

PMA.Evolution (model EV01Y1)

For use in zone 2 hazardous areas

Paint Mixing Scales



- One cable for power and data
- Integrated mixing module processes one formulation
- New high contrast display technology
- Glass sensor touch technology
- Multi-lock security function

Technical data

Capacity	7500 999,95 g
Readability	0,1 0,05 g
Tare range (subtractive)	-7500 g
Weighing pan (∅)	233 mm
Housing (WxDxH)	233 x 321 x 341 mm
Net weight	approx. 2.4 kg
External calibration weight	1, 2, 5 kg (class F2 or better)
Allowable operating temperature	5 ... 40 °C
Dust and water protection rating in accordance to EN 60529/IEC 60529:	IP40
Power supply	via USB 2.0 or higher or power supply YPS06-USB
Input voltage power consumption	5 V _{DC} 2 W
Approval	Ex zone 2
IECEX	Ex ic nA IIB T4 Gc
ATEX	II 3G Ex ic nA IIB T4 Gc
Data output	USB Type B
Format	7-bit ASCII, 1 Start-bit, 1 or 2 Stop-Bits
Parity	Even, odd, no parity
Transmission rates	1200 to 57,600 Bit/s
Handshake mode	Software, Hardware or None
Internal application	Integrated mixing module for one job Recalculation (correction of over-pours) Calculation by factor (0.25 – 5.0) Stand-alone recipe processing (absolute, additive) Dual range 0.05 g (up to 999.95 g) Hardware test and communication symbol

Standard accessories

USB cable, length 5 m (16.4 ft.)	YCC01-0040M5
----------------------------------	--------------

Optional accessories

In-use cover for display (set of 10)	YDC03PMA10
In-use cover for column (set of 10)	YDC03PMA-CO10
In-use cover for weighing pan (set of 10)	YDC03PMA-WP10

USB power plug 5V / 900 mA	YPS06-USB
AC-adapter set (US, EU, UK)	YAK01
AC- adapter set (AU, ZA, BR)	YAK02
AC- adapter set (IN, KR, CN)	YAK03

External calibration weight 1 kg (class F2)	YCW614-AC-00
External calibration weight 1 kg (class F2) with DAKKS certificate	YCW614-AC-02
External calibration weight 2 kg (class F2)	YCW624-AC-00
External calibration weight 2 kg (class F2) with DAKKS certificate	YCW624-AC-02
External calibration weight 5 kg (class F2)	YCW654-AC-00
External calibration weight 5 kg (class F2) with DAKKS certificate	YCW654-AC-02



EV01Y1 rear view

Dimension drawing

