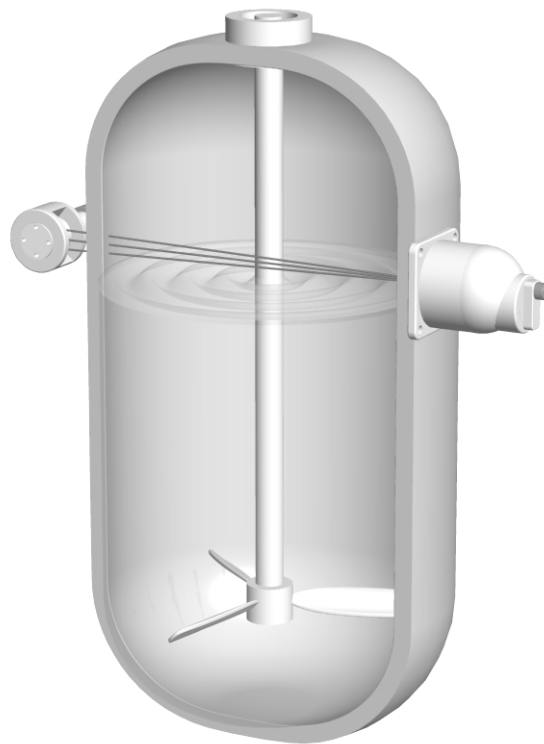


Technical Information

Level Switch LB 480

Level Switch
Grenzschalter

Field mounted components
Messstellen-Komponenten



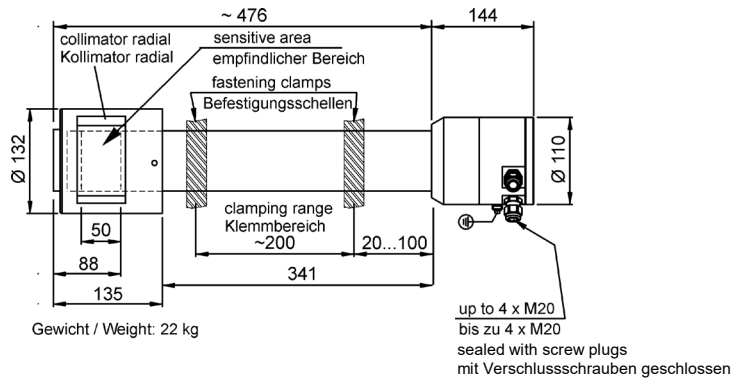
LB 480

Dimensions in mm
Abmessungen in mm

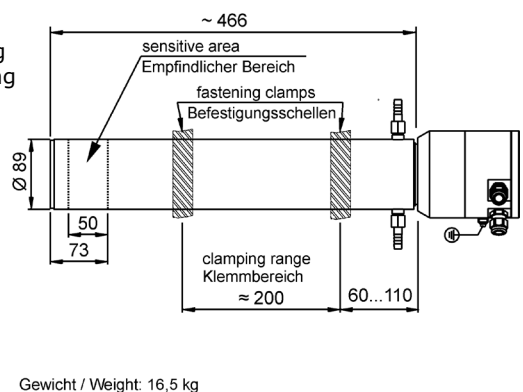
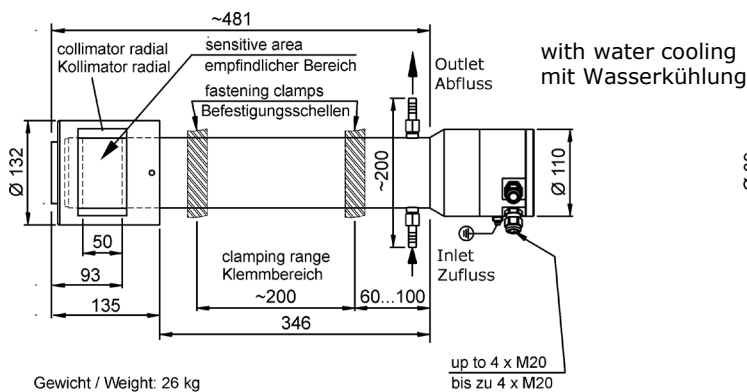
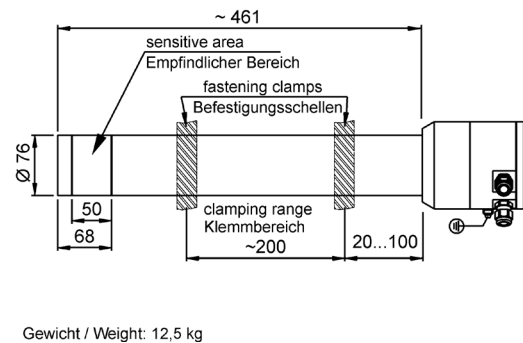
1.0 CrystaSENS (Version f. Zone 1/2)

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

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

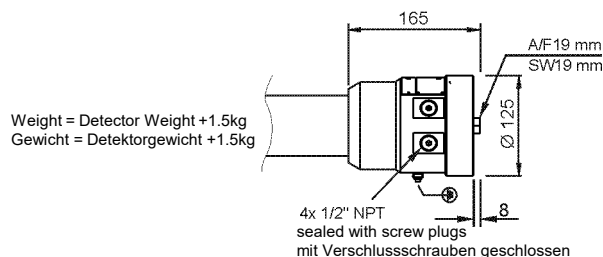


LB 480-1.-... -...-0.-



1.1 CrystaSENS (Version f. Divisions 1/2)

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



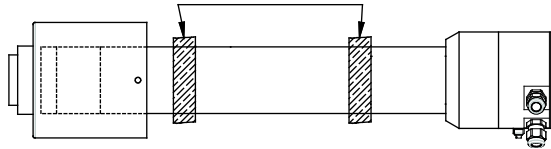
The detector version for divisions (NEC/CEC) differs only in the terminal housing, compared to the standard version illustrated above. The dimensions of this terminal housing are illustrated in this drawing.

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.

