# SARTURIUS

# PMA HD

Model SPO1S1 For Use in Non-hazardous Areas



### Benefits

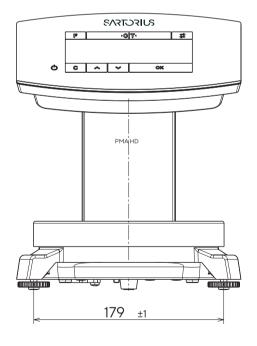
- One cable for power and data
- Integrated mixing module processes one job
- New high contrast display technology
- Glass sensor touch technology
- Multi-lock security function

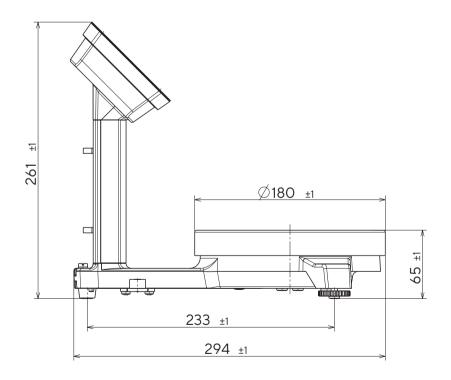
## Technical Specifications

Capacity	2,200 g	
Readability	0.01 g	
Tare range (subtractive)	-2,200 g	
Weighing pan (Ø)	180 mm	
Housing (W×D×H)	212×294×261 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 YEPS01-USB	

Input voltage   power consumption	5 V <sub>DC</sub>   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	1,200 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	

## Technical Drawings





### Accessories

### **Standard Accessories**

### Optional Accessories

In-use cover for display (set of 10)	YDC03PMA10
In-use cover for column (set of 10)	YDC04PMA-C010
In-use cover for weighing pan (set of 10)	YDC04PMA-WP10
USB power plug 5.2 V <sub>DC</sub>   1400 mA	YEPS01-USB
AC-adapter set (US, EU, UK)	YEPS01-PS1
AC-adapter set (AU, ZA, BR, AR)	YEPS01-PS6
AC-adapter set (IN, KR, CN)	YEPS01-PS7
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



SPO1S1 rear view

### Germany

Sartorius Lab Instruments GmbH & Co. KG Otto-Brenner-Strasse 20 37079 Goettingen Phone +49 551 308 0

⊕ For further information, visit

www.sartorius.com

#### **USA**

Sartorius Corporation 565 Johnson Avenue Bohemia, NY 11716 Phone +1 631 254 4249 Toll-free +1 800 635 2906