

GAMMAcast LB 6752

MOULD LEVEL DETECTOR

Fast response time

Fast changes of the level can be detected safely and reliably due to a cycle time of only 5 ms.

Microprocessor controlled

Constant monitoring and control of the detector status.

High sensitivity

Using our largest scintillator.

Temperature stable

Stable and reliable measurement even at high temperatures and significant temperature fluctuation.

Robust & water resistant

IP67, designed for harsh operating conditions such as high temperatures, vibration, shock, and water. Maximum robustness against mechanical stress.



TECHNICAL DATA & FACTS

GAMMAcast LB 6752

Detector

Application	Mould level
Power supply	5 ... 32 VDC, approximately 1.2 W
Temperature range	-20 ... +50 °C operation -20 ... +70 °C storage
Temperature stability	≤ 0.2 %/°C (-20 ... +50 °C)
Housing material	Stainless steel ISO 1.4301/AISI 304
IP protection	IP 67
Scintillator	Organic polymer 150 x 100 x 50 mm (W x H x L) or 150 x 100 x 25 mm (W x H x L)
Photomultiplier	2" Vacuum photomultiplier
Dimensions	~100 x 200 x 280 mm (W x H x L)
Weight	~19 kg
Water connection (optional)	R¼ with maximum pressure of 6 bar Water of drinking quality
Electrical connection	GAMMAcast female connector
Communication	RS485 and analog pulses to the castXpert LB 452 Control unit for redundancy
Electromagnetic flux tolerance	Handles up to 2 mT without signal degradation

Special Features

- Our most sensitive mould level detector
- Industry-leading measurement cycle time
- Industry-leading temperature stability and mechanical robustness
- Detector self-diagnostics and validation of communication
- Service and extensive spare parts program available