1.2 CrystaSENS Scintillator Size / Szintillatorgröße

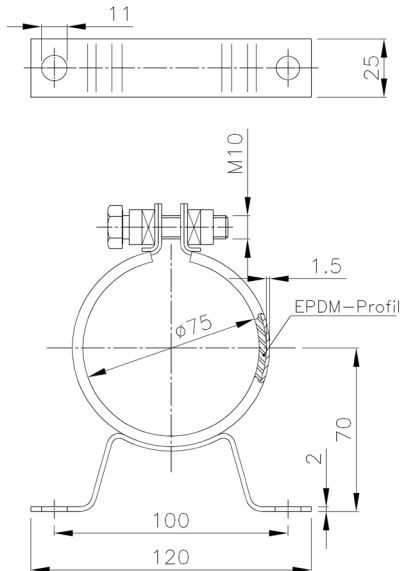
Type Typ	Scintillator Size Szintillatorgröße	($\varnothing$ /h)	Water cooling Wasserkühlung
LB 480-11	50/50		-
LB 480-12	50/50		✓
LB 480-13	40/35		-
LB 480-14	40/35		✓
LB 480-15	25/25		-
LB 480-16	25/25		✓

1.3 Mounting Clamps for Detector Befestigungsschellen für Detektor

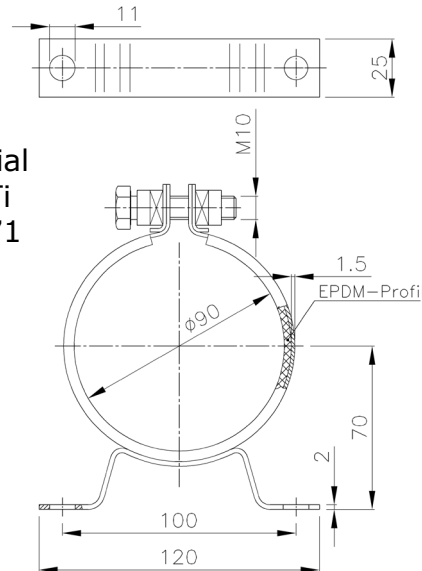


for Detectors without water cooling
für Detektoren ohne Wasserkühlung

for Detectors with water cooling
für Detektoren mit Wasserkühlung



Material
316Ti
1.4571



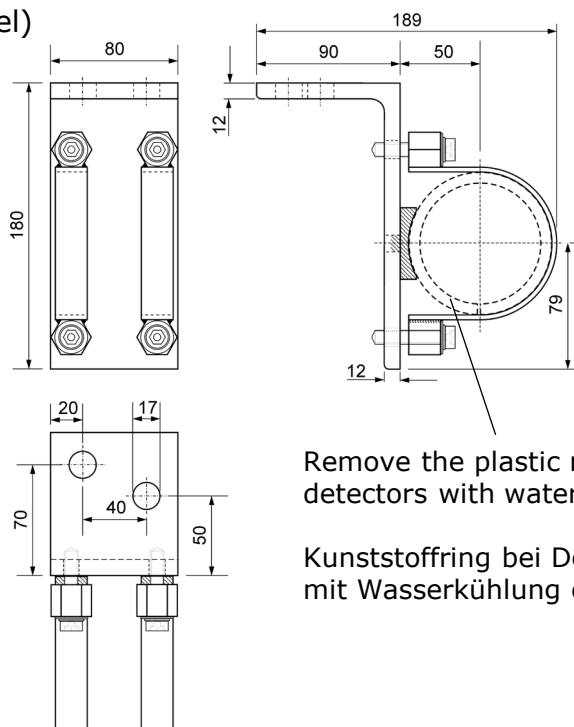
Id. Nr. 31346 (1 set = 2 clamps)
Id. Nr. 31345 (single clamp)

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

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)

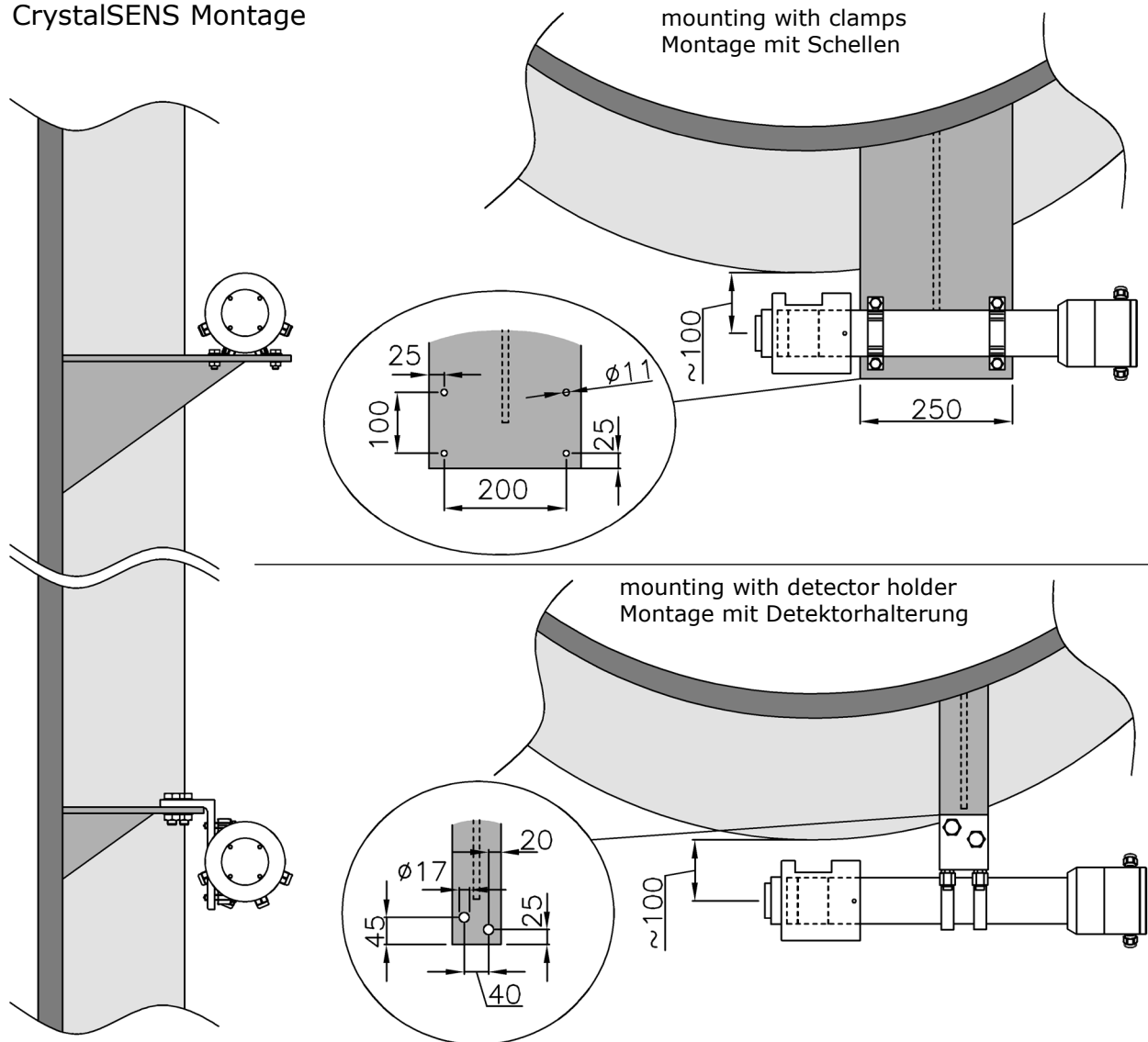
Part No. Id. Nr.	for Detector für Detektor
39246	without water cooling ohne Wasserkühlung
39247	with water cooling mit Wasserkühlung



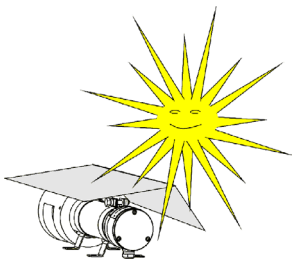
Remove the plastic ring for
detectors with water cooling.

