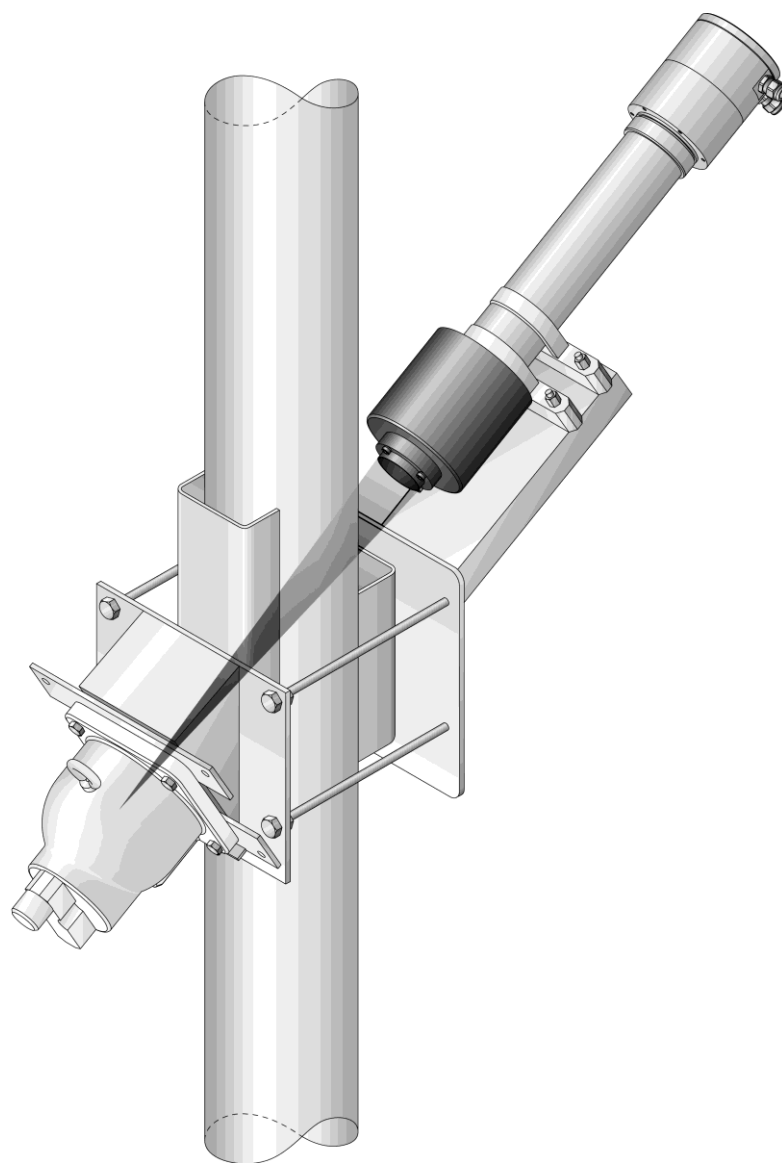


Technical Information

Density LB 480

Density Gauge
Dichte Messung

Field mounted components
Messstellen-Komponenten

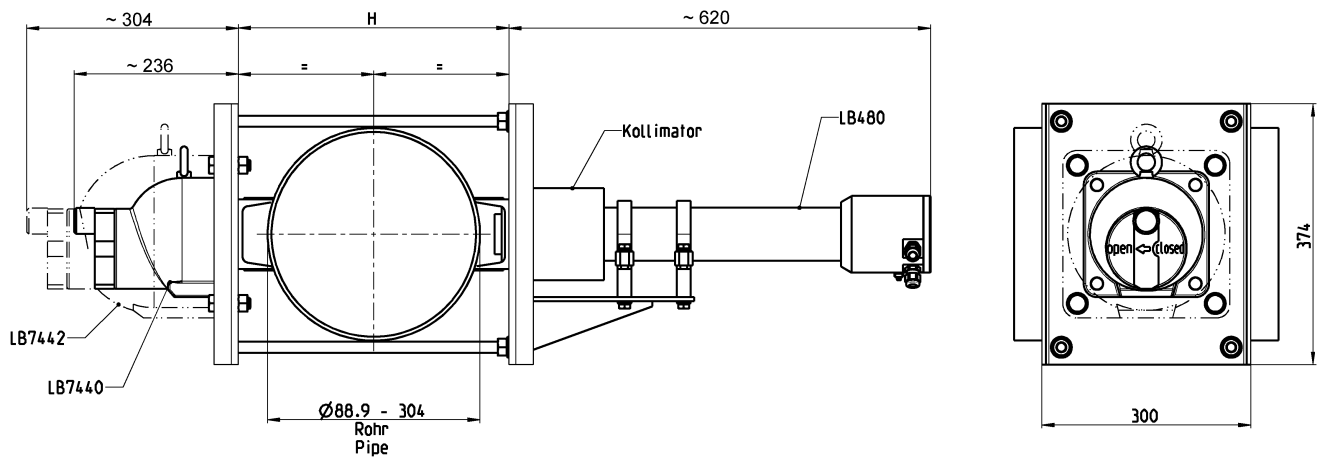
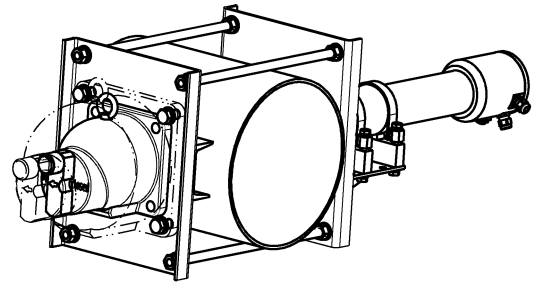


LB 480

Dimensions in mm
Abmessungen in mm

1.0 Clamping Device 90° for Pipe Diameter 88.9 ... 304 mm
Montagevorrichtung 90° für Rohrdurchmesser 88.9 ... 304 mm

Material	Carbon Steel Stahl St37
Painting Lackierung	Polyurethane, gray Polyurethan, grau



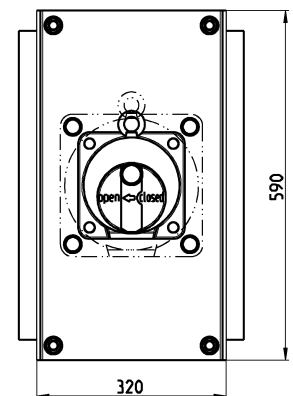
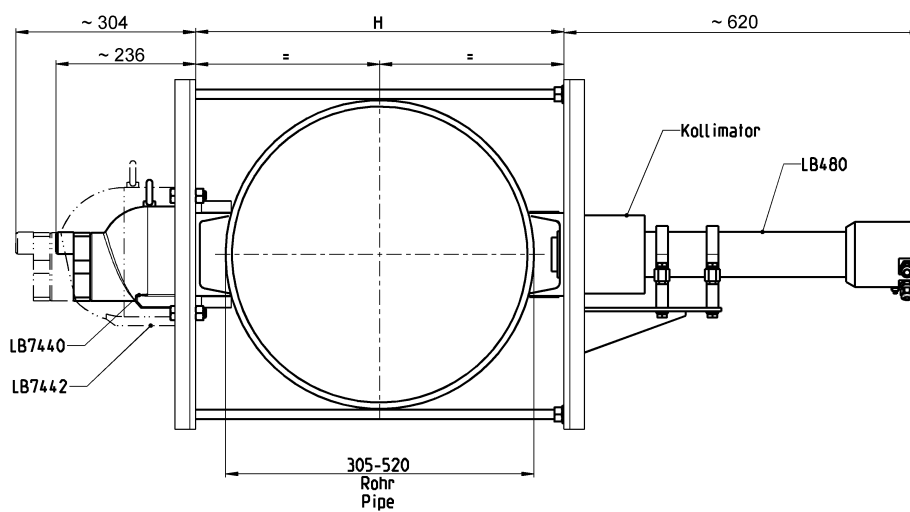
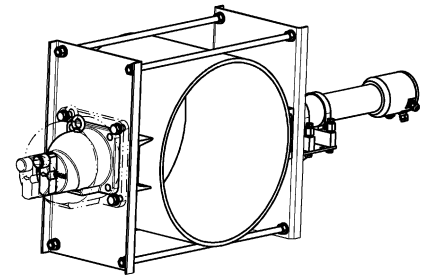
Part No. Id. Nr.	Pipe Diameter Rohrdurchmesser	H	Weight of Clamp. Device Gewicht der Montagevorrichtung
80795	88.9	120	23 kg
	101.6	156	
	114.3	180	
	141.3	218	
	168.3	250	
	219.1	310	
	273.0	368	
	304.0	402	

LB 480

Dimensions in mm
Abmessungen in mm

1.1 Clamping Device 90° for Pipe Diameter 305 ... 521 mm Montagevorrichtung 90° für Rohrdurchmesser 305 ... 521 mm

Material	Carbon Steel Stahl St37
Painting Lackierung	Polyurethane, gray Polyurethan, grau



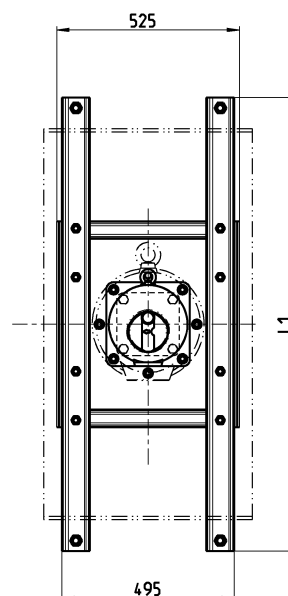
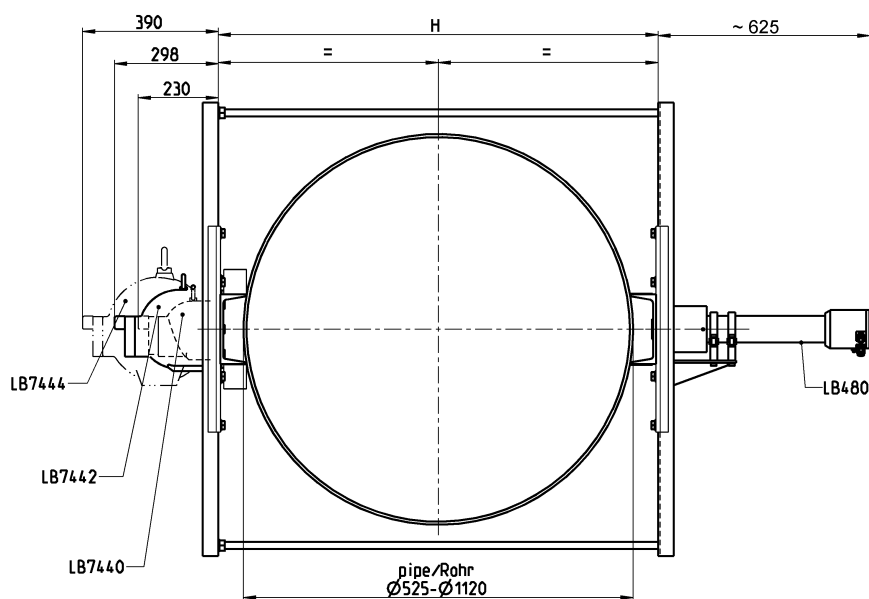
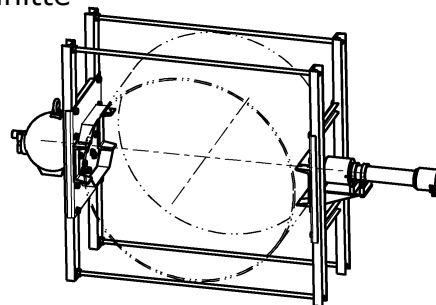
Part No. Id. Nr.	Pipe Diameter Rohrdurchmesser	H	Weight of Clamp. Device Gewicht der Montagevorrichtung
80796	305.0	400	34 kg
	318.0	413	
	323.8	419	
	355.6	451	
	406.4	501	
	457.2	552	
	508.0	603	
	521.0	616	

LB 480

Dimensions in mm
Abmessungen in mm

1.2 Clamping Device 90° for Large Pipe Diameters Montagevorrichtung 90° für große Rohrleitungsquerschnitte

Material	Carbon Steel Stahl St37
Painting Lackierung	Polyurethane, gray Polyurethan, grau



