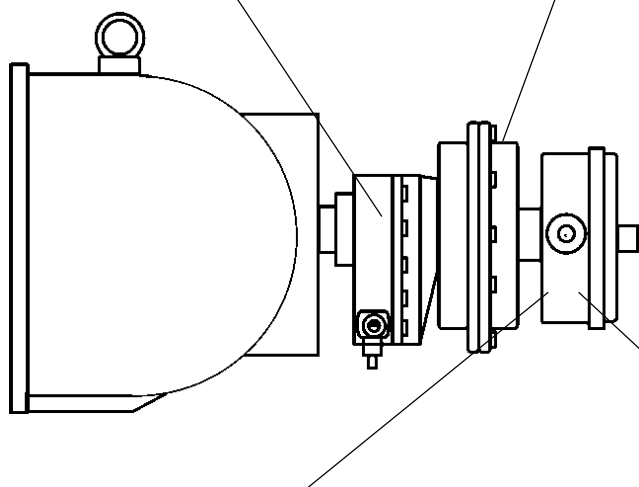


Return Spring
Rückstellfeder



male square spin-type handle
as position indicator, or for
manual operation

Außenvierkant zur Stel-
lungsanzeige, oder manueller
Betätigung



adjustable cam
verstellbare Nocken

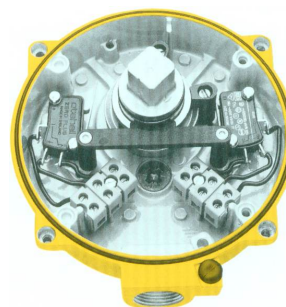
Cable Gland
Kabelverschraubung

Limit Switch in Ex de
Endschalterereinheit in Ex de



Cable Diameter 6 ... 12 mm
Kabeldurchmesser 6 ... 12 mm

Limit Switch in IP 65
Endschalterereinheit in IP 65



Cable Diameter 9 ... 12 mm
Kabeldurchmesser 9 ... 12 mm

Contact Rating
Kontaktbelastbarkeit

Volt		Load / Last (A)	
AC	DC	R	L
250		7	5
125		7	5
	12	7	5
	24	1	1
	48	0,5	0,06
	250	0,25	0,03

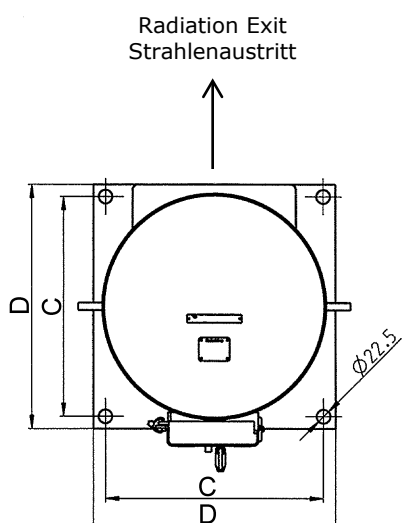
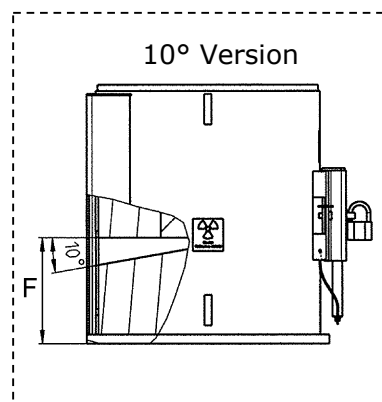
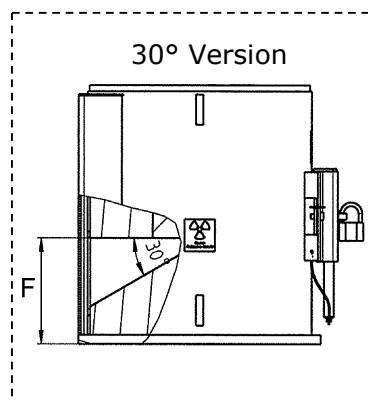
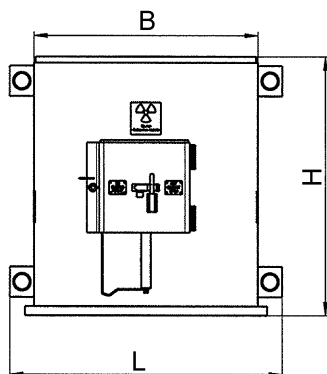
Contact Rating
Kontaktbelastbarkeit

