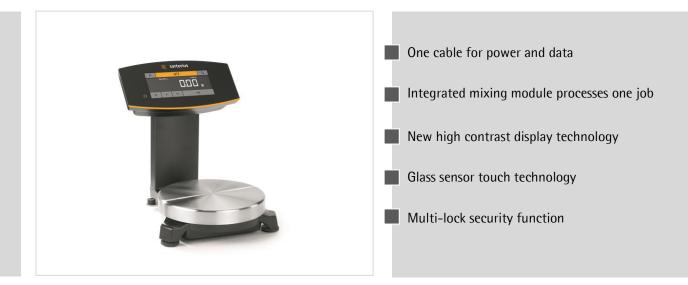


PMA.HD (model SP01S1)

For use in non-hazardous areas



Technical data

Canadity	2200 g
Capacity	5
Readability	0.01 g
Tare range (subtractive)	-2200 g
Weighing pan (Ø)	180 mm
Housing (WxDxH)	212 x 294 x 261 mm
Net weight	approx. 2.1 kg
External calibration weight	1, 2 kg (class F1 or better)
Allowable operating temperature	5 40 °C (41 104 °F)
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
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)
	Hardware test and communication symbol

Standard accessories

YCC01-0040M5

Optional accessories

In-use cover for display (set of 10)	YDC03PMA10
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 F1)	YCW613-AC-00
External calibration weight 1 kg (class F1) with DAKKS certificate	YCW613-AC-02
External calibration weight 2 kg (class F1)	YCW623-AC-00
External calibration weight 2 kg (class F1) with DAKKS certificate	YCW623-AC-02



SP01S1 rear view

Dimension drawing