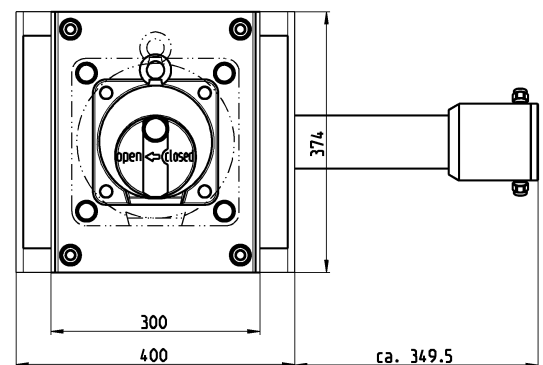
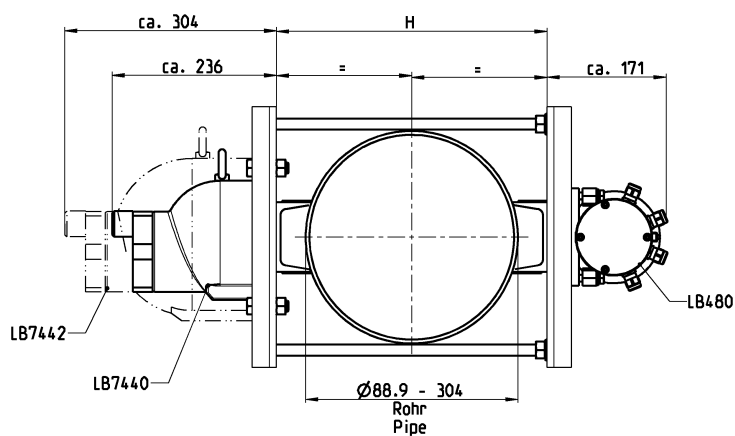
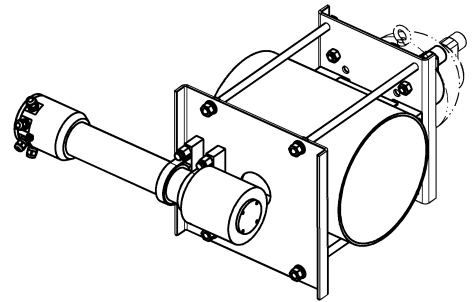
Part No. Id. Nr.	Pipe Ø Range Rohr Ø Bereich	H	Length of the thread bars Länge der Gewindestangen	L1	Weight of Clamp. Device Gewicht der Montagevorrichtung
51872-01	400 ... 519	643	735	680	76 kg
51872-02	520 ... 559	686	790	680	77 kg
51872-03	560 ... 659	792	890	740	79 kg
51872-04	660 ... 759	896	990	840	84 kg
51872-05	760 ... 869	1009	1100	950	89 kg
51872-06	870 ... 1020	1163	1250	1200	98 kg
51872-07	1020 ... 1120	1265	1350	1300	103 kg

LB 480

Dimensions in mm
Abmessungen in mm

2.0 Parallel Clamping Device 90° for Pipe Diameter 88.9 ... 304 mm
Parallele Montagevorrichtung 90° für Rohrdurchmesser 88.9 ... 304 mm

Material	Carbon Steel Stahl St37
Painting Lackierung	Polyurethane, gray Polyurethan, grau



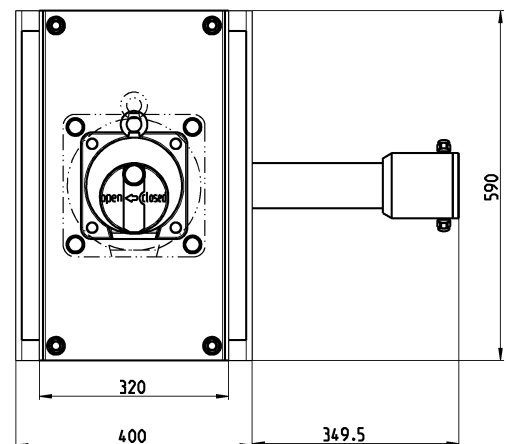
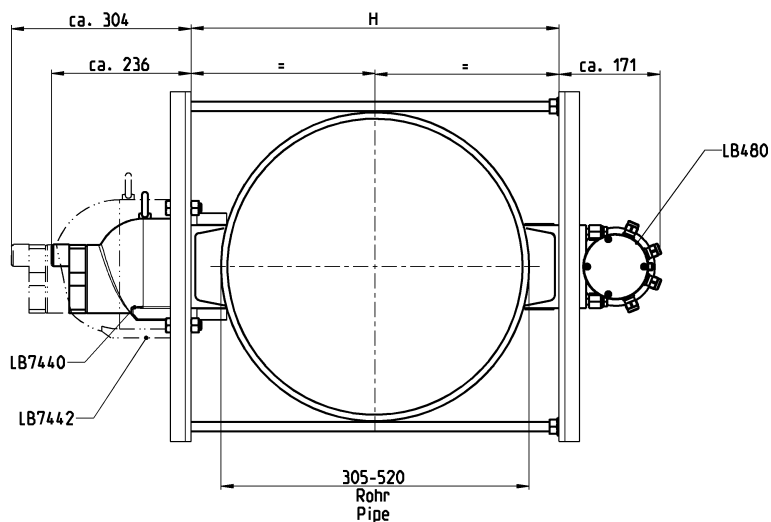
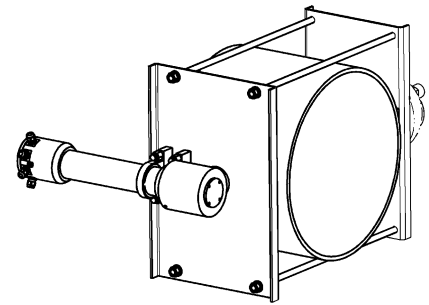
Part No. Id. Nr.	Pipe Diameter Rohrdurchmesser	H	Weight of Clamp. Device Gewicht der Montagevorrichtung
59296	88.9	120	22 kg
	101.6	156	
	114.3	180	
	141.3	218	
	168.3	250	
	219.1	310	
	273.0	368	
	304.0	402	

LB 480

Dimensions in mm
Abmessungen in mm

2.1 Parallel Clamping Device 90° for Pipe Diameter 305 ... 521 mm Parallele Montagevorrichtung 90° für Rohrdurchmesser 305 ... 521 mm

Material	Carbon Steel Stahl St37
Painting Lackierung	Polyurethane, gray Polyurethan, grau



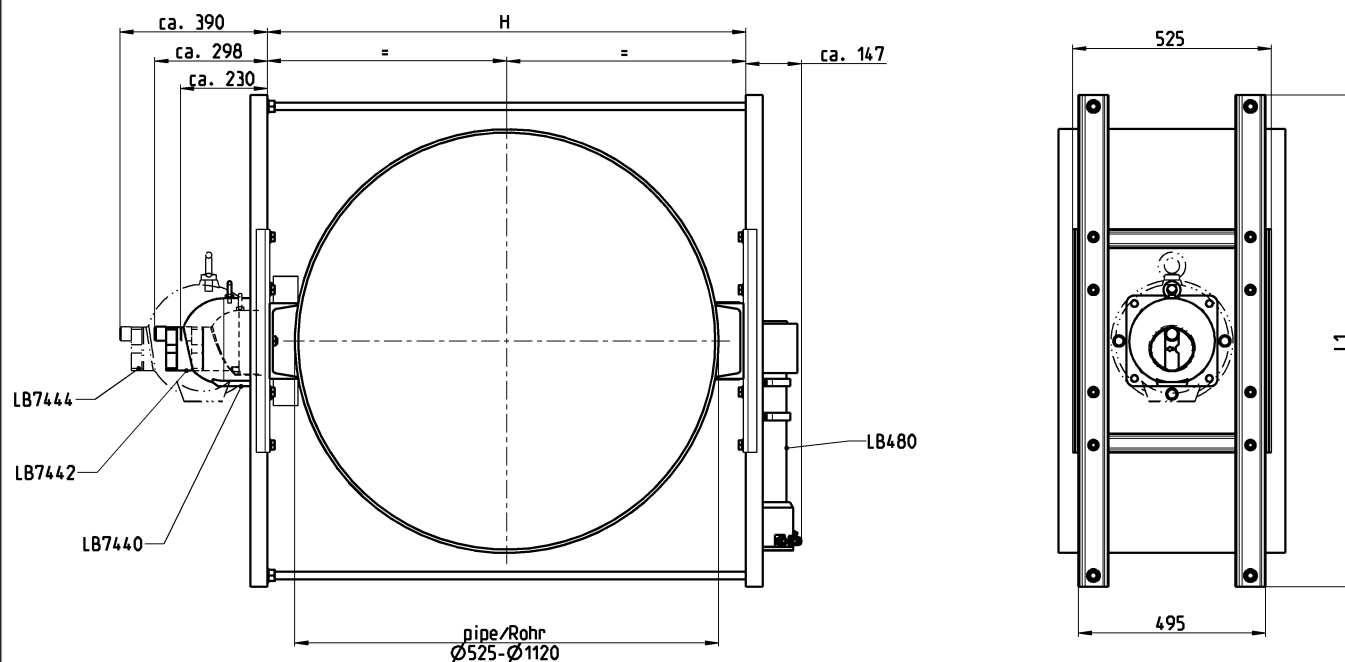
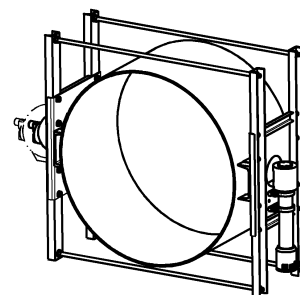
Part No. Id. Nr.	Pipe Diameter Rohrdurchmesser	H	Weight of Clamp. Device Gewicht der Montagevorrichtung
59293	305.0	400	34 kg
	318.0	413	
	323.8	419	
	355.6	451	
	406.4	501	
	457.2	552	
	508.0	603	
	521.0	616	

LB 480

Dimensions in mm
Abmessungen in mm

2.2 Clamping Device 90° for Large Pipe Diameters Montagevorrichtung 90° für große Rohrleitungsquerschnitte

Material	Carbon Steel Stahl St37
Painting Lackierung	Polyurethane, gray Polyurethan, grau



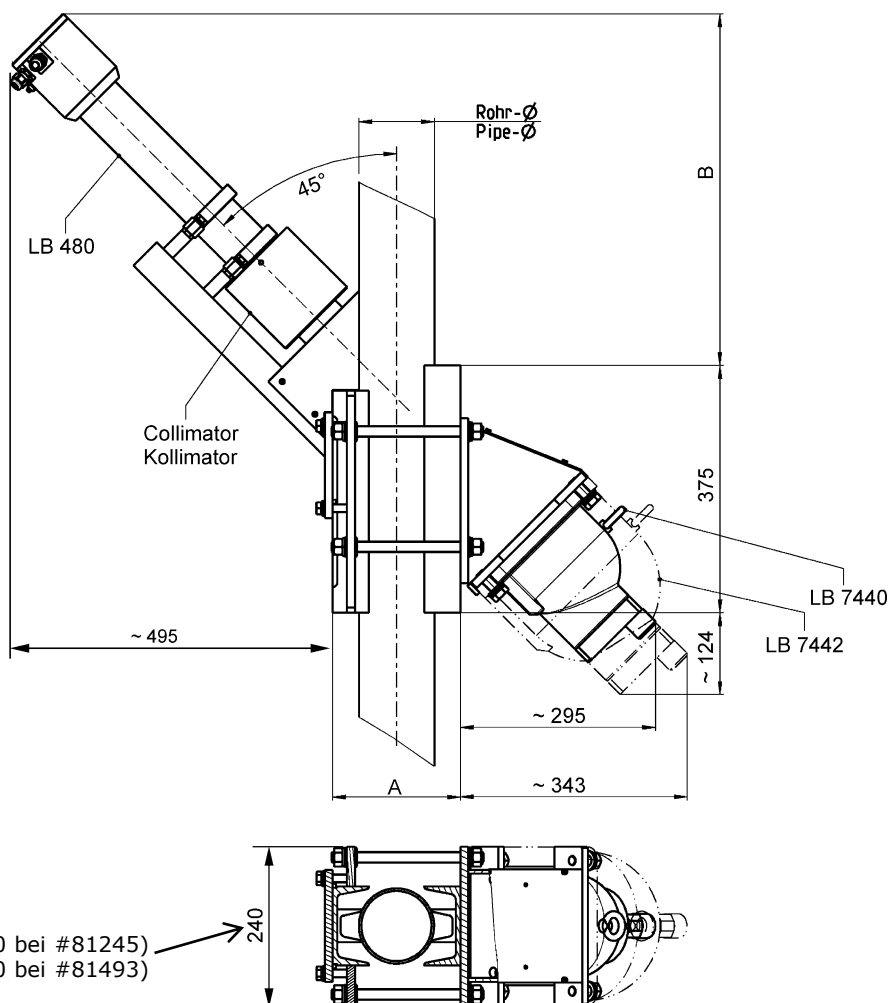
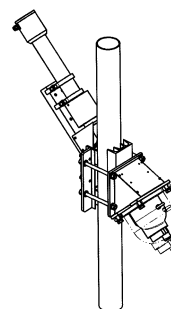
Part No. Id. Nr.	Pipe Ø Range Rohr Ø Bereich	H	Length of the thread bars Länge der Gewindestangen	L1	Weight of Clamp. Device Gewicht der Montagevorrichtung
81491	400 ... 519	643	735	680	73kg
	520 ... 559	686	790	680	74 kg
	560 ... 659	792	890	740	76 kg
	660 ... 759	896	990	840	79 kg
	760 ... 869	1009	1100	950	86 kg
	870 ...1020	1163	1250	1200	95 kg
	1020 ...1120	1265	1350	1300	100 kg