Kunststoffring bei Detektoren
mit Wasserkühlung entfernen.

1.4 CrystaSENS Mounting CrystaSENS Montage



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.

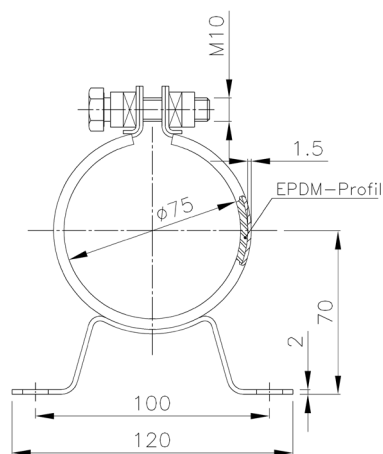
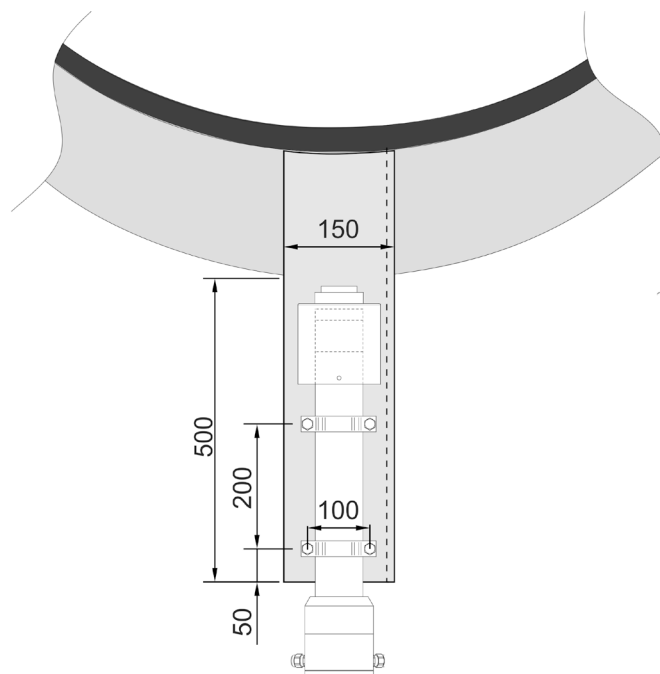
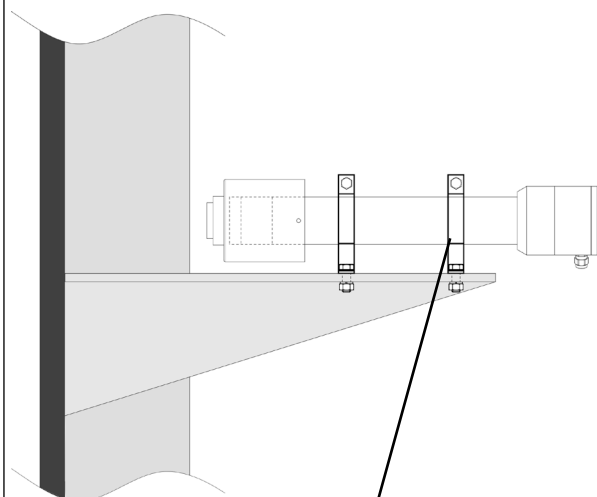
Wird durch Sonneneinstrahlung eine Detektortemperatur von über 50°C erreicht, so ist ein geeigneter Sonnenschutz zu montieren. Auch die Aufheizung des Detektors durch Wärmeabstrahlung vom Behälter kann durch ein dünnes Wärmeableitblech gemildert werden. Für jeden Detektor steht auch eine geeignete Wasserkühlung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

LB 480

Dimensions in mm
Abmessungen in mm

1.5 CrystaSENS Mounting Axial Collimator CrystaSENS Montage axialer Kollimator

mounting with clamps
Montage mit Schellen

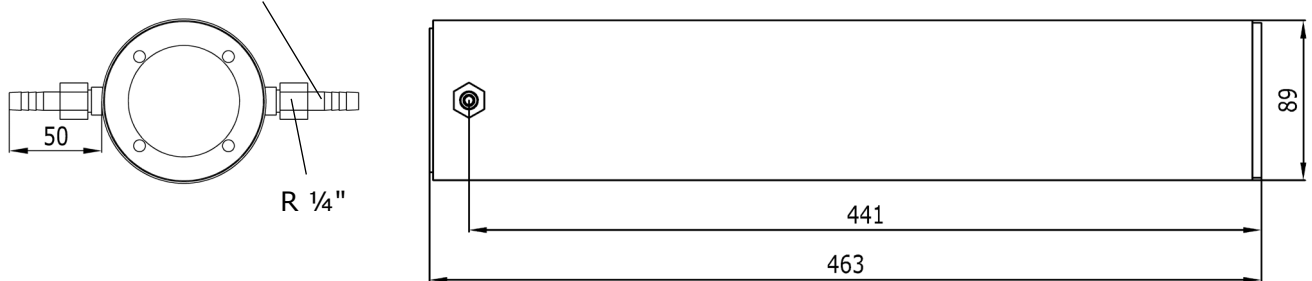


Id. Nr. 31346 (1 set = 2 clamps)

Id. Nr. 31345 (single clamp)

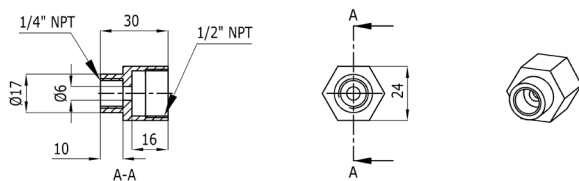
1.6 CrystaSENS Water Cooling Jacket and Adaptor Fittings CrystaSENS Wasserkühlung und Adapter Anschlussstücke

pipe connection diameter 10
Schlauchanschluss Ø 10



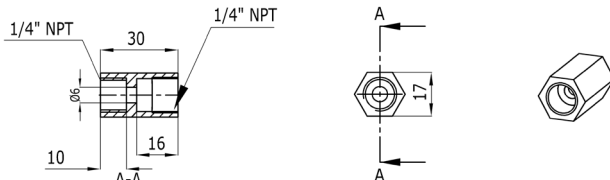
Fitting adaptor for standard water cooling Rp 1/4" → 1/2" NPT
stainless steel 304, part no: 47189

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



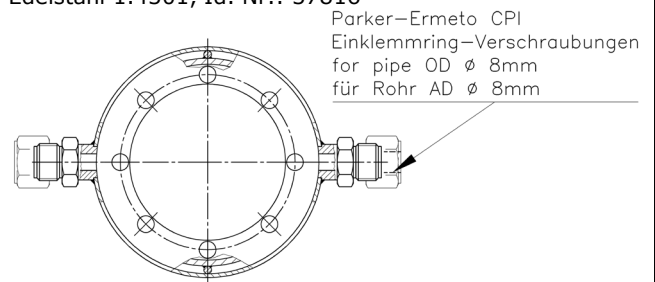
Fitting adaptor for standard water cooling Rp 1/4" > 1/4" NPT
stainless steel 304, part no: 46743

