

ROTATABLE POINT SOURCE SHIELD LB 8301 FOR DENSITY AND BULK FLOW

Reliable & safe

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

Corrosion resistant

Available in stainless steel

Flexible

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



TECHNICAL DATA & FACTS

Point source shields LB 8301

Shields	LB 8301, Type 100	LB 8301, Type 150	LB 8301, Type 200	LB 8301, Type 270
Application	Level, Bulk flow	Level, Bulk flow	Level, Bulk flow	Level, Bulk flow
Nuclide	Cs-137, Co-60	Cs-137, Co-60	Cs-137, Co-60	Cs-137, Co-60
Shielding material	Lead	Lead	Lead	Lead
Lead thickness	Approx. 42 mm	Approx. 67 mm	Approx. 94 mm	Approx. 130 mm
Exit angle	Standard is 45°, other angles upon request	Standard is 45°, other angles on request	Standard is 45°, other angles on request	Standard is 45°, other angles on request
Beam geometry	Fan-shaped collimated according to the beam exit angle	Fan-shaped collimated according to the beam exit angle	Fan-shaped collimated according to the beam exit angle	Fan-shaped collimated according to the beam exit angle
Outer material	Painted steel, stainless steel 1.4310, stainless steel 1.4571	Painted steel, stainless steel 1.4310, stainless steel 1.4571	Painted steel, stainless steel 1.4310, stainless steel 1.4404 (Super Duplex)	Painted steel, stainless steel 1.4404 (Super Duplex)
Fire resistance class acc. to IEC 62598:2011	K	K	K	K
Mounting	Base mounting / bulk flow measuring frame	Base mounting / bulk flow measuring frame	Base mounting / bulk flow measuring frame	Base mounting
Weight	Approx. 32 kg	Approx. 70 kg	Approx. 150 kg	Approx. 370 kg
Pneumatic shutter control possible	Yes	Yes	Yes	Yes
Ambient, storage and transport temperature	−40 ... +100 °C	−40 ... +100 °C	−40 ... +100 °C	−40 ... +100 °C

Special Features

- Multiple exit angles: suitable for your application
- Various sizes: for compliance with country-specific radiation protection requirements for low to very high source activities
- Outer material: painted steel or various grades of stainless steel available for use in demanding environments
- Nuclides: Cs-137 and Co-60 point sources can be used
- Remote-controlled pneumatically-driven shutter: can be fitted to all variants if required