Volt		Load / Last (A)		
AC	DC	R	L	Lamp / Lampe
250		15	3	1,5
125		15	3	1,5
	12	15	3	1,5
	24	10	2	1
	48	3	0,6	0,3
	250	0,25	0,05	0,025

LB 480

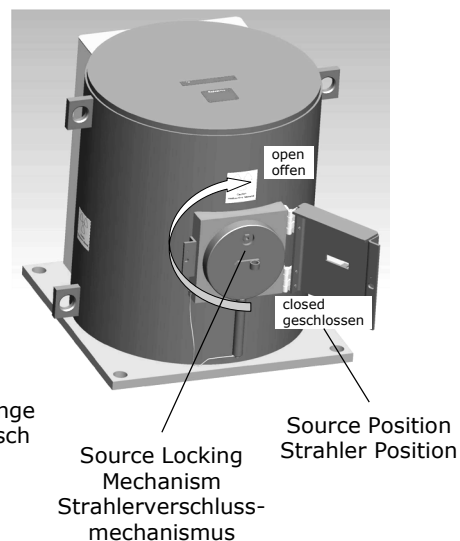
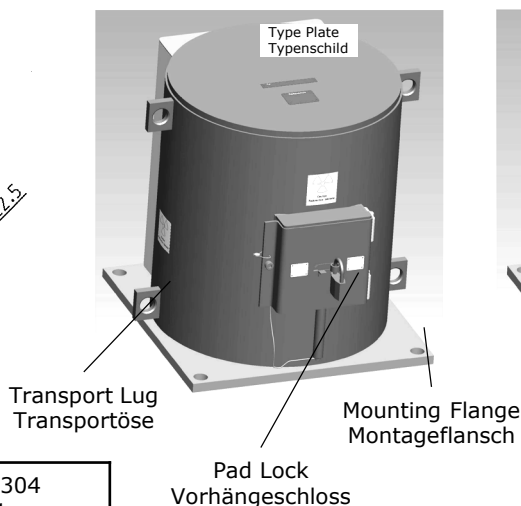
Dimensions in mm
Abmessungen in mm

9. Point Source Shielding LB 8030/8040 Punktstrahler-Abschirmbehälter LB 8030/8040



Shutter mechanism
locked
Strahlermechanismus
abgesperrt

Shutter mechanism
accessible
Strahlermechanismus
zugänglich

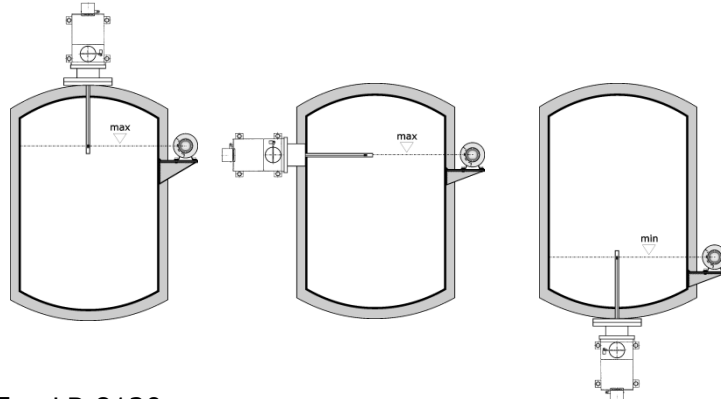


Material	Stainless Steel 304 Edelstahl 1.4301
Painting Lackierung	Polyurethane, yellow Polyurethan, gelb

