

GAMMAcast LB 6739 / 6760

MOULD LEVEL DETECTORS

SiPM Technology

The world's first and most robust, mould level detector based on Silicon Photomultiplier (SiPM) technology. (LB 6760)

Cooling option

Available with water cooling.

Robust & water resistant

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

Microprocessor controlled

Constant monitoring and control of the detector status.

Compatible with electromagnetic stirrers

Specially designed for use in the direct environment of electromagnetic stirrers and brakes. (LB 6760)

Long established system

The industry standard and our best-selling mould level detector with well over 3,000 strands in operation worldwide.

Fast response time

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

Temperature stability

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



TECHNICAL DATA & FACTS

GAMMAcast LB 6739 / 6760

Detector	LB 6739	LB 6760
Application	Mould level	Mould level
Power supply	5 ... 32 V DC, approximately 1.2 W	5 ... 32 V DC, approximately 1.2 W
Temperature range	-20 ... +50 °C operation -20 ... +70 °C storage	-20 ... +60 °C operation -20 ... +70 °C storage
Temperature stability	≤ 0.002 %/°C (-20 ... +50 °C)	≤ 0.002 %/°C (-20 ... +60 °C)
Housing material	Stainless steel ISO 1.4301/AISI 304	Stainless steel ISO 1.4301/AISI 304
IP protection	IP 67	IP 67
Scintillator	CsI crystal; Ø 40 mm or Ø 25 mm, H: 50 mm	Nal crystal; Ø 40 mm x 50 mm
Photomultiplier	1" Vacuum photomultiplier	Silicone based photomultiplier array
Dimensions	Ø 50 mm (Ø 60 mm with water cooling) 327 mm length	Ø 50 mm (Ø 60 mm with water cooling) 327 mm or 264 mm length
Weight	~2 kg and ~3 kg with water cooling	~2 kg and ~3 kg with water cooling
Water connection (optional)	R¼ with maximum pressure of 6 bar Water of drinking quality	R¼ with maximum pressure of 6 bar Water of drinking quality
Electrical connection	GAMMAcast female connector	GAMMAcast female connector
Communication	RS485 and analog pulses to the castXpert LB 452 Control unit for redundancy	RS485 and analog pulses to the castXpert LB 452 Control unit for redundancy
Electromagnetic flux tolerance	Handles up to 2 mT without signal degradation	Tested up to to 150 mT without signal degradation

Special Features

- For use with electrometric stirrer and brakes without signal deterioration (LB 6760)
- Automatic voltage control
- Industry-leading measurement cycle time
- Industry-leading temperature stability and mechanical robustness
- Detector self-diagnostics and communications verification
- Service and extensive spare parts program available