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



Water cooling jacket with Parker Ermeto Fittings
stainless steel 304, part no: 37816

Wasserkühlung mit Parker Ermeto Anschlüssen
Edelstahl 1.4301, Id. Nr.: 37816



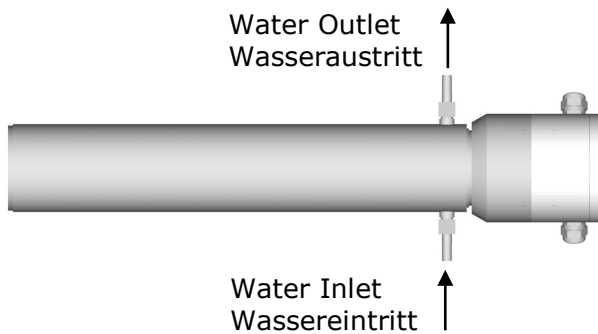
Further fitting adaptors for standard water cooling jacket:
Rp 1/4" > 1/2" NPT male, stainless steel 304, part no: 06352
Rp 1/4" > 1/4" NPT male, stainless steel 304, part no: 06349
Weitere Adapter für die Standard-Wasserkühlung:
Rp 1/4" > 1/2" NPT Außengewinde, 1.4301, Id. Nr.: 06352
Rp 1/4" > 1/4" 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)

1.7 Water Cooling Installation Instruction Anweisung zur Installation der Wasserkühlung

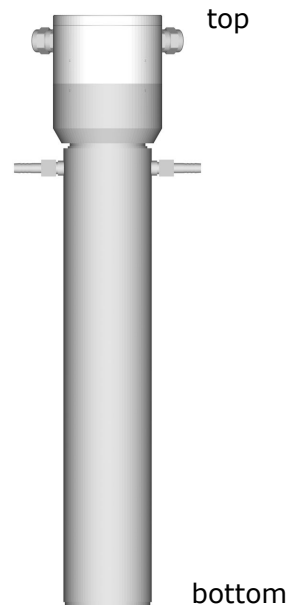
Horizontal Detector Installation Horizontale Detektor Installation



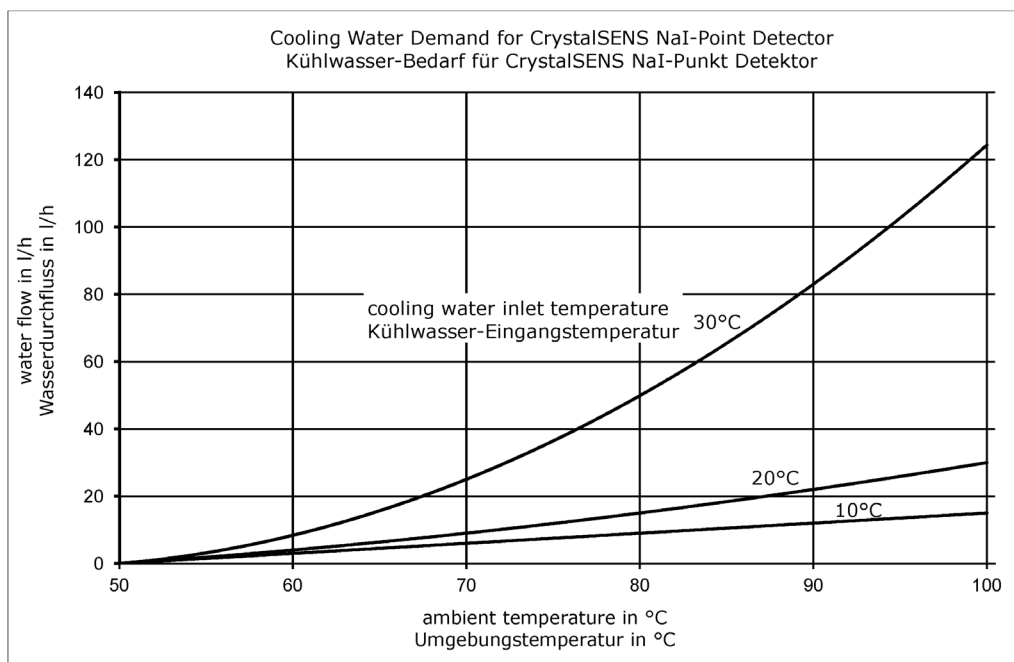
In order to fill the entire water cooling jacket, incoming water must enter from the bottom. Damit sich die Wasserkühlung vollständig mit Wasser füllt, muss der Wasserzufluss von unten erfolgen.