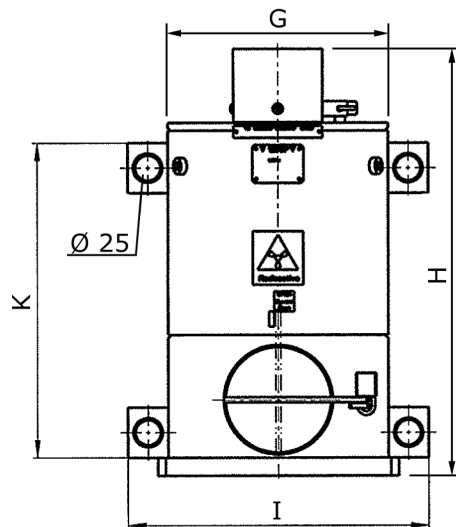
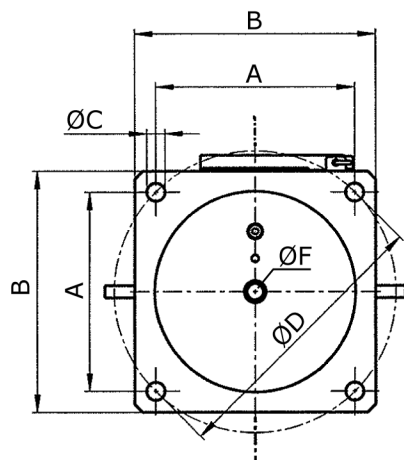
Type Typ	B	C	D	F	H	L		kg
LB 8030-01	370	360	400	174	425	450	30°	516
LB 8030-02							10°	
LB 8040-01	456	460	500	305	567	536	30°	1014
LB 8040-02							10°	

 Radiation Angle of the Shielding / Abstrahlwinkel der Abschirmung

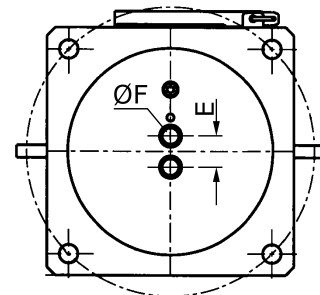
10. Point Source Flange Shielding LB 81xx for Dip Pipe Installations Punktstrahler-Flansch-Abschirmbehälter LB 81xx für Tauchrohr-Applikationen



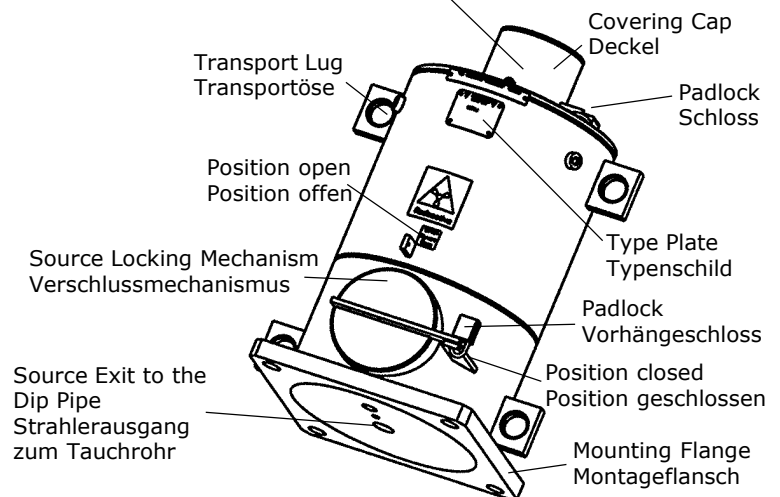
Type/Typ LB 8115 + LB 8120



Version for 2 Sources LB 81xx-20
Version für 2 Strahler LB 81xx-20



Junction Box for Shaft Core or Steel Cable
Anschlussraum für Wellenseele oder Seil

