

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.

Durable

Two options for protective sleeve: standard and heavy duty.

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 to +125 °C operation –20 to + 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

| | Red cable | Black cable |
|-------------------|---|--|
| Length | 5 m or according to cable length | 5 m or according to cable length |
| Material | Silicone coated fiberglass | Galvanized steel core with plastic sheath |
| Robustness | Handles splashes of liquid steel and iron | Additional mechanical robustness in comparison to red hose |
| Temperature range | –55 to +260 °C operation +1090 °C for 20 minutes +1640 °C for 15 – 30 seconds | –60 to +150 °C for black plastic sheath; melting temperature of galvanized steel core: 1400 °C |

Electrical cable

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

GAMMAcast connector

| | |
|-------------------|--------------------------|
| Material | Stainless steel |
| Temperature range | –20 to +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