Vertical Detector Installation Vertikale Detektor Installation

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



1.8 Detector Cooling Water Demand Detektor Kühlwasserbedarf

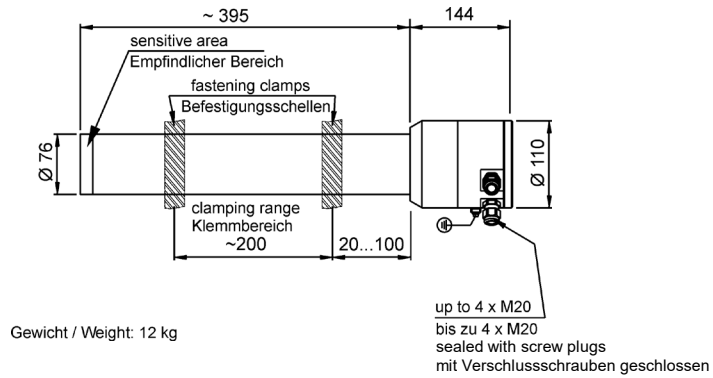


LB 480

Dimensions in mm
Abmessungen in mm

1.8 CrystalSENS 44/5 f. Am-241
(none Ex, nicht-Ex)

LB 480-17-00 -..-0.-
LB 480-17-Z0 -..-0.-



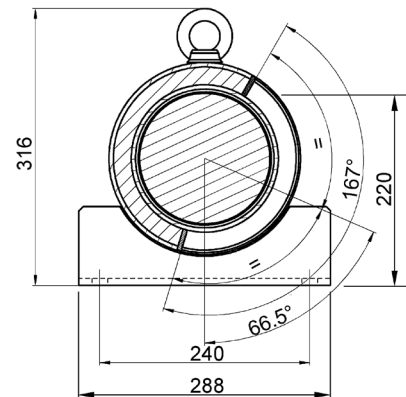
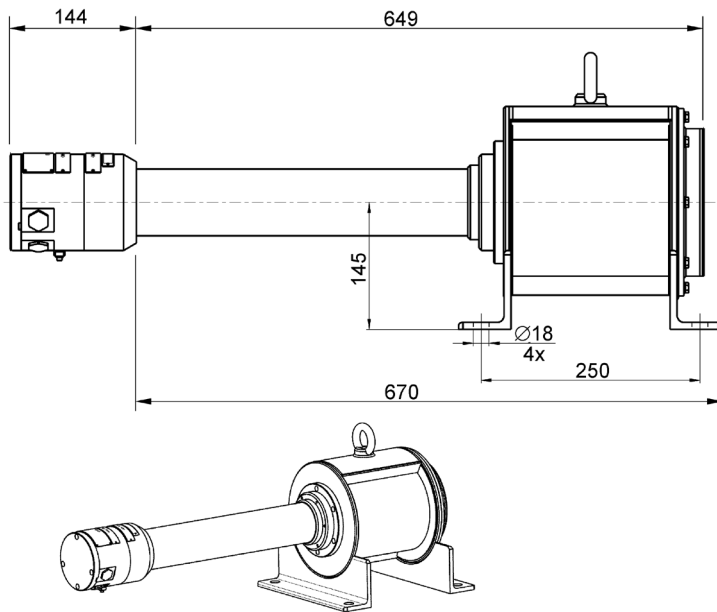
LB 480

Dimensions in mm
Abmessungen in mm

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

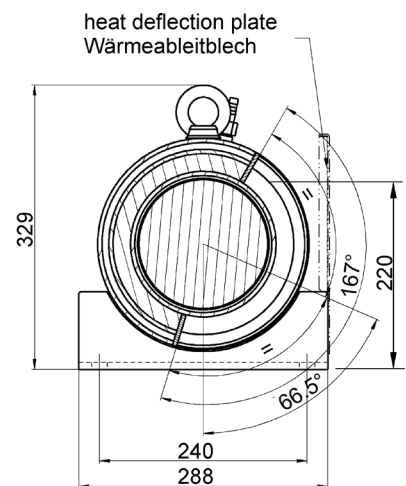
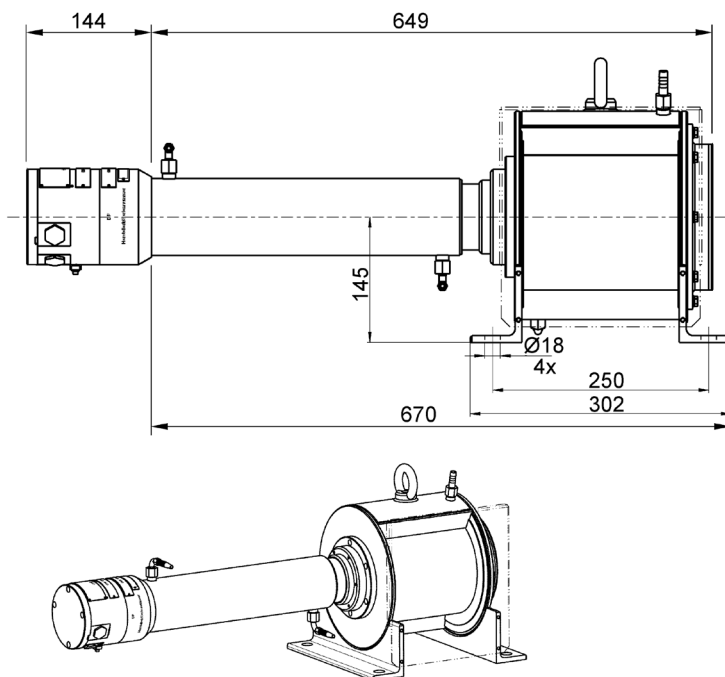
LB 480-3.-0.
LB 480-3.-1.
LB 480-3.-2.
LB 480-3.-Z.

