

GAMMAcast Cable FOR MOULD LEVEL MEASUREMENT

Secure & reliable

Proven, screw-locked industrial stainless-steel plug.

Universal

Compatible with all GAMMAcast detectors.

Flexible

Available in four different lengths and with straight or angled connector.

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 cable

Complete cable assembly

Application	Mould level
Temperature range	–20 ... +125 °C operation –20 ... + 70 °C storage
IP protection	IP67
Length	Available in 5, 10, 15, and 20 meters
Detector connection	Standard GAMMAcast female connector – straight or angled
Junction box connection	Open ends or industrial connector

Protective hose

Length	5 m for all cable lengths
Material	Silicone coated fiberglass
Robustness	Handles splashes of liquid steel and iron
Temperature range	–55 ... +260 °C operation +1090 °C for 20 minutes +1640 °C for 15 – 30 seconds

Electrical cable

Length	Available in 5, 10, 15, and 20 meters
Material	Silicone coated
Diameter	13 mm
Temperature range	–25 ... +180 °C operation
Minimum bend radius	~ 130 mm dynamic installation ~ 80 mm static installation

GAMMAcast connector

Material	Stainless steel
Temperature range	–20 ... +125 °C operation
Variants	Straight or angled

Special Features

- Strain relief spring
- Stainless-steel screw connection on detector side
- Backward compatible with existing detectors and junction boxes
- Developed in cooperation with our customers