LB 480

Dimensions in mm
Abmessungen in mm

3.0 Clamping Device 45° Montagevorrichtung 45°

Material	Carbon Steel Stahl St37
Painting Lackierung	Polyurethane, gray Polyurethan, grau



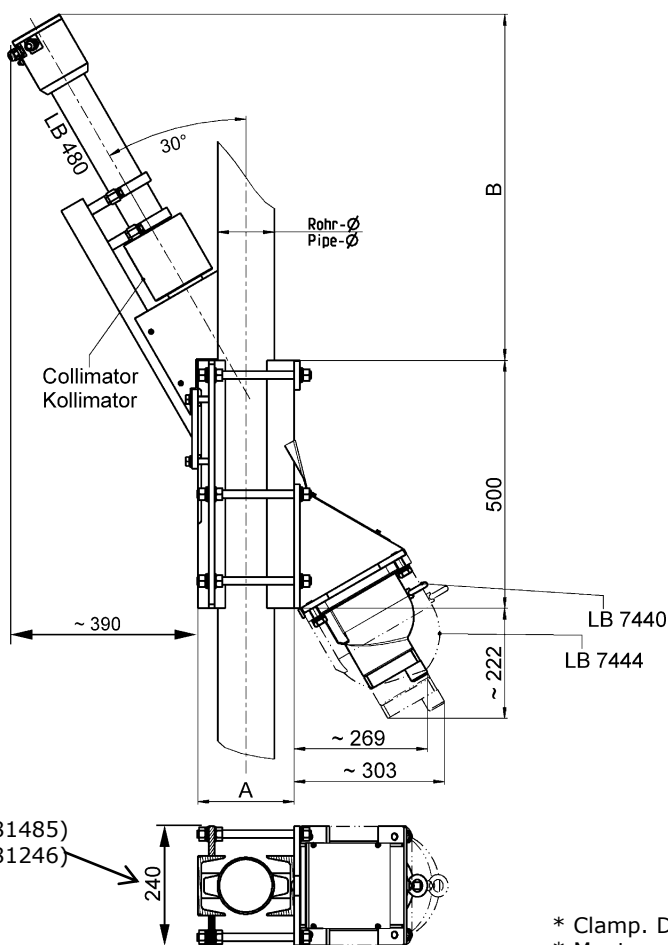
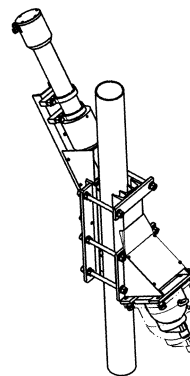
Part No. Id. Nr.	Pipe Ø	A	B	Weight approx Gewicht ca		Part No. Id. Nr.	Pipe Ø	A	B	Weight approx Gewicht ca.
27249	48.3	128	480	31 kg		80793	127.0	165	523	30 kg
27248	60.3	128	480			26997	133.0	177	534	
27250	63.5	133	485			26998	139.7	189	545	
27251	70.0	142	494			26999	146.0	199	555	
27252	76.1	149	500			27000	152.4	209	563	
27253	82.5	157	508			27001	159.0	219	573	
27254	88.9	165	516			27002	165.1	227	581	
27255	95.0	172	522			27003	168.3	232	585	
27256	101.6	179	529			27004	171.0	235	589	
27257	108.0	187	536			27005	177.8	244	598	
27258	114.3	194	543			27006	191.0	261	614	
80794	121.0	201	550			81245	220.0	296	646	
						81493	273.0	356	706	

LB 480

Dimensions in mm
Abmessungen in mm

3.1 Clamping Device 30° Montagevorrichtung 30°

Material	Carbon Steel Stahl St37
Painting Lackierung	Polyurethane, gray Polyurethan, grau



(332 → #81485)
(280 → #81246)

* Clamp. Device Weight (kg)
* Montagevorrichtung Gewicht (kg)

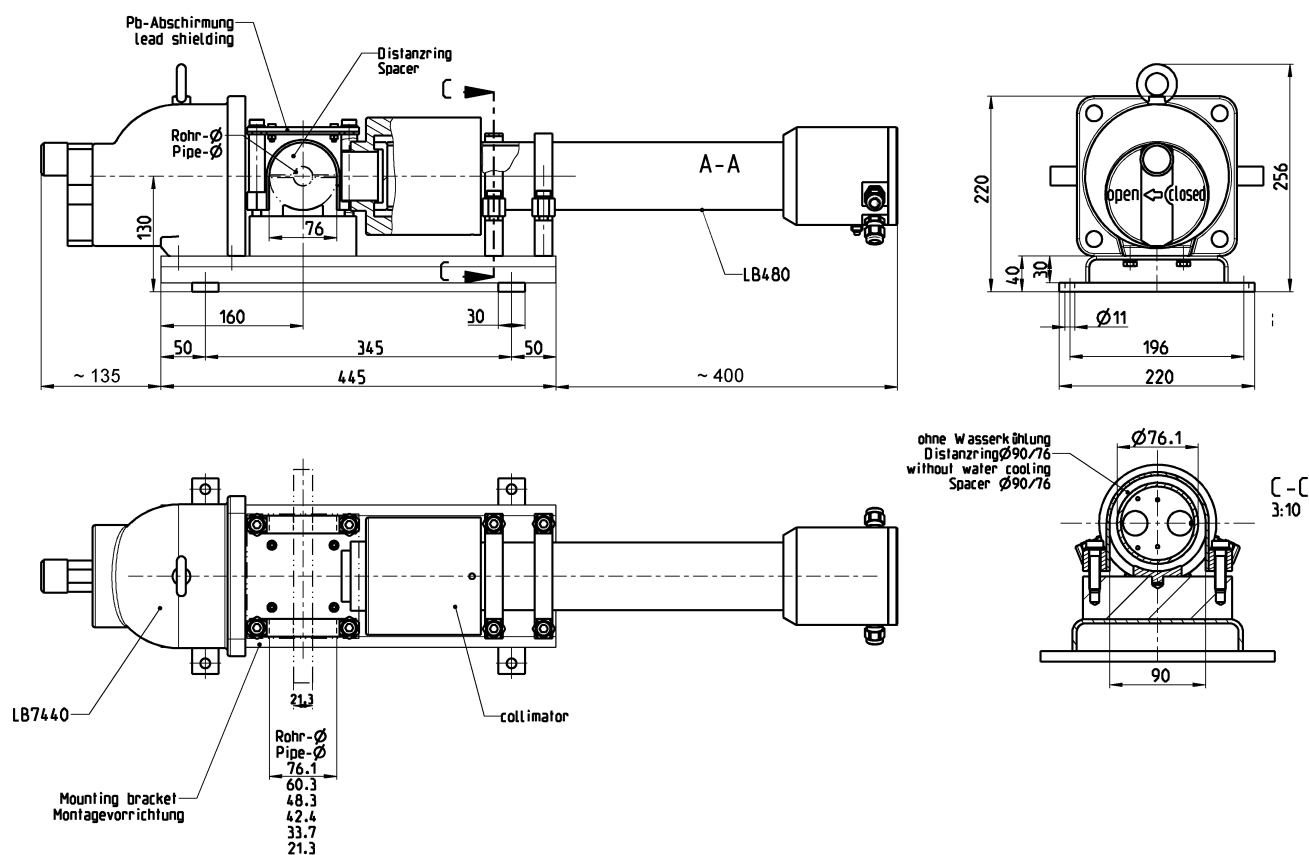
Part No. Id. Nr.	Pipe Ø	A	B	Weight* Gewicht*	Part No. Id. Nr.	Pipe Ø	A	B	Weight* Gewicht*
25964	48.3	128	602	39 kg	80791	127.0	165	674	39 kg
80792	60.3	128	602		26655	133.0	177	692	
25971	63.5	133	610		26656	139.7	189	712	
25972	70.0	142	625		26657	146.0	199	728	
25973	76.1	149	636		26658	152.4	209	741	
25974	82.5	157	650		26659	159.0	219	759	
25975	88.9	165	663		26660	165.1	227	773	
25976	95.0	172	674		26661	168.3	232	780	
25977	101.6	179	686		26662	171.0	235	787	
25978	108.0	187	698		26663	177.8	244	803	
25979	114.3	194	710		26664	191.0	261	830	
25980	121.0	201	723		81246	220.0	296	884	40 kg
					81485	273.0	356	988	

LB 480

Dimensions in mm
Abmessungen in mm

4.0 Clamping Device 90° for Small Pipe Diameters Montagevorrichtung 90° für kleine Rohrleitungsquerschnitte

Material	Carbon Steel Stahl St37
Painting Lackierung	Polyurethane, gray Polyurethan, grau



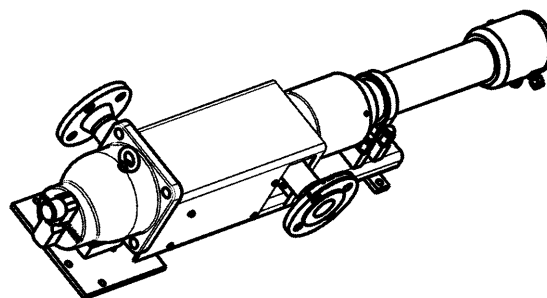
Part No. Id. Nr.	Pipe Ø	Weight of Clamp. Device Gewicht der Montagevorrichtung
47292-01	21.3	70 kg
47292-02	33.7	
47292-03	42.4	
47292-04	48.3	
47292-05	60.3	
47292-06	76.1	

LB 480

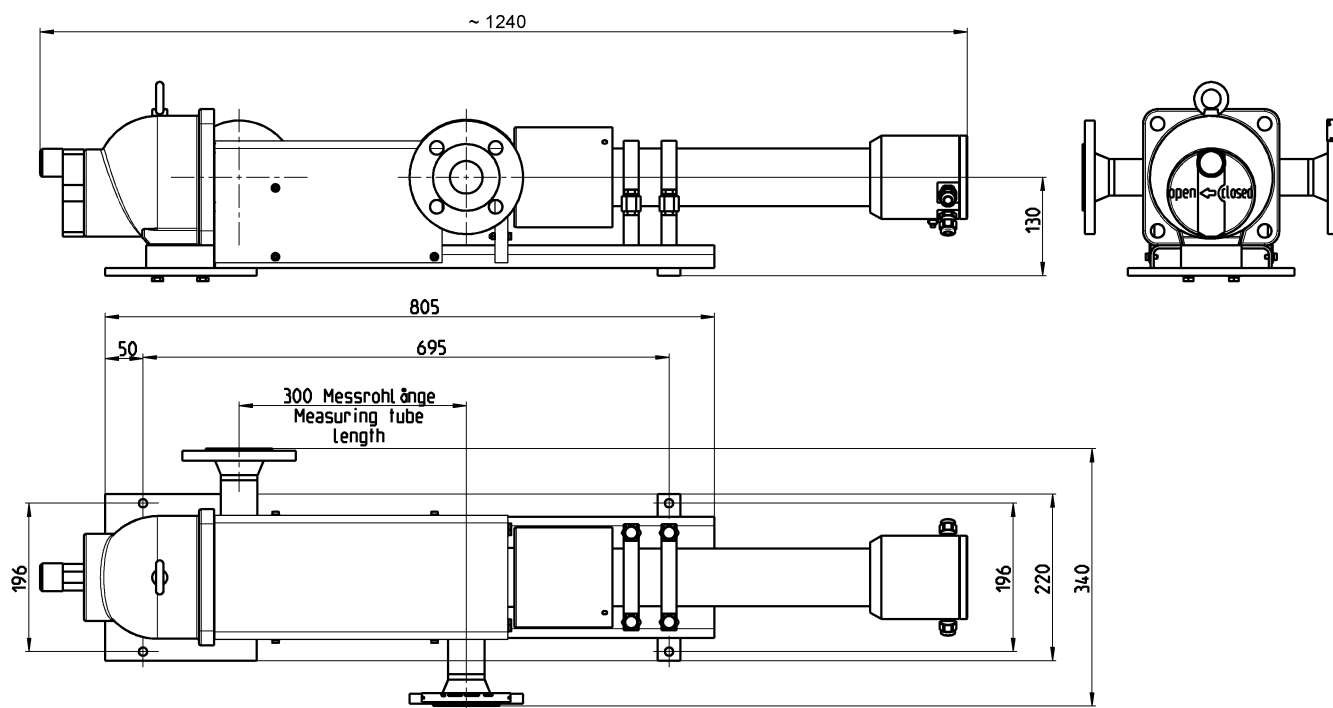
Dimensions in mm
Abmessungen in mm

4.1 S + U Pipe Clamping Device S + U förmige Montagevorrichtung

Material	Carbon Steel Stahl St37
Painting Lackierung	Polyurethane, gray Polyurethan, grau



Part No. / Id. Nr. 21087



Verschiedene Versionen erhältlich.