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



Weight/Gewicht 52 kg

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

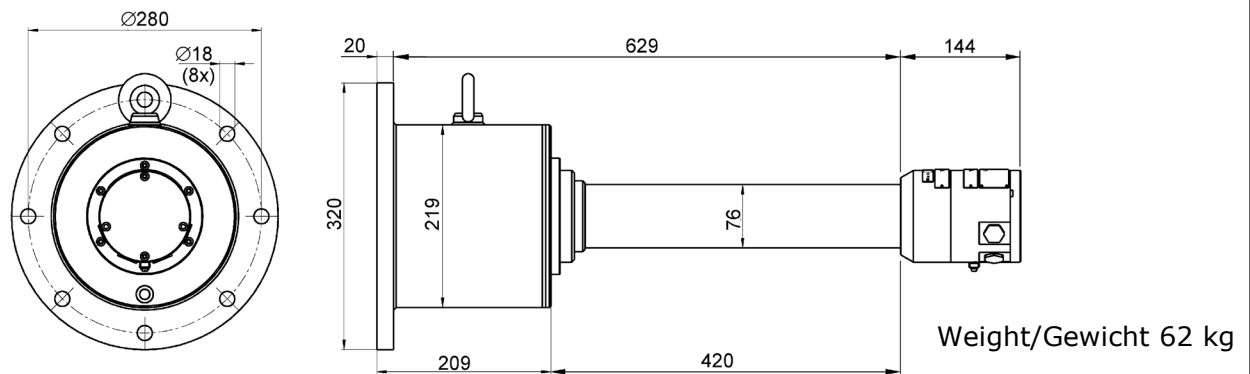


Weight/Gewicht 59 kg

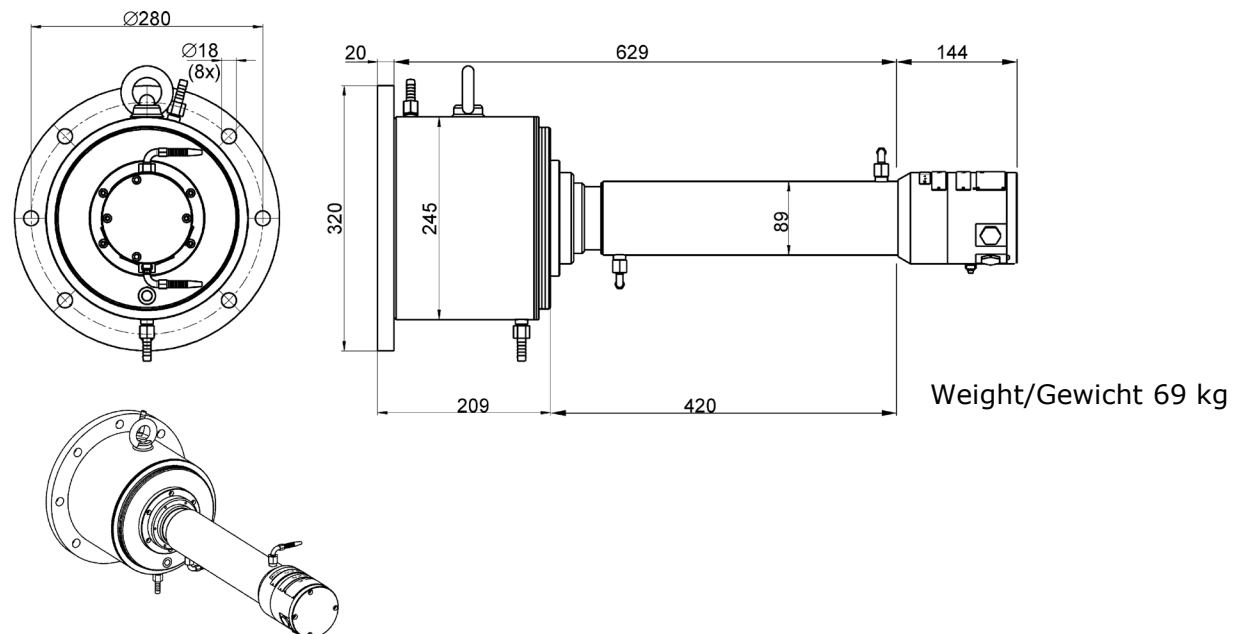
LB 480

Dimensions in mm
Abmessungen in mm

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

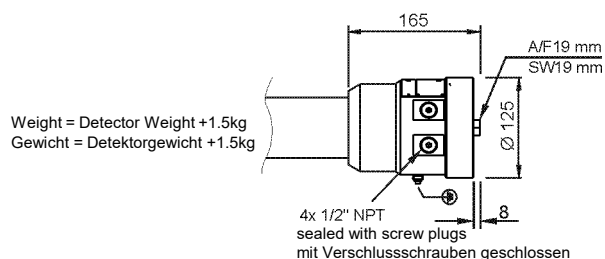


2.4 with Axial Irradiation and Water Cooling mit frontaler Einstrahlung und Wasserkühlung LB 480-32-...-a.



2.5 SuperSENS (Version f. Divisions 1 + 2)

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



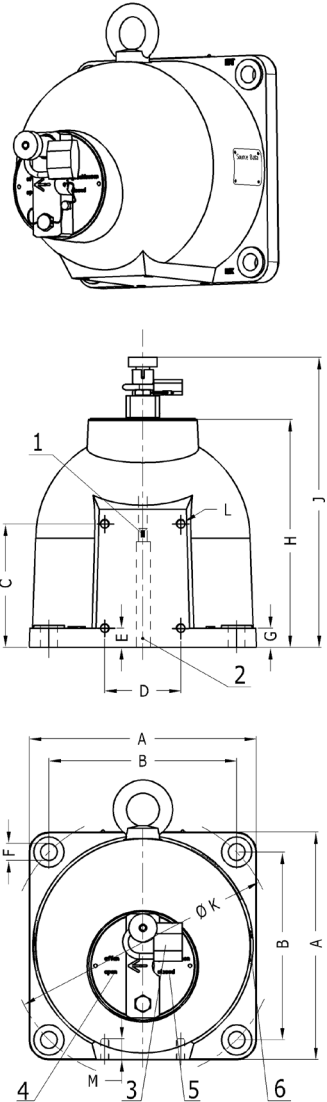
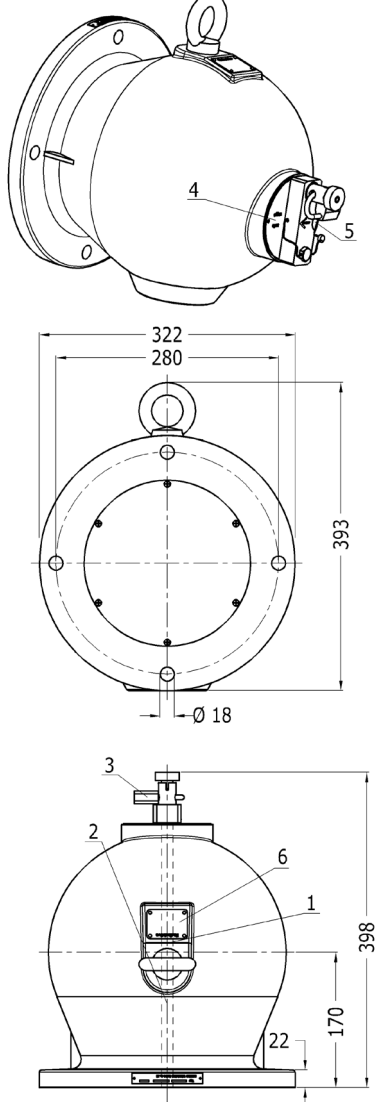
The detector version for divisions (NEC/CEC) differs only in the terminal housing, compared to the standard version illustrated above. The dimensions of this terminal housing are illustrated in this drawing.

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

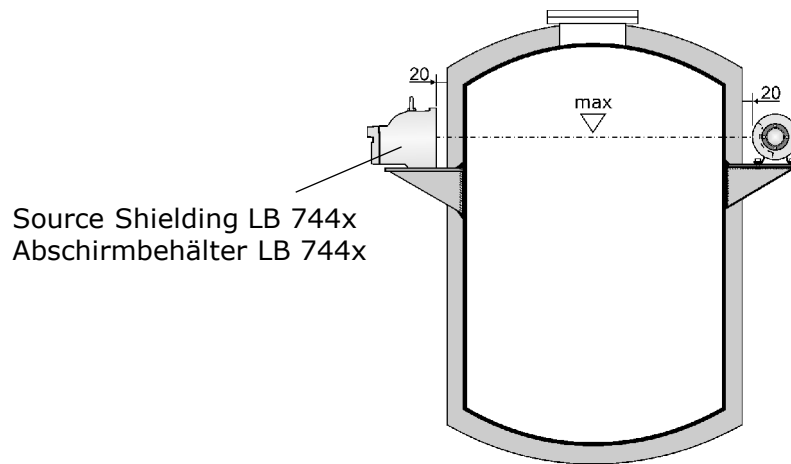
3.0 Point Source Shielding LB 744x Punktstrahler-Abschirmbehälter LB 744x

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

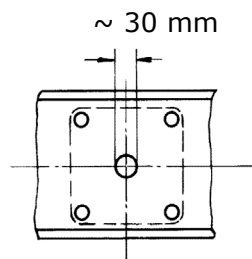
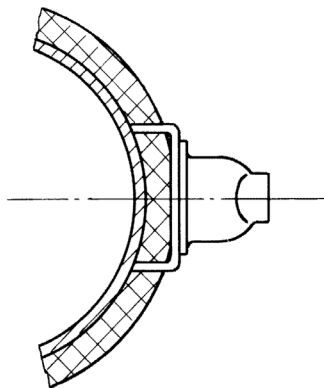
Housing Gehäuse	Type Typ	A	B	C	D	E	FØ	G	H	J	KØ	L	M		kg
Carbon Steel St 37	LB 7440-F-CR LB 7445-F-CR	180	142	75	60	15	18	20	173	238	200	M8	12	16°	31
Super Duplex UNS 32750 SAF 2507 1.4410	LB 7440-FE-CR LB 7445-FE-CR														
	LB 7442-FE-CR	240	198	130	80	20	18	20	242	306	280	M10	14	9°	81
Carbon Steel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