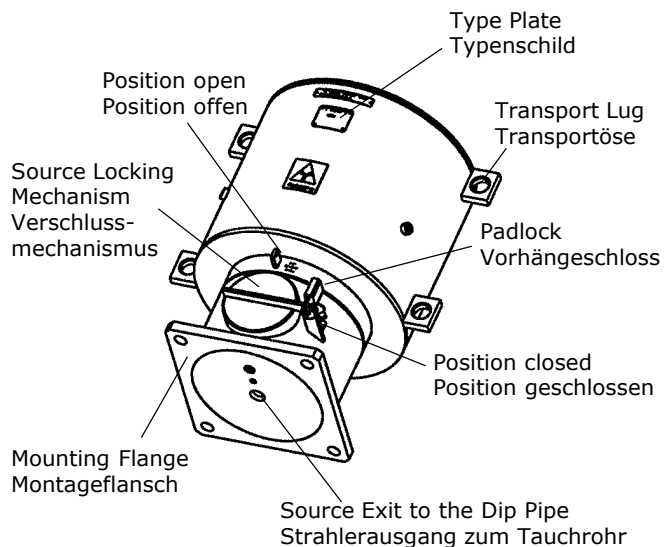
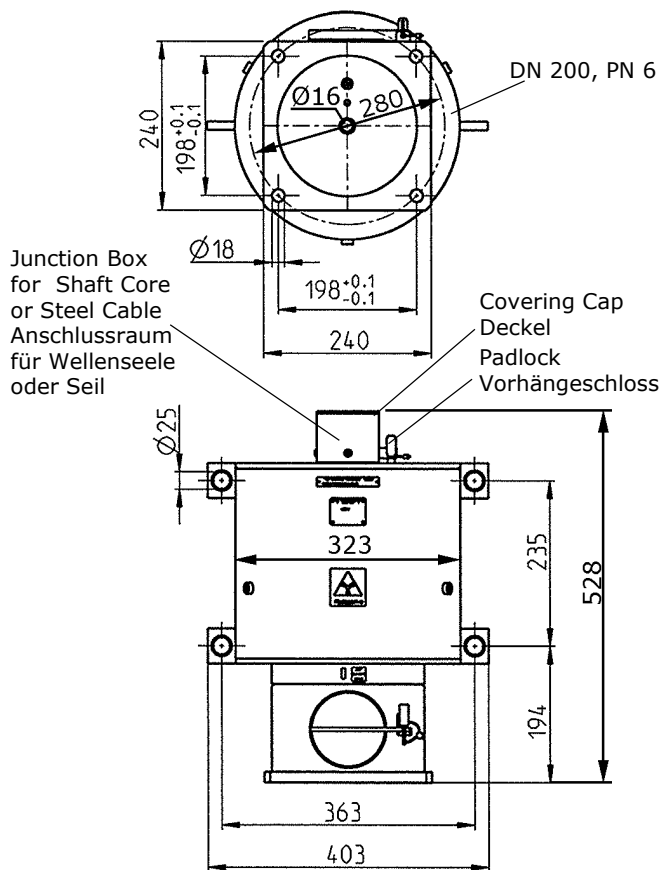


Type Typ	A	B	ØC	ØD	E	ØF	G	H	I	K	kg
LB 8115-01 LB 8115-20	141,5	180	18	200	- 30	16	159	350	239	237	60 kg
LB 8120-01 LB 8120-20	198	240	18	280	- 30	16	219	425 429	300	313	137 kg

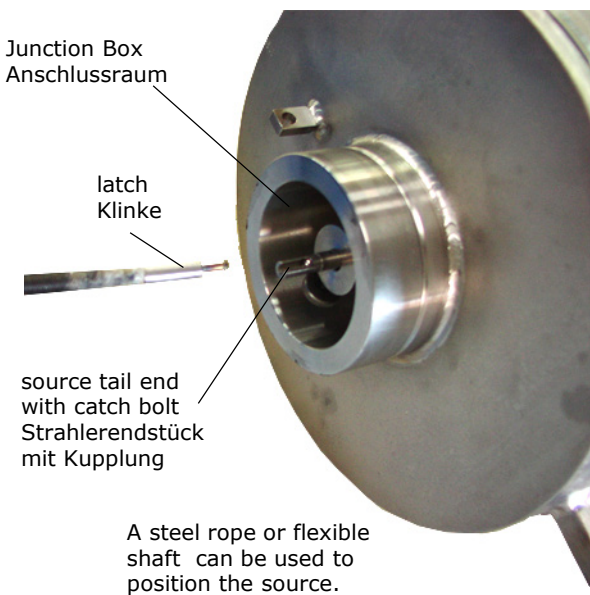
LB 480

Dimensions in mm
Abmessungen in mm

Type/Typ LB 8125-01

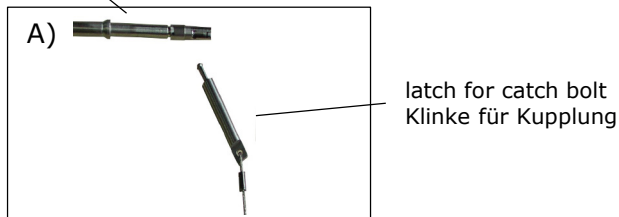
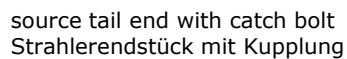


Operation of Flange Shielding Bedienung der Flanschabschirmung



A steel rope or flexible shaft can be used to position the source.

