

CONGAUGE LB 6755

MOULD LEVEL DETECTOR FOR STEEL & POWDER

Backwards compatible

Mechanically backward compatible with LB 6752 / LB 6652 detectors.

Superior redundancy

Automatic detection and compensation if a scintillation failure should occur.

Fast response time

Fast changes of the level can be detected due to a cycle time of only 5 ms and 500 ms for powder level.

Compatible with electromagnetic stirrers

Specially designed for use in the direct environment of electro-magnetic stirrers and brakes.



in cooperation with

SMS  **group**
SMS Concast AG



TECHNICAL DATA & FACTS

CONGAUGE LB 6755

Detector

Application	Mould level (steel and powder level simultaneously)
Power supply	15...32 V DC, approximately 6 W
Temperature range	-20 ... +55 °C (-4...131 °F) operation -20 ... +70 °C (-4...158 °F) storage
Temperature stability	≤0.2 %/°C (-20 ... +55 °C)
Housing material	Stainless steel ISO 1.4301 / AISI 304
IP protection	IP66/67
Scintillator	10 x NaI(Tl)-crystals; 40 mm x 45 mm x 14 mm
Photomultiplier	Silicon based photomultiplier arrays
Dimensions	Variant 1: ~100 x 200 x 157 mm (W x H x L) Variant 2: ~100 x 200 x 280 mm (W x H x L)
Weight	Variant 1: ~10 kg Variant 2: ~13 kg
Water connection	R ¼ with maximum pressure of 6 bar Water of drinking quality
Electrical connection	GAMMAcast male connector
Communication	RS485 and analog pulses to the castXpert LB 452 Control unit for redundancy
Internal cycle time	5 ms standard steel level 500 ms for powder level
Redundancy	A scintillation module failure is automatically detected and compensated for
Electromagnetic flux tolerance	Handles up to 150 mT without signal degradation

Special Features

- Measures the steel level and the powder level with one technology simultaneously and accurately
- For use with electrometric stirrer and brakes without signal deterioration
- Industry-leading measurement cycle time
- Industry-leading temperature stability and mechanical robustness
- Available in two variants, one of which is fully mechanically compatible with the LB 6752 / LB 6652 detector
- Superior safety with automatic detection and compensation if a scintillation failure should occur
- Service and extensive spare parts program available