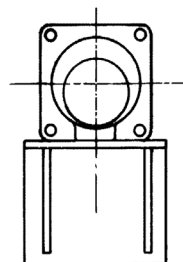
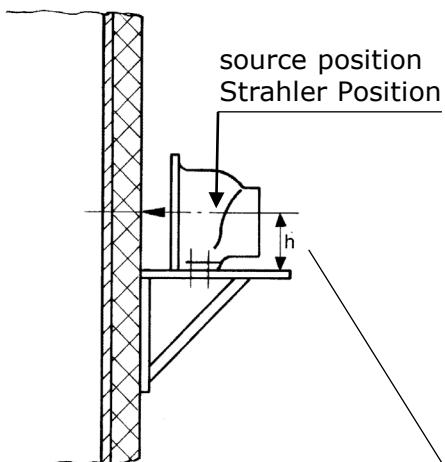
3.1 Mounting Proposal for Source Shielding LB 744x Montagevorschlag für Abschirmbehälter LB 744x



Source Shielding LB 744x
Abschirmbehälter LB 744x



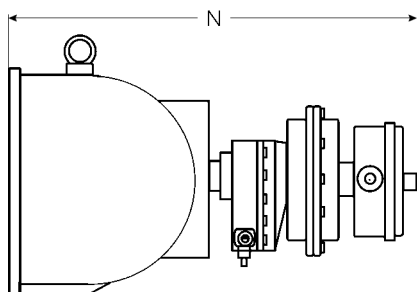
Flange Installation
Flanschinstallation



Pedestal Installation
Sockelinstallation

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

3.2 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 Pneumatischer Verschlussantrieb mit Endschalter
Ex de IIC T6	80919	Pneumatic Actuator with Limit Switch with ATEX Pneumatischer Verschlussantrieb mit Endschalter with ATEX

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 for intrinsically safe power supply 2 Näherungssensoren für Eigensichere Speisung

3.3 Components for Pneumatic Actuator Einzelteile für Pneumatischen Verschlussantrieb

Pneumatic Actuator
Pneumatischer Antrieb

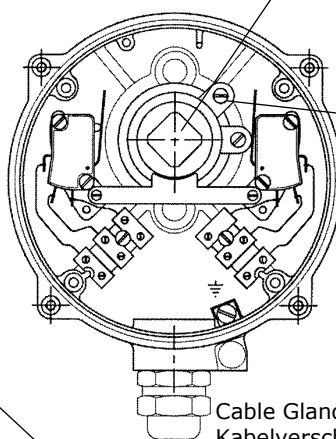
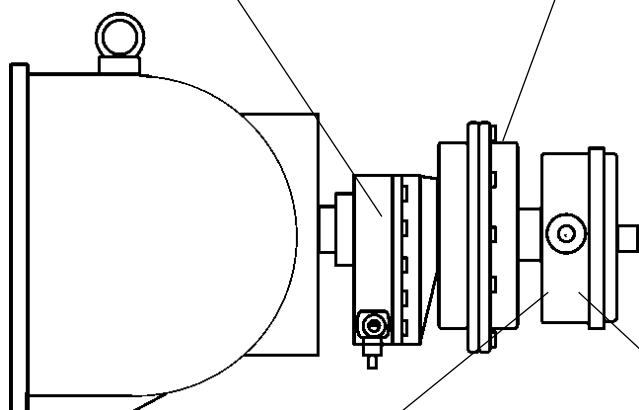


Return Spring
Rückstellfeder



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

Außenvierkant zur Stel-
lungsanzeige, oder manueller
Betätigung



adjustable cam
verstellbare Nocken

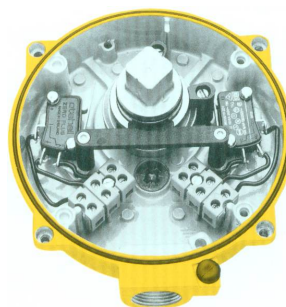
Cable Gland
Kabelverschraubung

Limit Switch in Ex de
Endschaltereinheit in Ex de



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

Limit Switch in IP 65
Endschaltereinheit 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
	30	7	5
	75	1	1
	125	0,5	0,06
	250	0,25	0,03

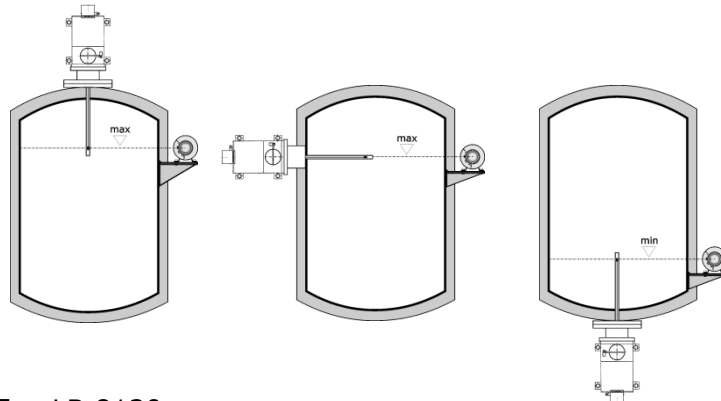
Contact Rating
Kontaktbelastbarkeit

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

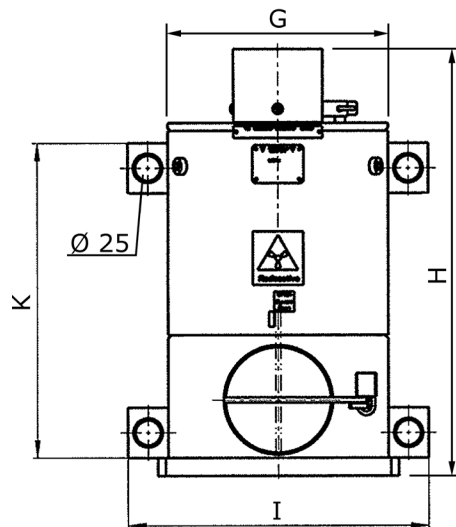
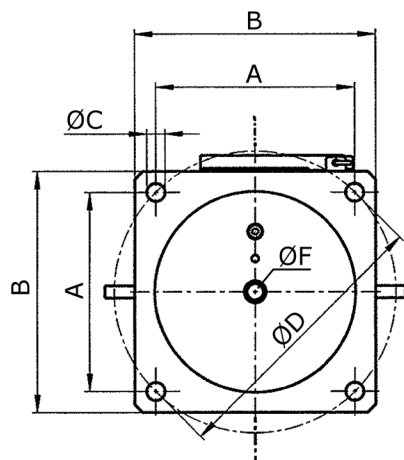
LB 480