Zur Positionierung kann ein Stahlseil oder eine Wellenseele genutzt werden.



plug the latch into the catch bolt
Klinke in Kupplung einklinken



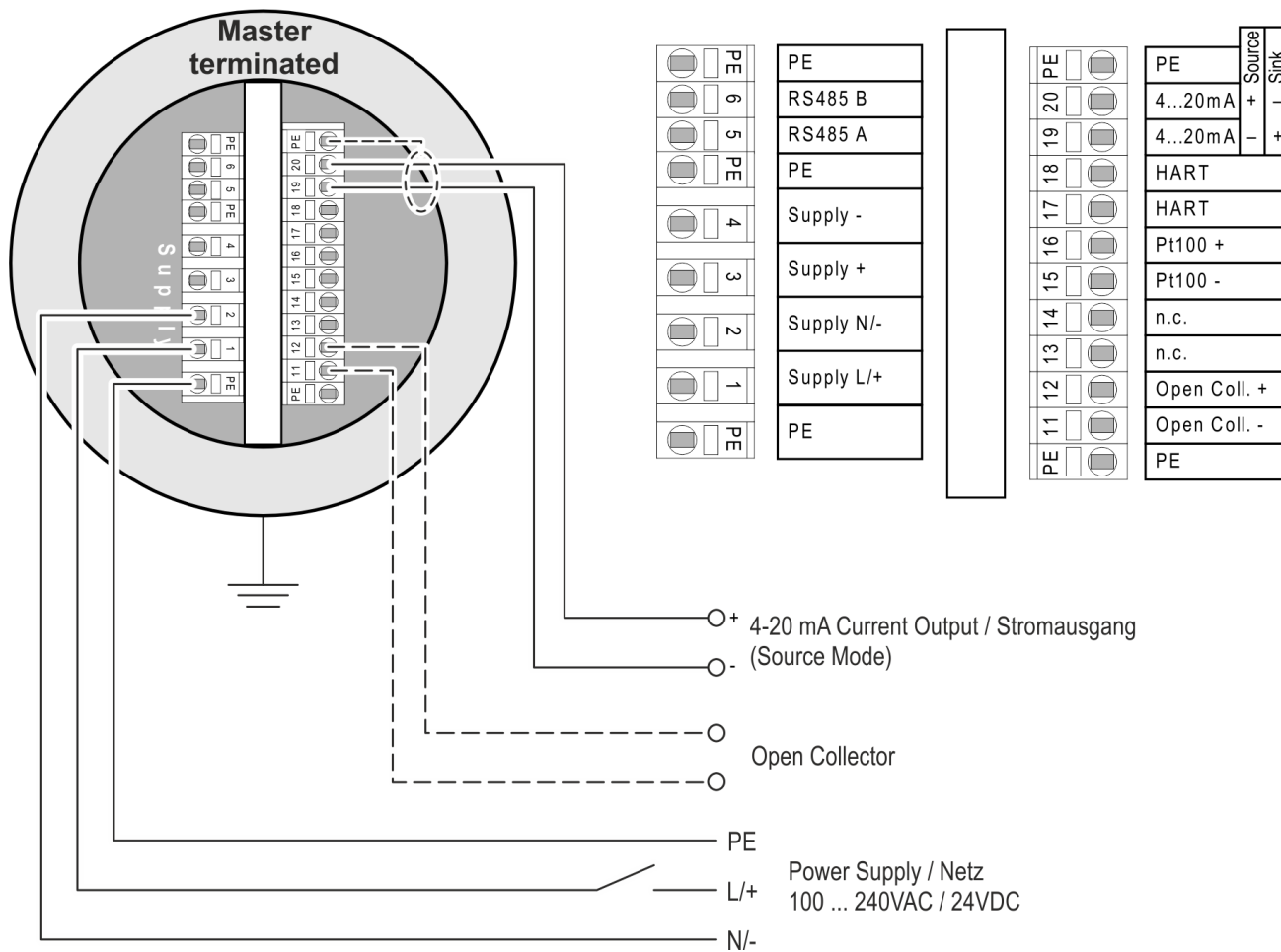
latch plugged into the catch bolt
Klinke in Kupplung einklinken