Obige beispielhafte Abbildung zeigt die Messstrecke in:

- S-Form
- DN40
- Messrohrlänge 300
- PN16
- ohne Temp.-Fühler

Various versions available.

The drawing above shows exemplarily a clamping device with:

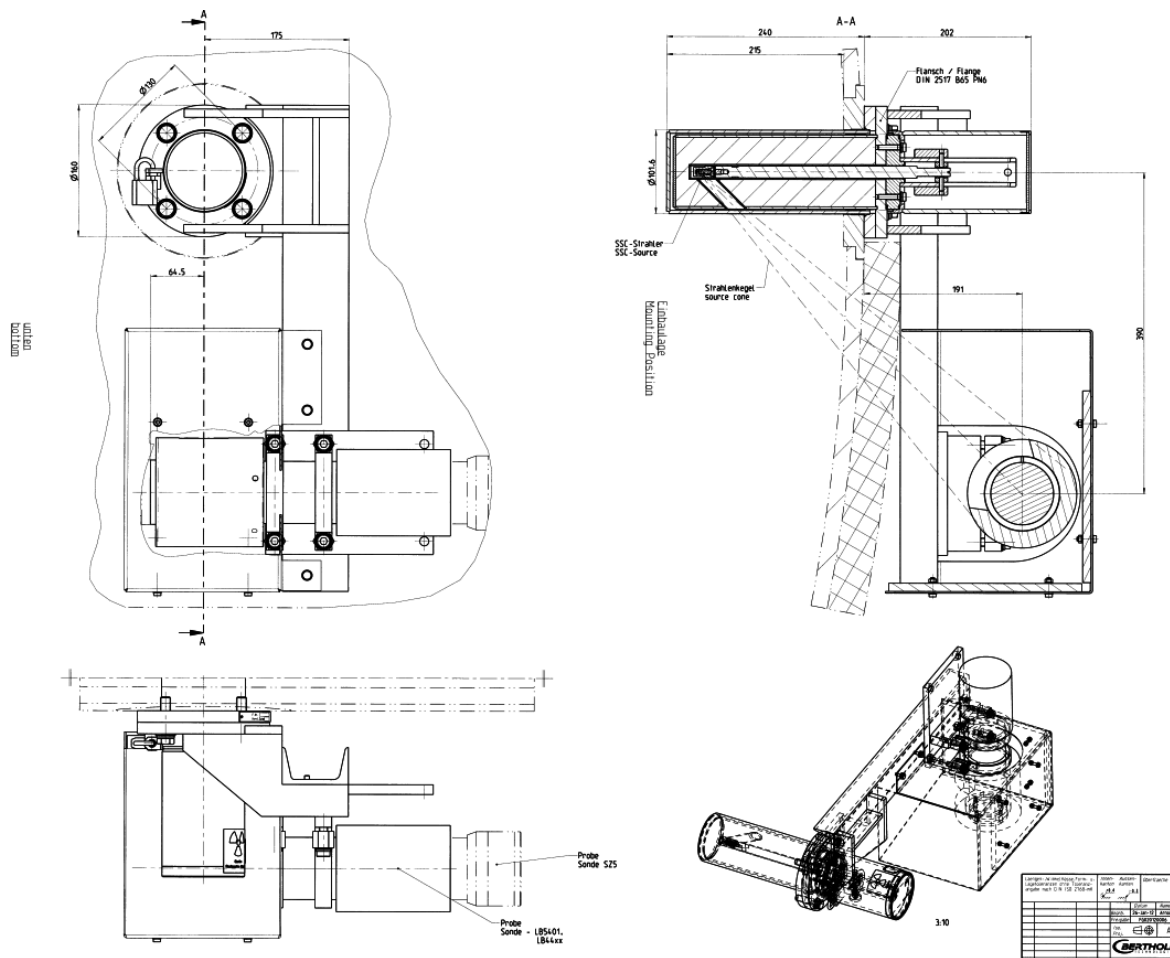
- S-shaped
- DN40
- pipe length 300
- PN16
- without temp.-sensor

LB 480

Dimensions in mm
Abmessungen in mm

5.0 Immersion Shielding Tauchabschirmung

Material Dip Pipe Material Tauchrohr	Stainless Steel AISI 316Ti 1.4571
Painting Lackierung	Polyurethane, gray Polyurethan, grau



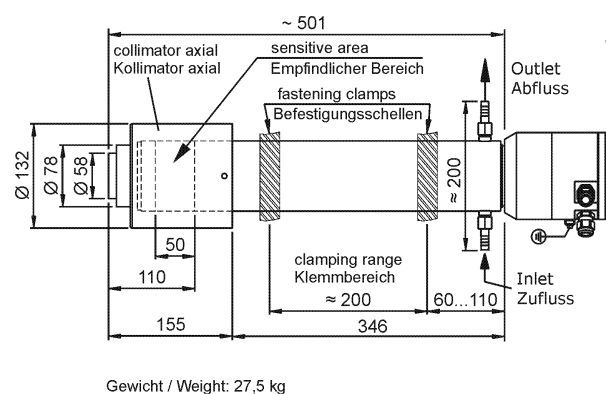
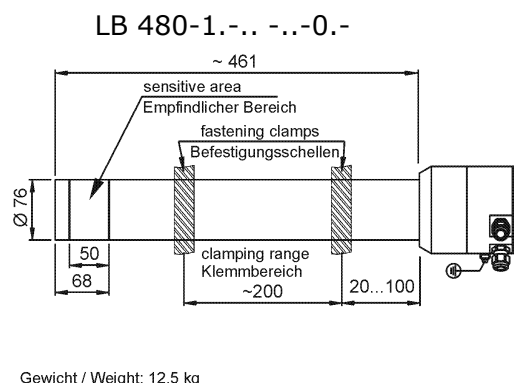
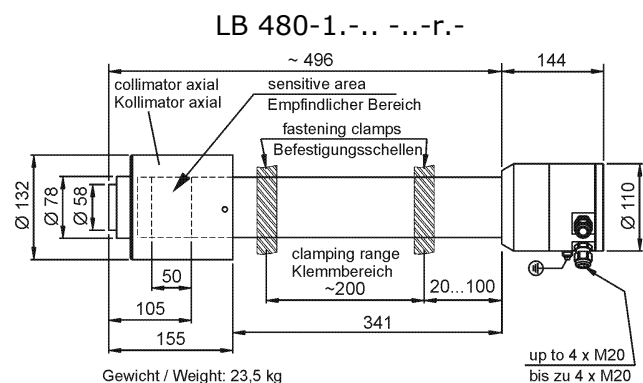
Part No. Id. Nr.	Measuring Range Messbereich
58959-01	depending on der vessel diameter

LB 480

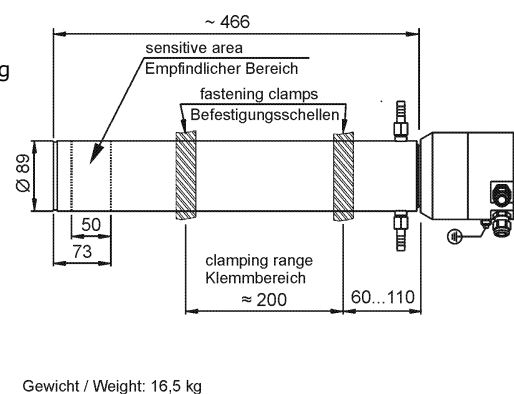
Dimensions in mm
Abmessungen in mm

6.0 CrystaSENS (Version f. Zone 1/2)

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

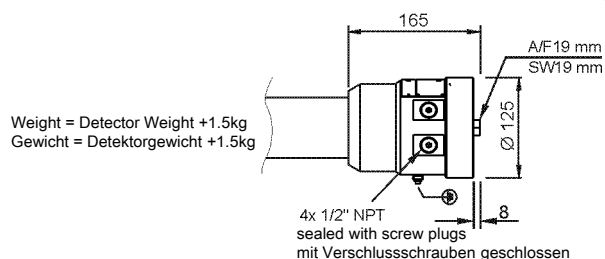


with water cooling
mit Wasserkühlung



6.1 CrystaSENS (Version f. Divisions 1/2)

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



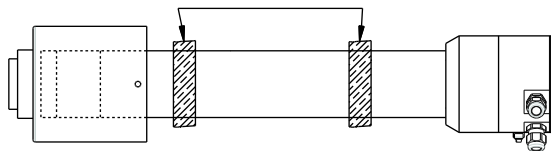
The detector version for divisions (NEC/CEC) differs only in the terminal housing, compared to the standard version illustrated above. The dimensions of this terminal housing are illustrated in this drawing.

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.

6.2 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		✓

6.3 Mounting Clamps for Detector Befestigungsschellen für Detektor

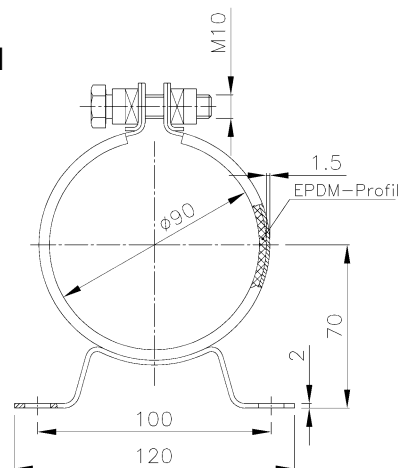
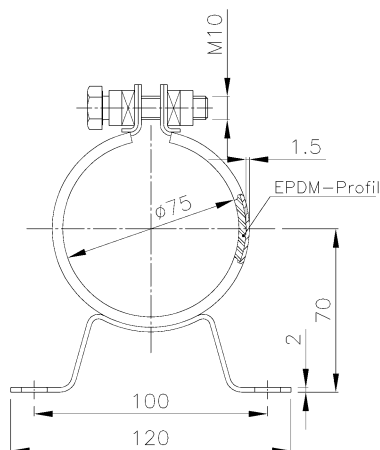


for Detectors without water cooling
für Detektoren ohne Wasserkühlung

for Detectors with water cooling
für Detektoren mit Wasserkühlung



Material
316Ti
1.4571



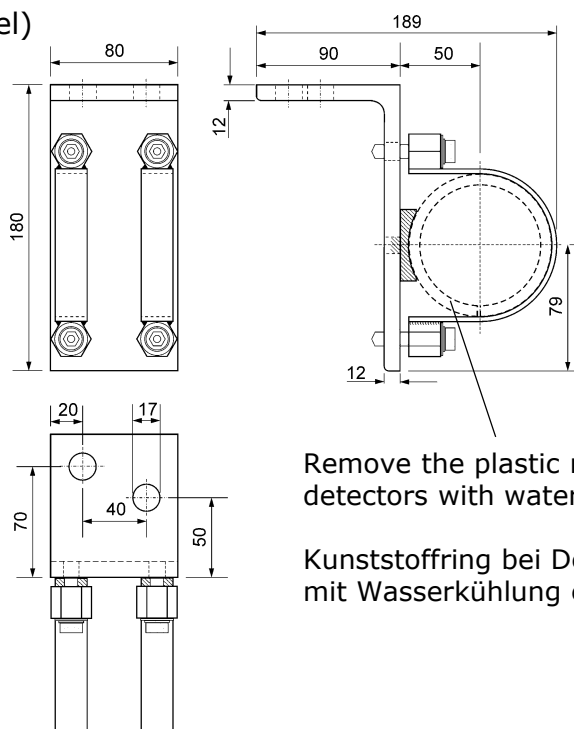
Id. Nr. 31346 (1 set = 2 clamps)
Id. Nr. 31345 (single clamp)

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)