Dimensions in mm
Abmessungen in mm

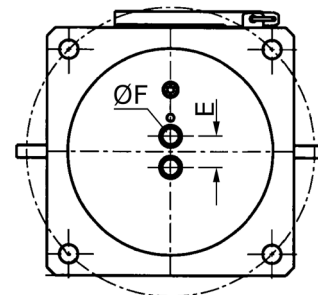
4.0 Point Source Flange Shielding LB 81xx for Dip Pipe Installations Punktstrahler-Flansch-Abschirmbehälter LB 81xx für Tauchrohr-Applikationen



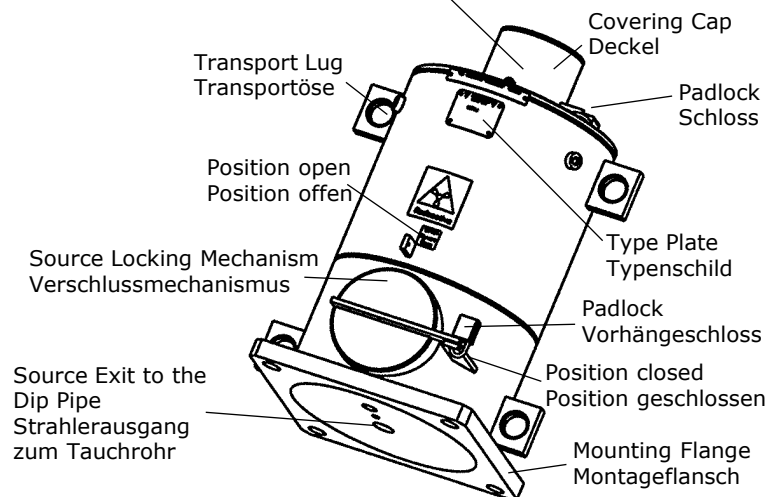
Type/Typ LB 8115 + LB 8120



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

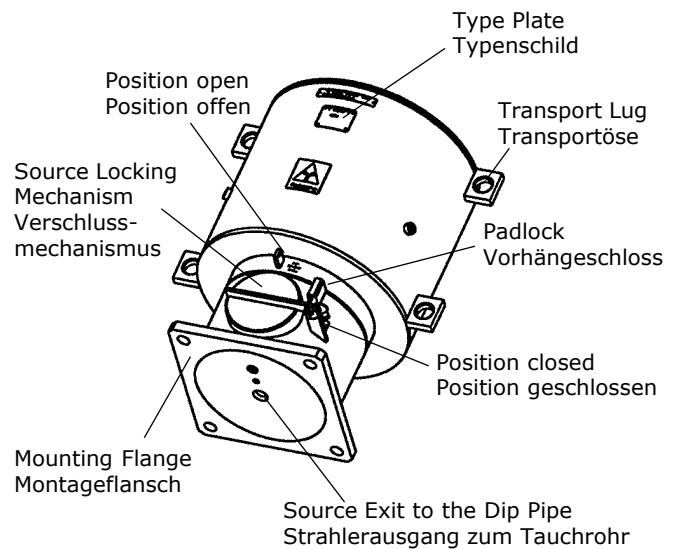
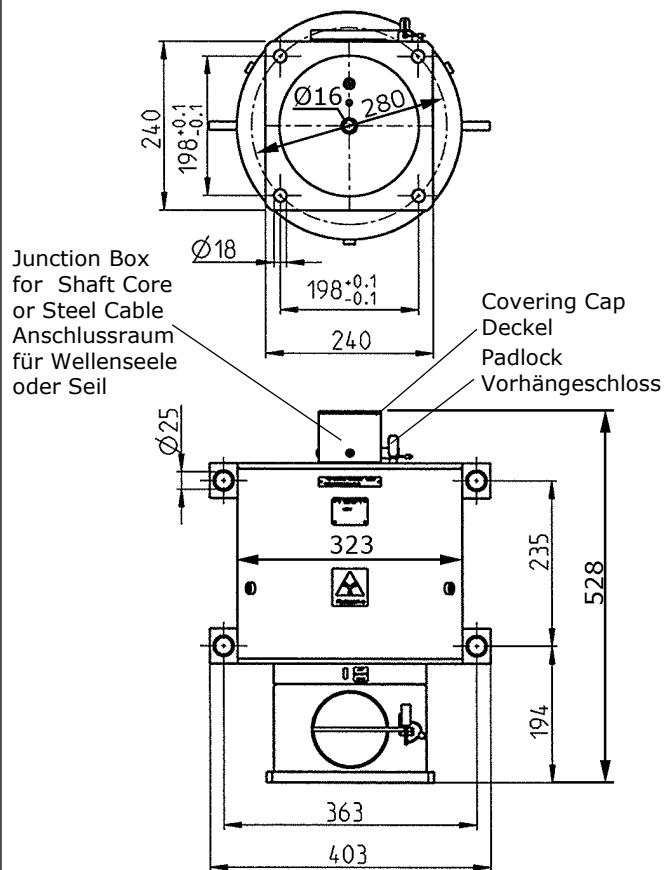


Type Typ	A	B	ØC	ØD	E	ØF	G	H	I	K	kg
LB 8115-01 LB 8115-20	141,5	180	18	200	- 30	16	159	350	239	237	60 kg
LB 8120-01 LB 8120-20	198	240	18	280	- 30	16	219	425 429	300	313	137 kg

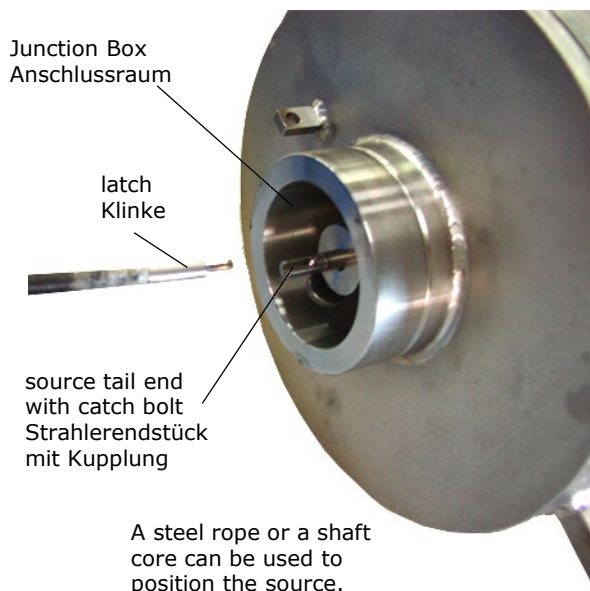
LB 480

Dimensions in mm
Abmessungen in mm

Type/Typ LB 8125



4.1 Operation of Flange Shielding Bedienung der Flanschabschirmung



Zur Positionierung kann ein Stahlseil oder eine Wellenseele genutzt werden.

source tail end with catch bolt
Strahlerendstück mit Kupplung

A)

latch for catch bolt
Klinke für Kupplung

plug the latch into the catch bolt
Klinke in Kupplung einklinken

B)

latch plugged into the catch bolt
Klinke in Kupplung einklinken

C)

LB 480

Dimensions in mm
Abmessungen in mm

5.0 Terminal Connection Elektrischer Anschlussplan

