

# POINT SOURCE SHIELDS LB 744x FOR DENSITY AND LEVEL SWITCH

## Reliable & safe

Complies with international standards for constructional requirements and classification of radiometric gauges.

## Corrosion resistant

Available in stainless steel (Super Duplex)

## Fire resistant

Fireproof version available

## Flexible

Available in three different sizes for a wide range of activities.



# TECHNICAL DATA & FACTS

## Point source shields LB 744x

Shields	LB 7440	LB 7442	LB 7444
Application	Density, Level switch	Density, Level switch	Density, Level switch
Nuclide	Punktstrahler	Punktstrahler	Punktstrahler
Shielding material	Lead	Lead	Lead
Lead thickness	67 mm	94 mm	132 mm
Exit angle	9°, 14 °	6°, 10°	6°
Beam geometry	Collimated round	Collimated round	Collimated round
Outer material	Painted steel, stainless steel 1.4410 (Super Duplex)	Painted steel, stainless steel 1.4410 (Super Duplex)	Painted steel
Fire resistance class acc. to IEC 62598:2011	Class F (optional)	Class F (optional)	N
Mounting	Flange, socket	Flange, socket	Flange
Weight	32 kg	85 kg	177 kg
Pneumatic shutter control possible	Yes	Yes	Yes
Ambient, storage and transport temperature	-40 ... +100 °C	-40 ... +100 °C	-40 ... +100 °C

### Special Features

- Various exit angles: suitable for your application
- Various sizes gradations: for compliance with country-specific radiation protection requirements for low and very high source activities
- Outer material: painted steel or various grades of stainless steel available for use under demanding environmental conditions
- Nuclides: Cs-137 and Co-60
- Remote-controlled pneumatically-driven shutter: Can be fitted to all variants