LB 480

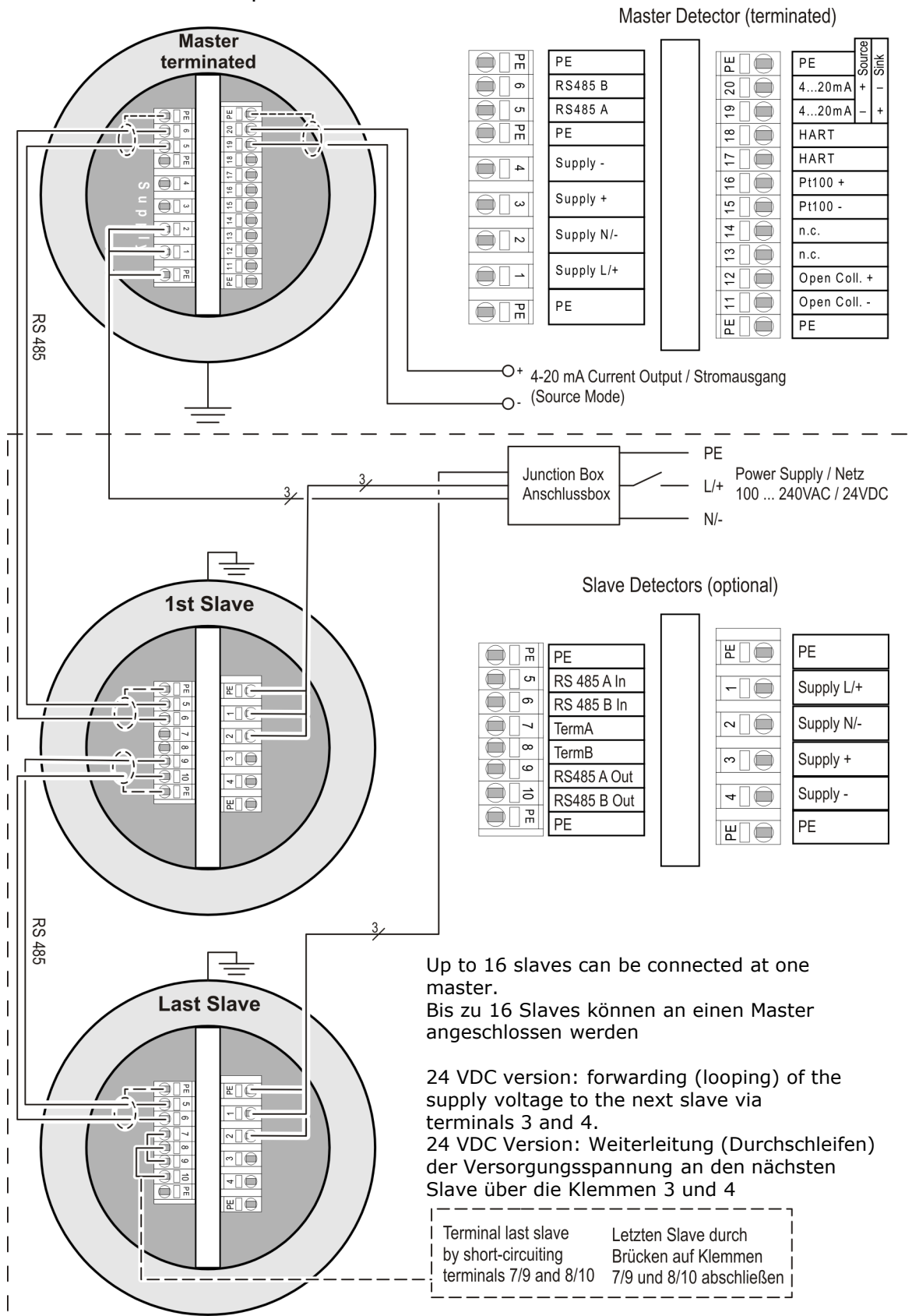
11. Terminal Connections Elektrische Anschlusspläne