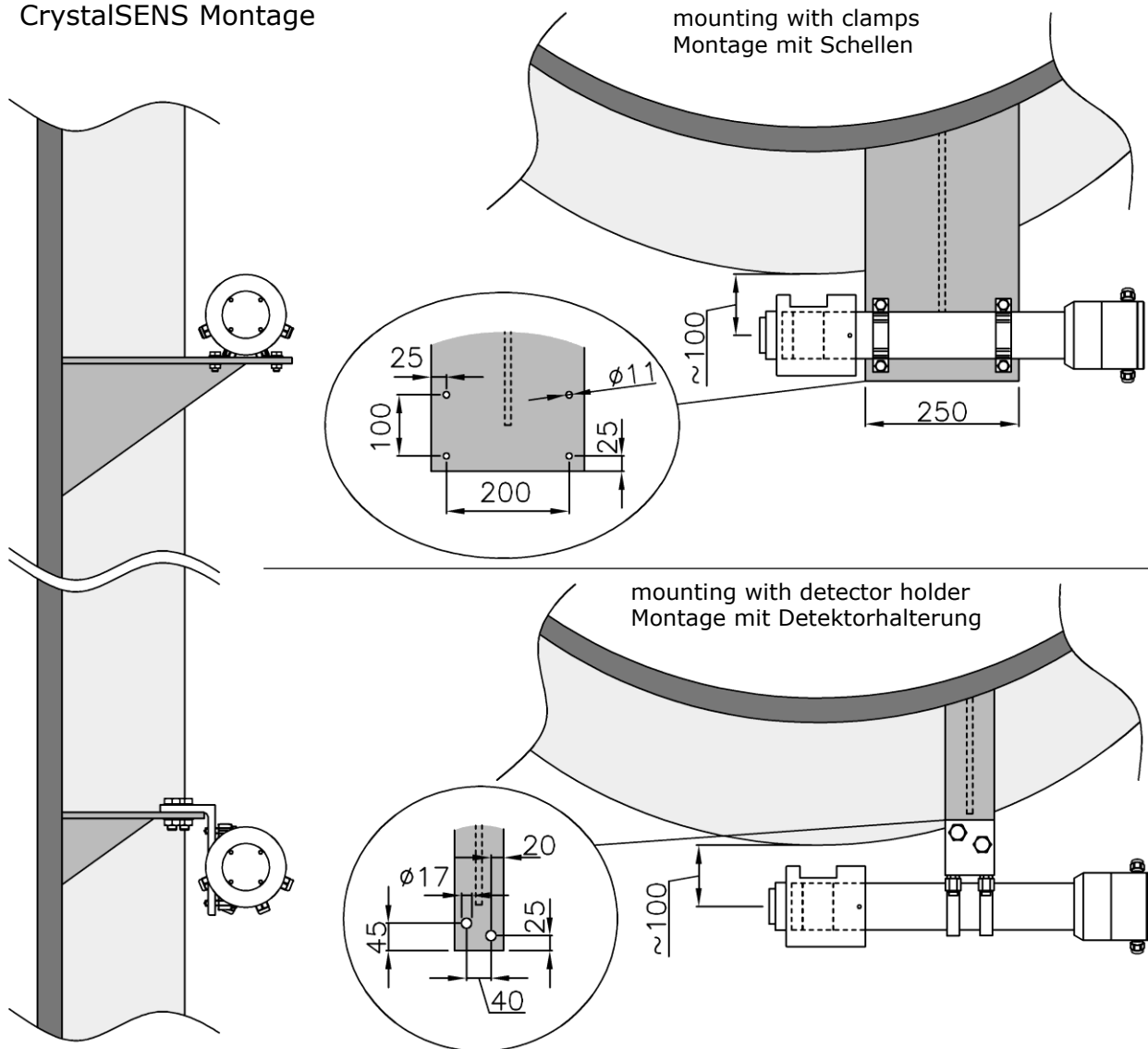
Part No. Id. Nr.	for Detector für Detektor
39246	without water cooling ohne Wasserkühlung
39247	with water cooling mit Wasserkühlung



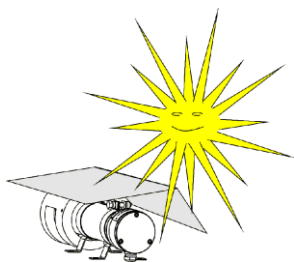
Remove the plastic ring for
detectors with water cooling.

Kunststoffring bei Detektoren
mit Wasserkühlung entfernen.

6.4 CrystalSENS Mounting CrystalSENS Montage



6.5 Sun Roof against Strong Sun Radiation Sonnendach gegen starke Sonneneinstrahlung

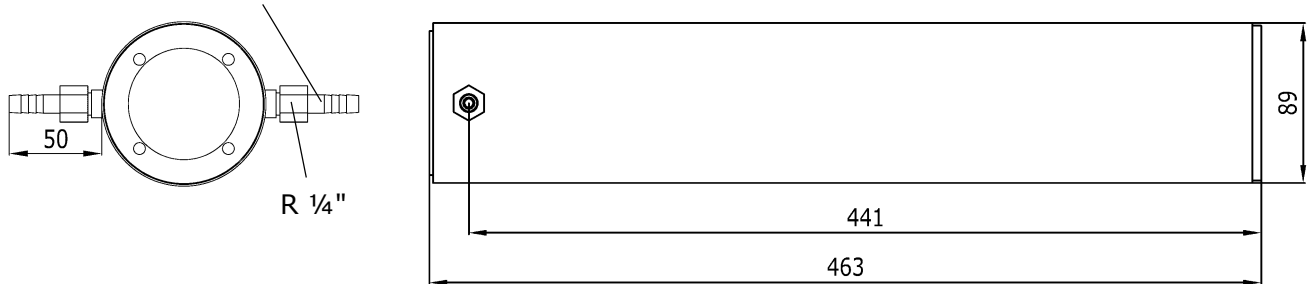


Direct sun radiation can overheat the detector. If the detector temperature can reach more than 50°C, a suitable sun roof must be installed. The heating of the detector by thermal radiation from the vessel can also be moderated by a thermal sheet, e.g. by a thin metal plate. For each detector a water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

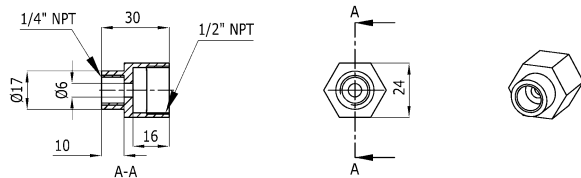
6.6 CrystaSENS Water Cooling Jacket and Adaptor Fittings CrystaSENS Wasserkühlung und Adapter Anschlussstücke

pipe connection diameter 10
Schlauchanschluss Ø 10



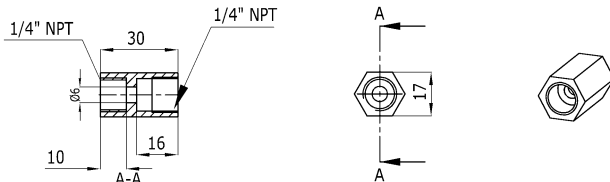
Fitting adaptor for standard water cooling Rp 1/4" → 1/2" NPT
stainless steel 304, part no: 47189

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



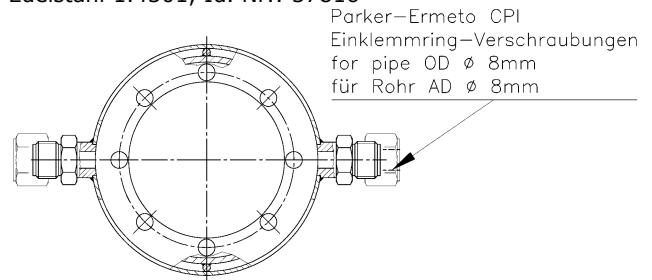
Fitting adaptor for standard water cooling Rp 1/4" > 1/4" NPT
stainless steel 304, part no: 46743

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



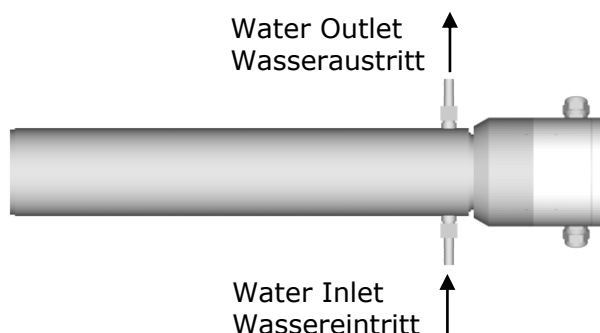
Further fitting adaptors for standard water cooling jacket:
Rp 1/4" > 1/2" NPT male, stainless steel 304, part no: 06352
Rp 1/4" > 1/4" NPT male, stainless steel 304, part no: 06349
Weitere Adapter für die Standard-Wasserkühlung:
Rp 1/4" > 1/2" NPT Außengewinde, 1.4301, Id. Nr.: 06352
Rp 1/4" > 1/4"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

6.7 Water Cooling Installation Instruction Anweisung zur Installation der Wasserkühlung

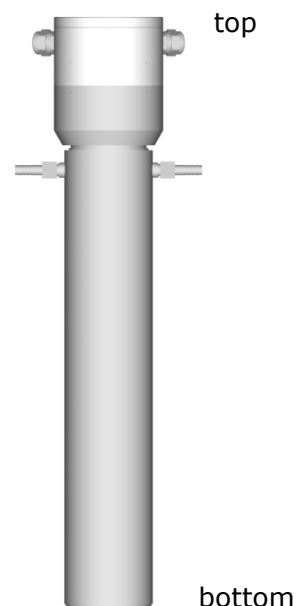
Horizontal Detector Installation Horizontale Detektor Installation



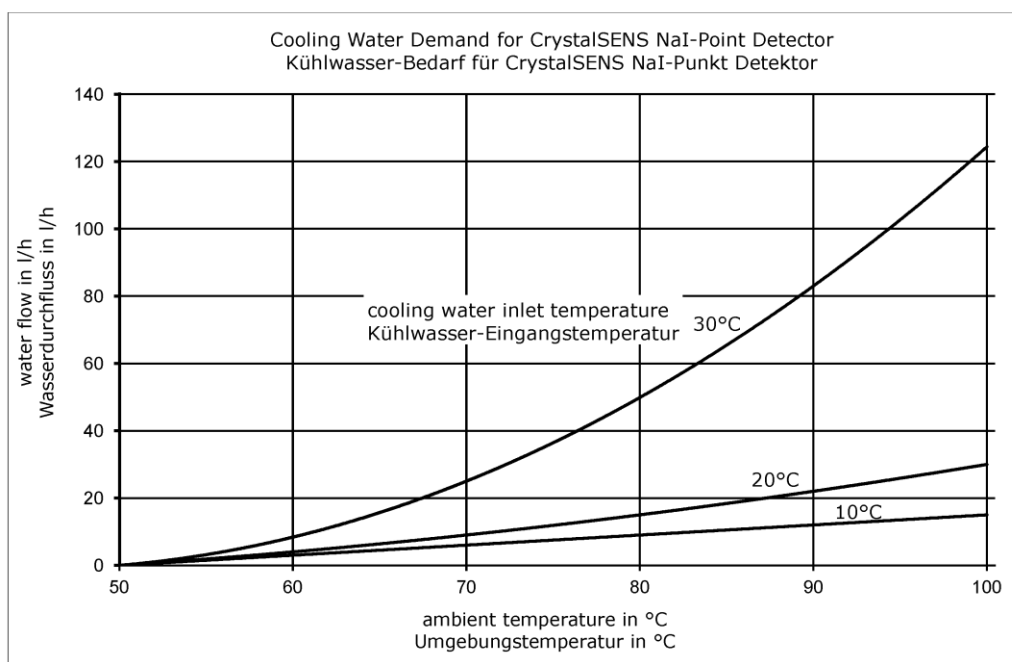
In order to fill the entire water cooling jacket, incoming water must enter from the bottom. Damit sich die Wasserkühlung vollständig mit Wasser füllt, muss der Wasserzufluss von unten erfolgen.

