

castXpert LB 452

MOULD LEVEL CONTROL UNIT

Highest measurability & repeatability

Built-in filter for suppression of mould oscillation. Display of measurement accuracy in the system.

Flexibility

Quick and easy format change by storing several calibration curves.

User-friendly

Intuitive and multilingual user interface via touchscreen.

Safety

Independent measuring channels with separate power supply.



Modularly expandable

Expansion to up to 4 measuring channels and additional signal input and output cards possible.

Tried and tested

Top customer satisfaction with over 1,000 systems in use worldwide.

TECHNICAL DATA & FACTS

castXpert LB 452

Control unit (handles up to 4 active channels)

Application	Mould level
Dimension	19" Rack, 3 HE (7" colour touch screen)
Power supply	90 ... 264 VAC, 50/60 Hz
Temperature range	-20 ... +50 °C operation -20 ... +70 °C storage
IP protection	IP 20
Weight	4 – 6 kg depending on number of boards installed
Connection	USB port for of USB storage devices or external keyboard or mouse Ethernet for remote control and logging

Measurement board (required per active channel)

Each active channel works separately with own power supply unit and connections.

Power supply	90 ... 264 VAC, 50/60 Hz
Inputs	Detector connection 2 digital inputs for remote empty and full adjustment
Outputs	1x 4 ... 20 mA current output for mould level Relay output for error signal
Configuration	Stores up to 8 calibration curves with up to 21 calibration points each

Extension boards (optional per active channel)

Inputs	2 digital inputs for remote selection of up to 4 calibration curves
Outputs	1x 4 ... 20 mA current output for mould level, detector temperature, or detector counts 2 relay outputs for alarm signal

Bus boards (optional per active channel)

Interfaces	Profibus DP or ProfiNet for cyclic data output and system selection
------------	---