11.1 Terminal Connections for Single Detector (terminated) Elektrische Anschlussplan für Einzeldetektor (terminiert)



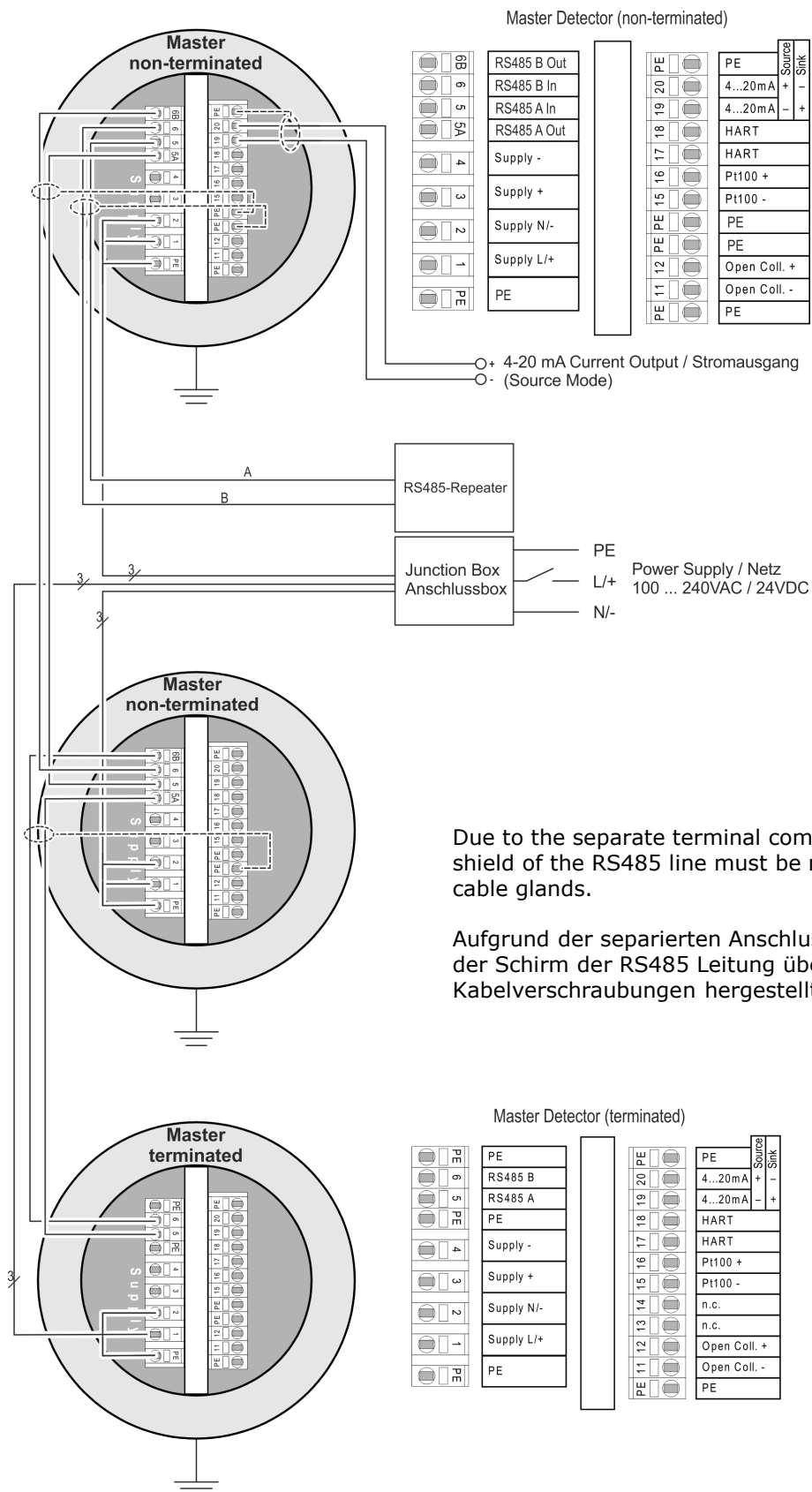
LB 480

11.2 Multi Detector Connection Diagram Master-Slave Detectors Multidetektor-Anschlussplan Master-Slave Detektoren



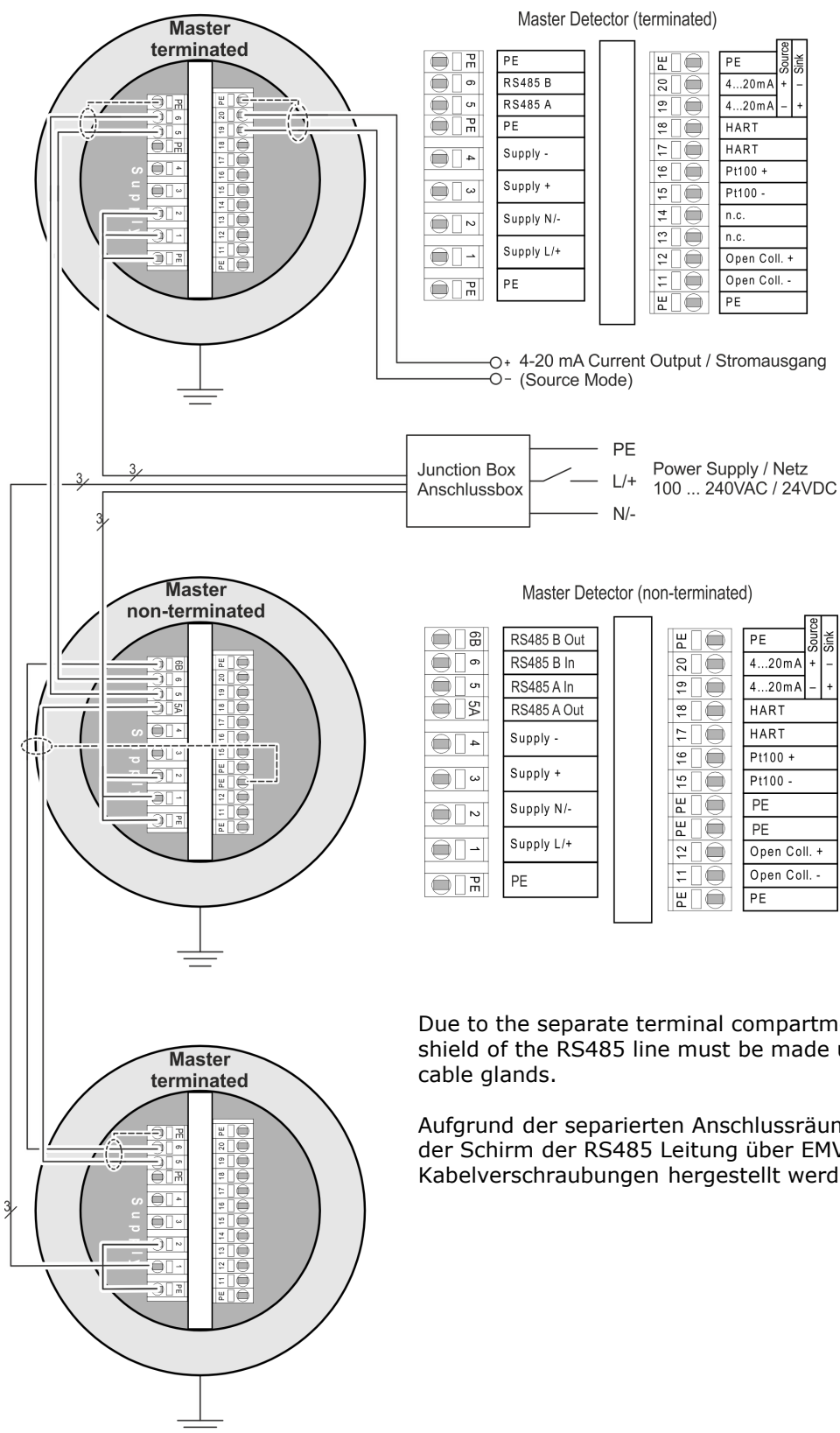
LB 480

11.3 Multi Detector Connection Diagram Master-Master Detectors at Repeater Module Multidetektor-Anschlussplan Master-Master Detektoren an Repeater Modul



LB 480

11.4 Multi Detector Connection Diagram Master-Master Detectors at cascaded Arrangement Multidetektor-Anschlussplan Master-Master Detektoren bei kaskadierter Anordnung



LB 480

[illegible]