Vertical Detector Installation Vertikale Detektor Installation

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

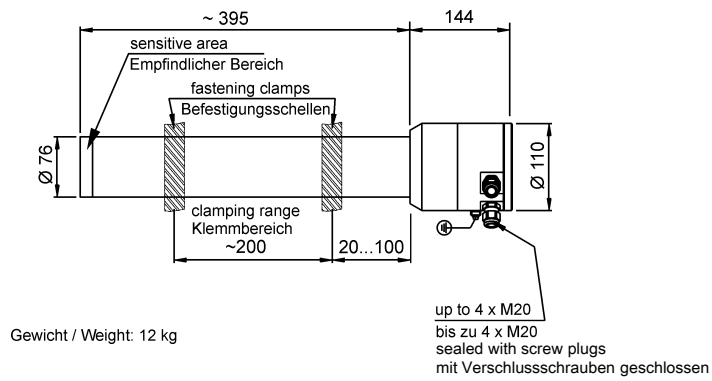


6.8 Detector Cooling Water Demand Detektor Kühlwasserbedarf



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

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

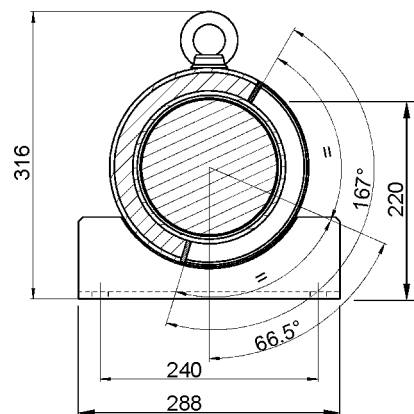
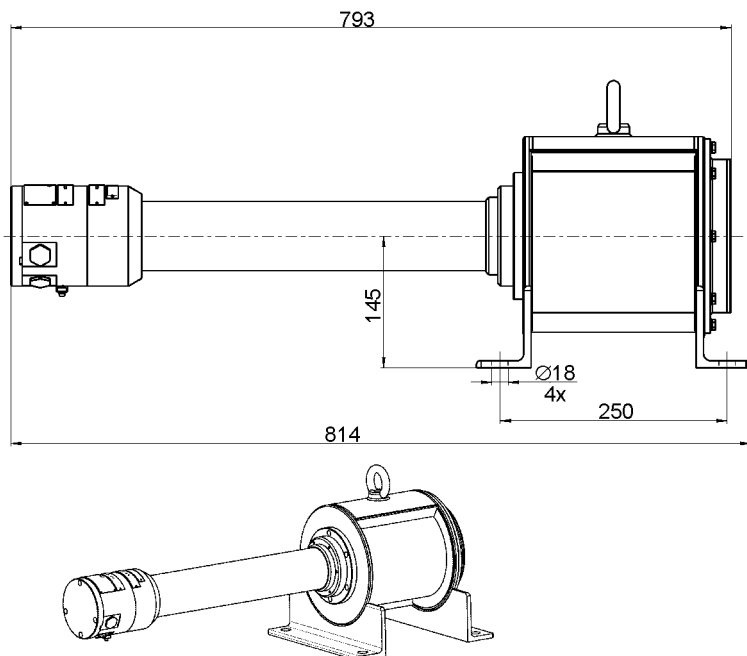


LB 480

Dimensions in mm
Abmessungen in mm

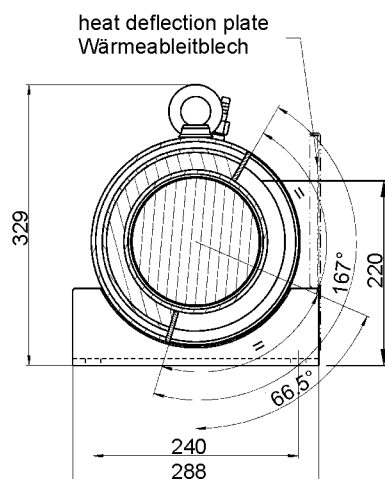
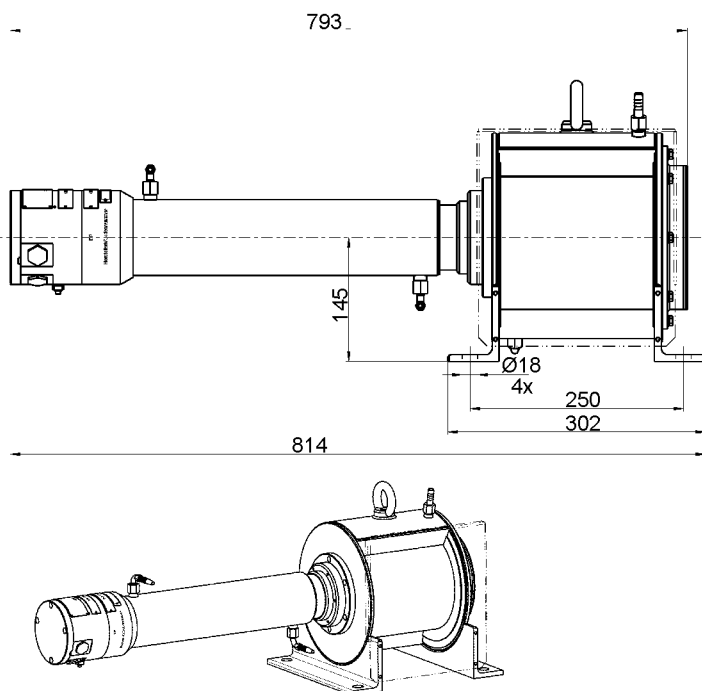
7. SuperSENS

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



Weight/Gewicht 52 kg

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

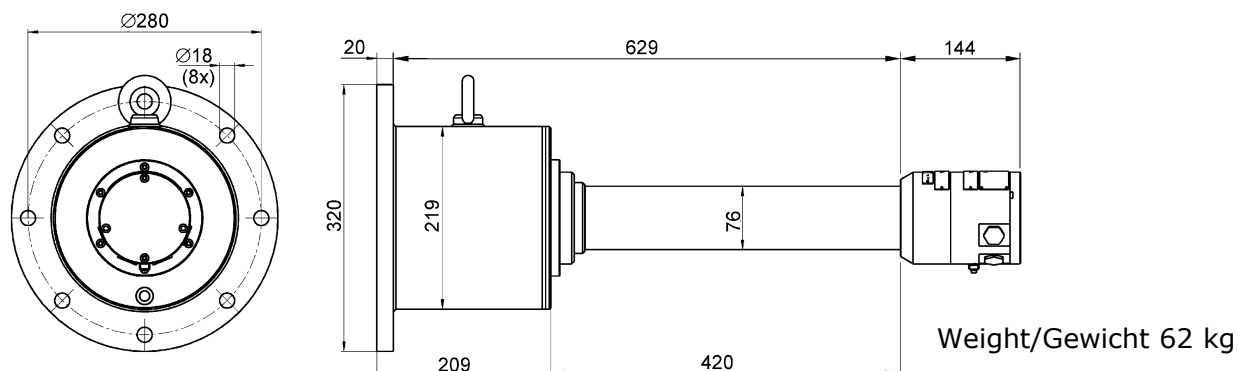


Weight/Gewicht 59 kg

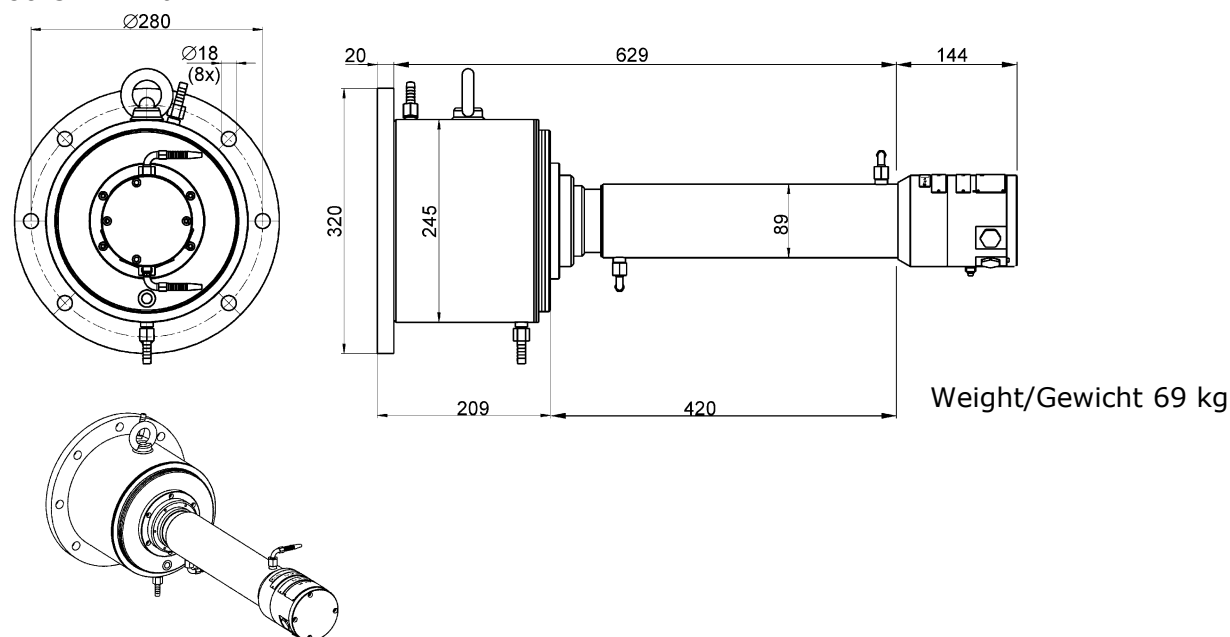
LB 480

Dimensions in mm
Abmessungen in mm

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

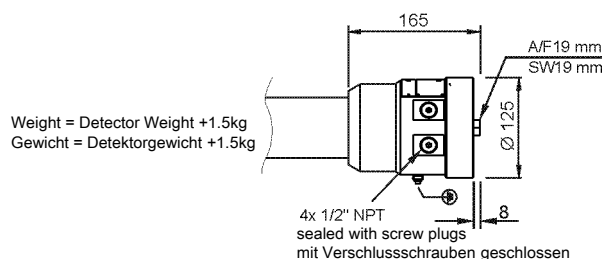


7.4 with Axial Irradiation and Water Cooling mit frontaler Einstrahlung und Wasserkühlung LB 480-32-...-a.



7.5 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.

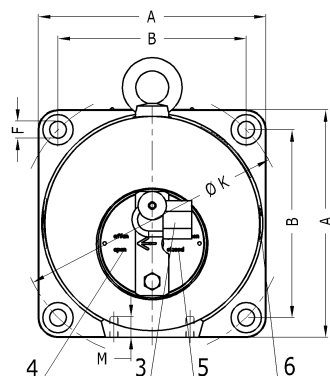
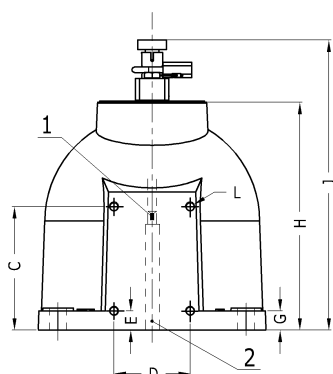
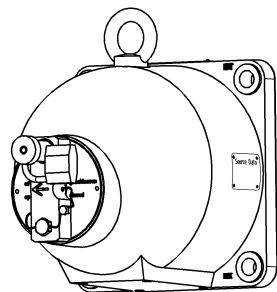
Die Detektor-Version mit Ex-Zulassung für Divisions (NEC/CEC) unterscheidet sich ausschließlich im Anschlusskopf, verglichen mit der oben aufgeführten Standardvariante. Die Abmessungen für den Anschlusskopf sind aus dieser Zeichnung zu entnehmen.

LB 480

Dimensions in mm
Abmessungen in mm

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

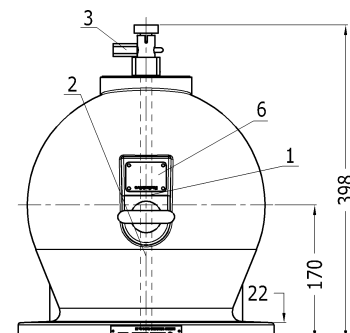
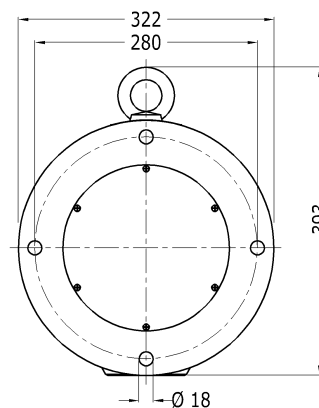
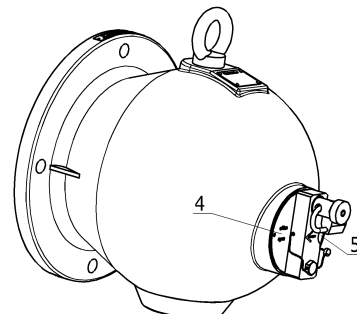
LB 7440 / 7442 / 7445 / 7446



- 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

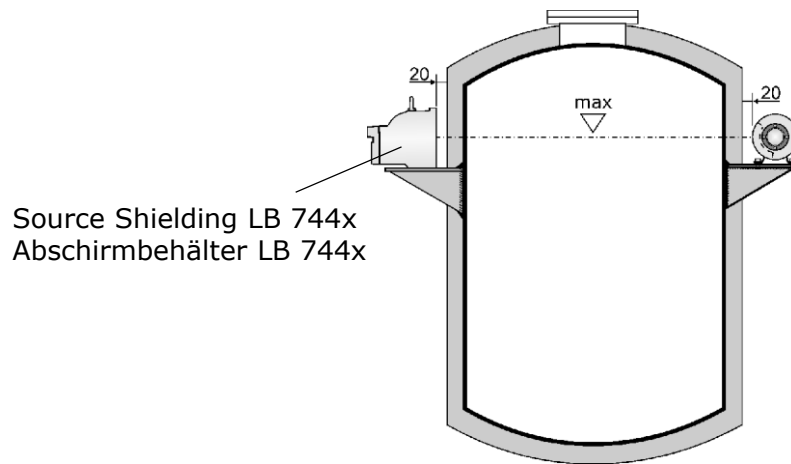
LB 7444



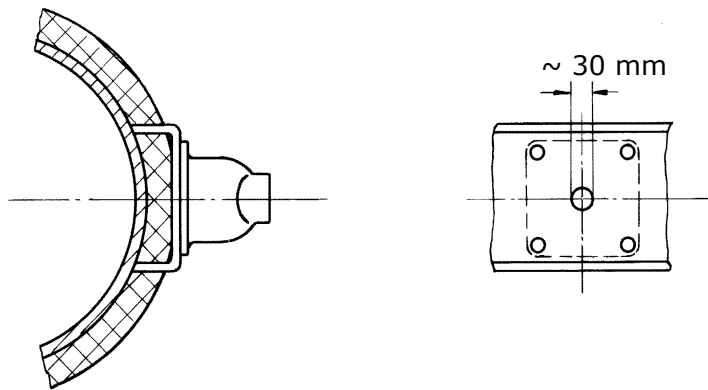
Housing Gehäuse	Type Typ	A	B	C	D	E	FØ	G	H	J	KØ	L	M	Flange Flansch		kg
Carbon Steel St 37	LB 7440-D-CR LB 7445-D-CR	180	142	75	60	15	18	20	173	238	200	M8	12	ND 125 PN 6	11°	31
Super Duplex UNS 32750 SAF 2507 1.4410	LB 7440-DE-CR LB 7445-DE-CR															
	LB 7442-DE-CR	240	198	130	80	20	18	20	242	306	280	M10	14	ND 200 PN 6	7°	81
Carbon Steel St 37	LB 7442-D-CR LB 7446-D-CR															
	LB 7444-CR	Dimensions in drawing / Abmessungen in Zeichnung												ND 200 PN 6	6°	170

 Radiation Angle of the Shielding / Abstrahlwinkel der Abschirmung

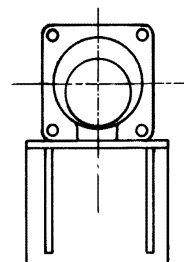
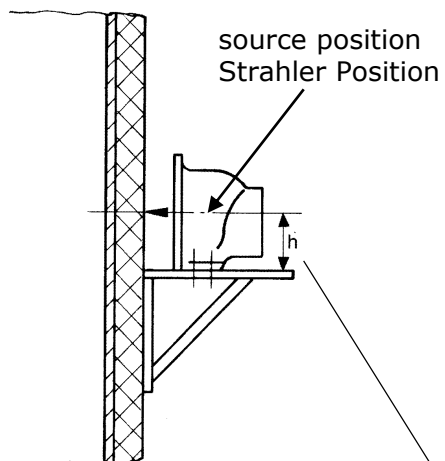
8.1 Mounting Proposal for Source Shielding LB 744x Montagevorschlag für Abschirmbehälter LB 744x



Source Shielding LB 744x
Abschirmbehälter LB 744x



Flange Installation
Flanschinstallation

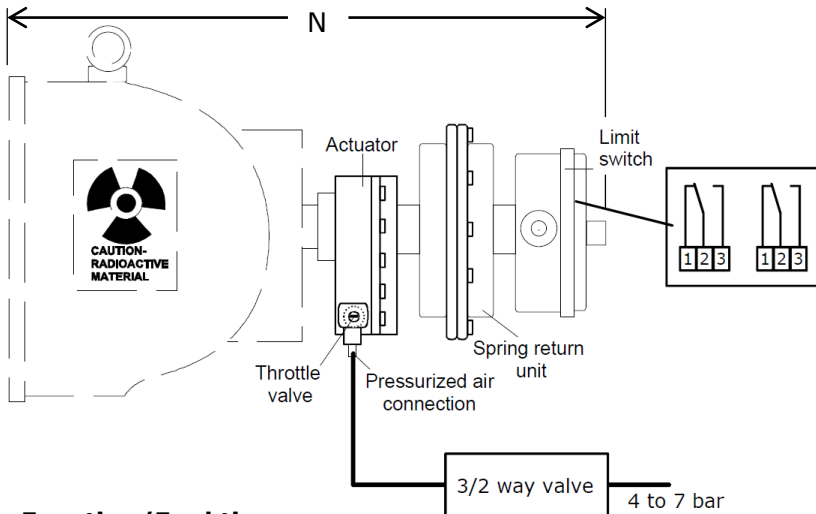


Pedestal Installation
Sockelinstallation

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

8.2 Pneumatic Actuator for Source Shielding LB 744x

Pneumatischer Verschlussantrieb für Abschirmbehälter LB 744x



N approx N 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

Function/Funktion:

The pressurized air moves the shutter to the OPEN position. If the pressurized air is turned off or in case of failure the moveable shutter is turned back to the CLOSED position by the spring return unit.
Die Druckluft bewegt den Verschlusskern in die Position AUF. Wenn der Druck abfällt, bewegt sich der Verschlusskern, mit Hilfe der Rückzugfeder, zurück in die Position ZU (Fail Safe).

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 OPEN/CLOSED

Endschaltereinheit, Optionen für Signalisierung AUF / ZU

Option I:	IP 65, 2 contacts (OPEN/CLOSED) IP 65, 2 Kontakte (AUF/ZU)
Option II:	2 contacts (OPEN/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 Näherungsinitiatoren für Eigensichere Speisung

The pneumatic actuator is equipped with a throttle valve. The factory setting of the valve is set such that the shield opening and closing process takes at least 2 s; otherwise the shield may get damaged!

Der pneumatische Antrieb hat ein Drosselventil. Diese Ventil ist werkseitig so eingestellt, dass die Bewegung für AUF/ZU mindestens 2 s dauert. Im anderen Fall könnte die Abschirmung beschädigt werden.

8.3 Components for Pneumatic Actuator Einzelteile für Pneumatischen Verschlussantrieb

Pneumatic Actuator
Pneumatischer Antrieb

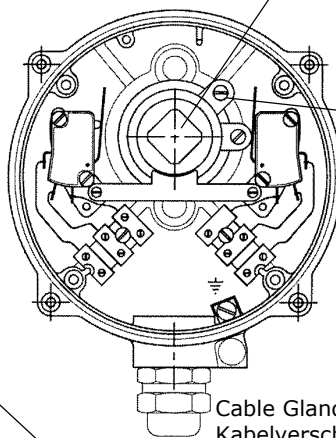
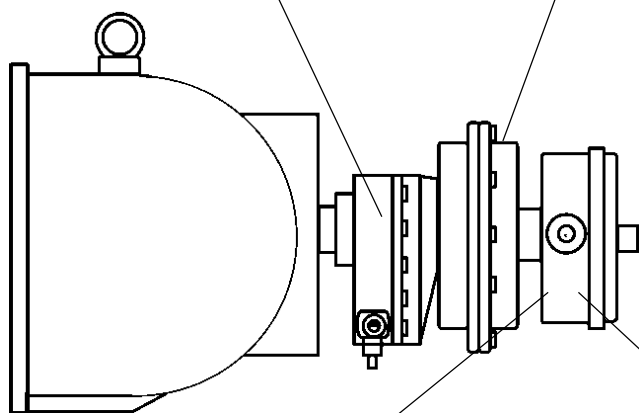


Return Spring
Rückstellfeder



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

Außenvierkant zur Stel-
lungsanzeige, oder manueller
Betätigung



adjustable cam
verstellbare Nocken

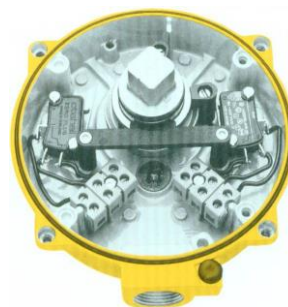
Cable Gland
Kabelverschraubung

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
	30	7	5
	75	1	1
	125	0,5	0,06
	250	0,25	0,03

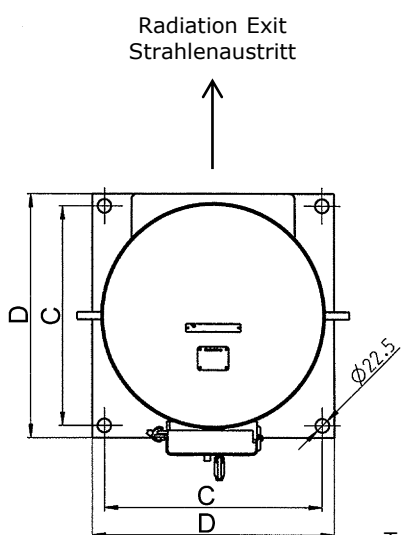
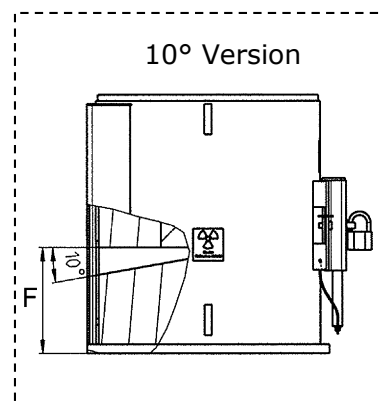
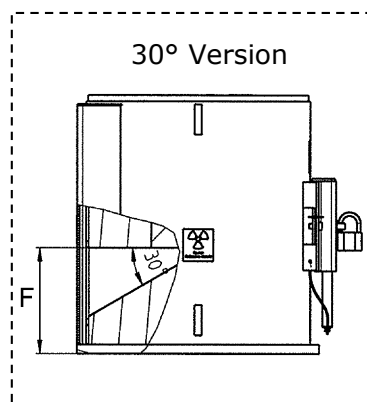
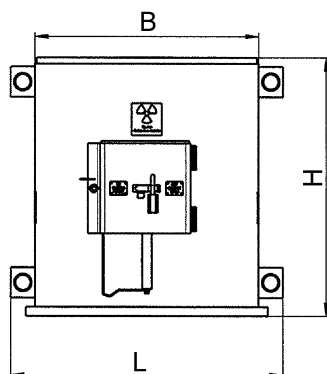
Contact Rating
Kontaktbelastbarkeit

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

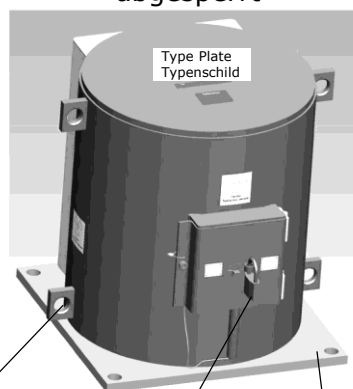
LB 480

Dimensions in mm
Abmessungen in mm

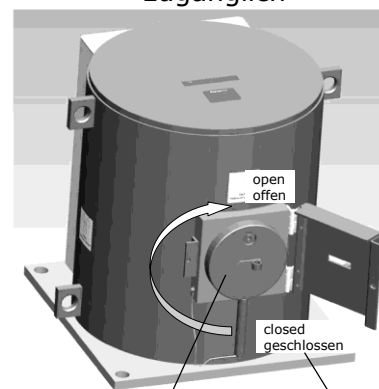
8.4 Point Source Shielding LB 8030/8040 Punktstrahler-Abschirmbehälter LB 8030/8040



Shutter mechanism
locked
Strahlermechanismus
abgesperrt



Shutter mechanism
accessible
Strahlermechanismus
zugänglich



Transport Lug
Transportöse

Mounting Flange
Montageflansch

Pad Lock
Vorhängeschloss

Source Locking
Mechanism
Strahlerverschluss-
mechanismus

Source Position
Strahler Position

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

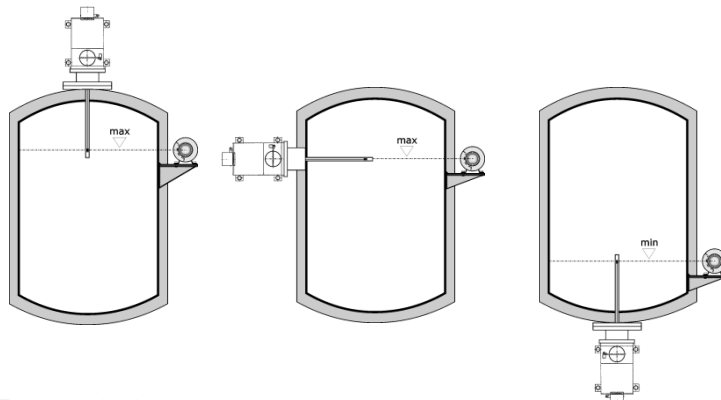
Type Typ	B	C	D	F	H	L		kg
LB 8030-01	370	360	400	174	425	450	30°	516
LB 8030-02							10°	
LB 8040-01	456	460	500	305	567	536	30°	1014
LB 8040-02							10°	

Radiation Angle of the Shielding / Abstrahlwinkel der Abschirmung

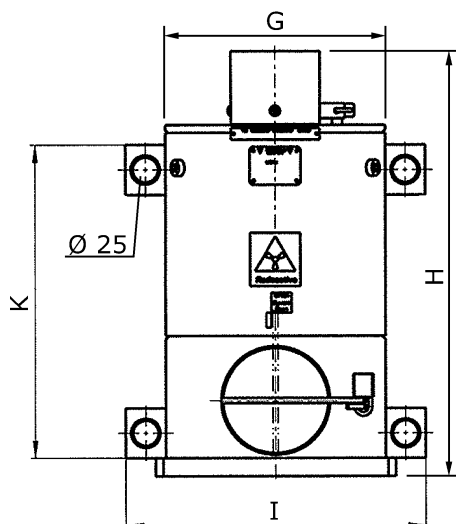
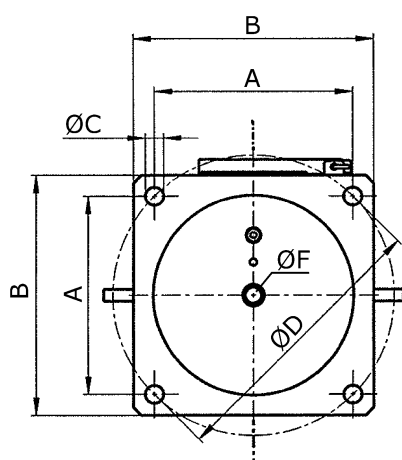
LB 480

Dimensions in mm
Abmessungen in mm

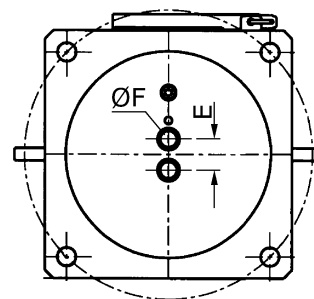
8.5 Point Source Flange Shielding LB 81xx for Dip Pipe Installations Punktstrahler-Flansch-Abschirmbehälter LB 81xx für Tauchrohr-Applikationen



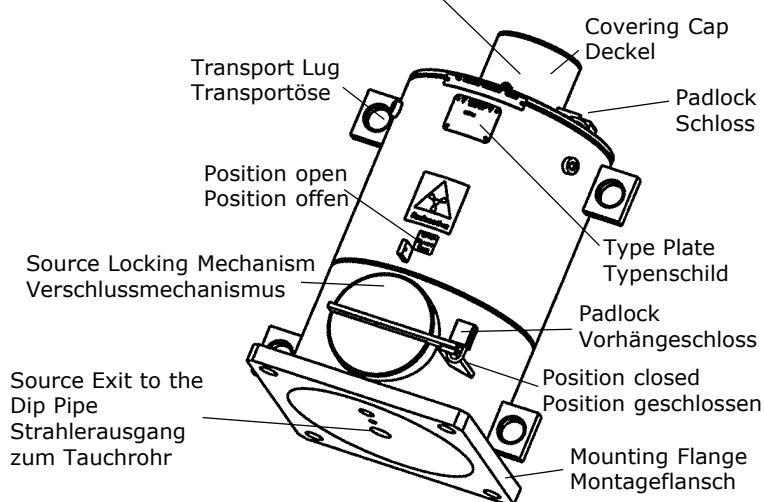
Type/Typ LB 8115 + LB 8120



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

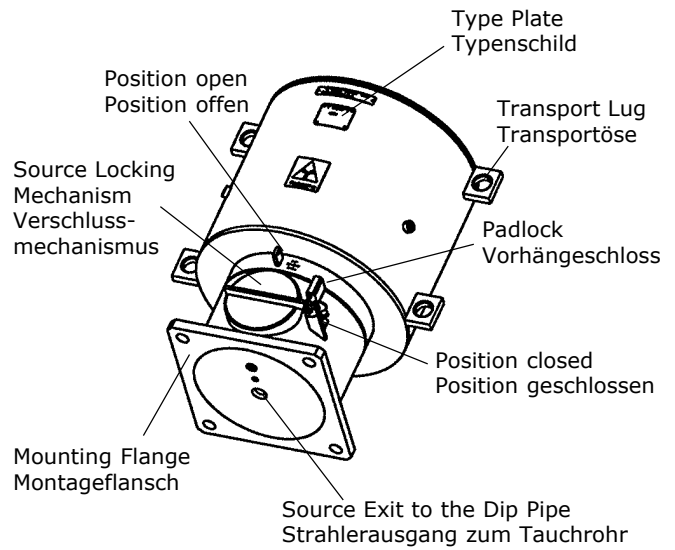
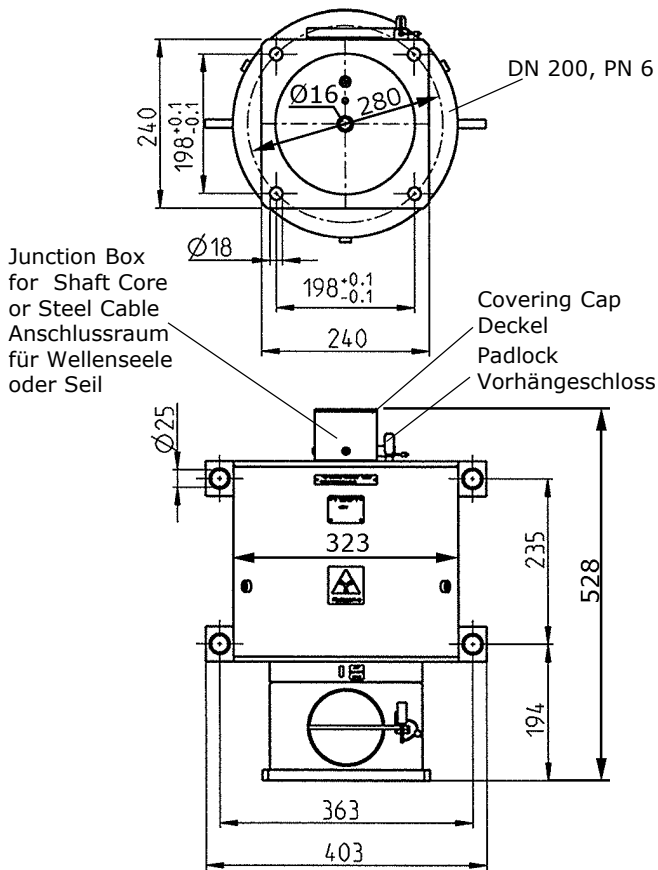


Type Typ	A	B	ØC	ØD	E	ØF	G	H	I	K	kg
LB 8115-01 LB 8115-20	141,5	180	18	200	- 30	16	159	350	239	237	60 kg
LB 8120-01 LB 8120-20	198	240	18	280	- 30	16	219	425 429	300	313	137 kg

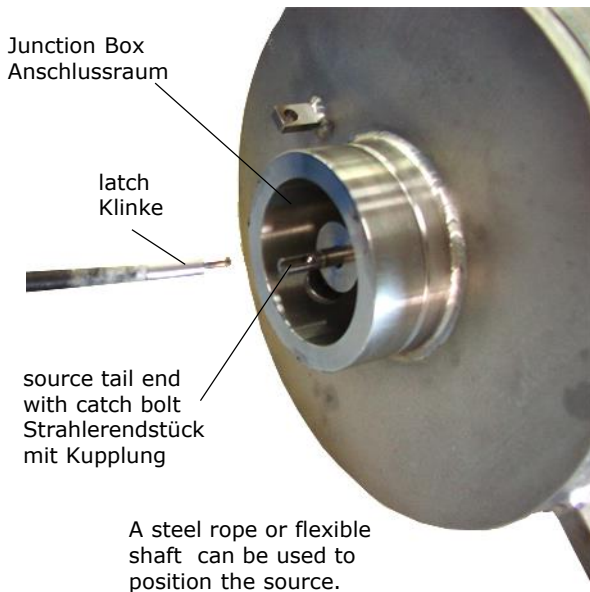
LB 480

Dimensions in mm
Abmessungen in mm

Type/Typ LB 8125-01

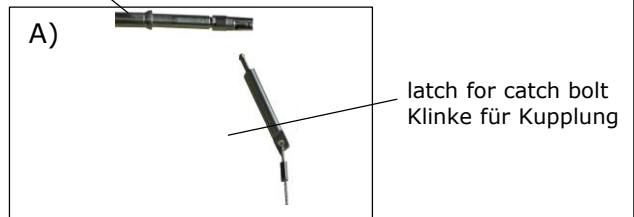
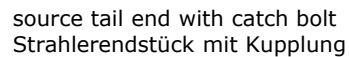


8.6 Operation of Flange Shielding Bedienung der Flanschabschirmung



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

Zur Positionierung kann ein Stahlseil oder eine Wellenseele genutzt werden.



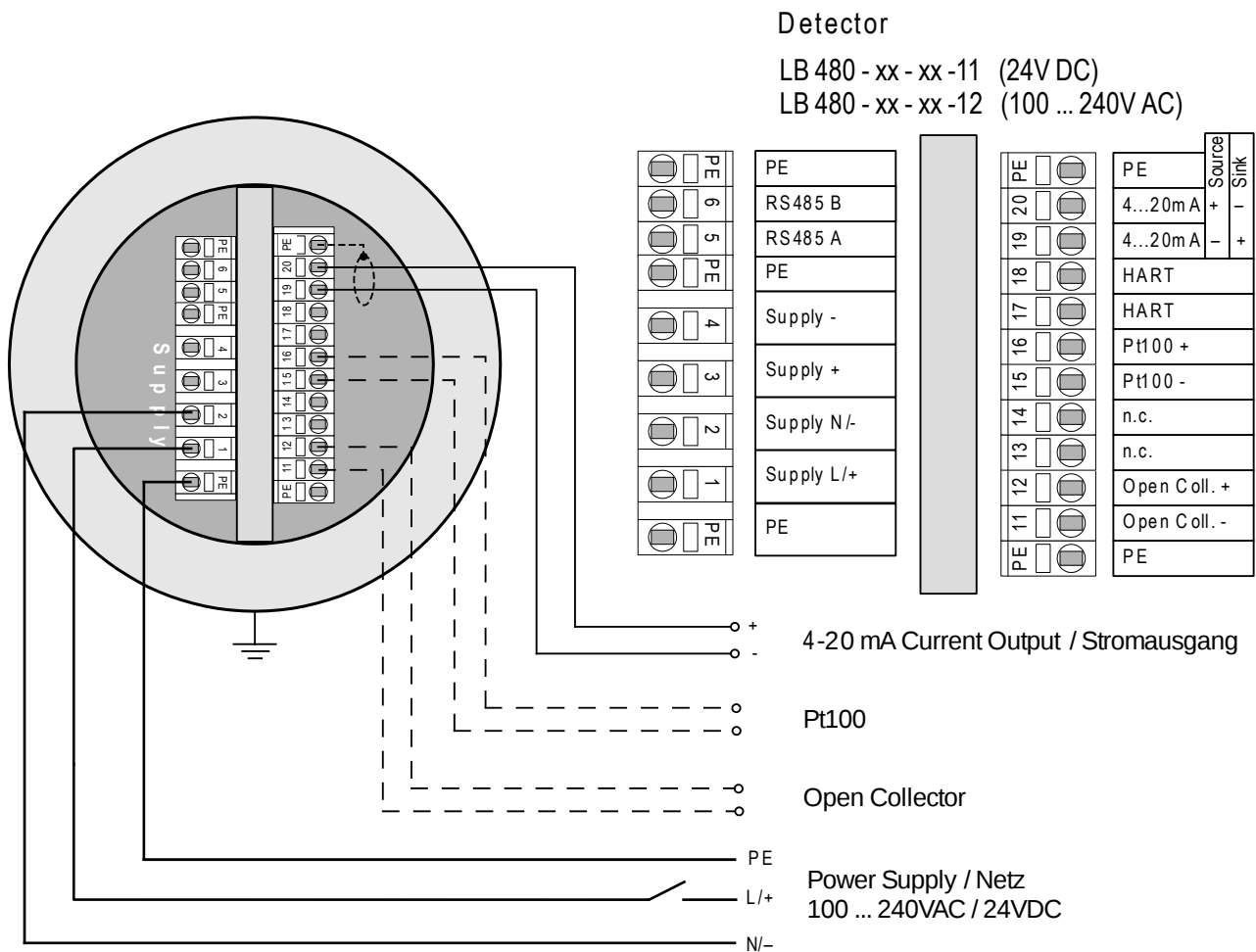
plug the latch into the catch bolt
Klinke in Kupplung einklinken



latch plugged into the catch bolt
Klinke in Kupplung einklinken



9.0 Terminal Connection Elektrischer Anschlussplan



LB 480

Dimensions in mm
Abmessungen in mm

[illegible]