

SARTORIUS

PMA Power

Awesome Accuracy up to 35 kg

Your Advantages:

- Bench or wall mount flexibility
- Built-In overpour recalculation feature
- Fractional mixing for different can sizes
- Solvent-resistant housing components
- Standard RS-232 interface port with flexible connection options
- Multi-language menu



Learn More:



PMA Power



PMA 35001-X

The PMA Power is the ideal solution if you don't want to make any compromises when mixing large amounts of paint. Its monolithic weigh cell featuring a resolution of 0.1 g and a weighing capacity extending up to 35 kg enables you to pour components quickly and accurately into relatively large paint cans.

Thanks to its high resolution, you will also find it easy to pour in small amounts with total accuracy on the PMA Power scale. This lets you save money by not having to buy a second scale.

The PMA Power comes standard with an interface port. This is how it works:

- As a stand-alone scale
- Connected to a PC
- In conjunction with paint-mixing systems

Its enhanced, large 14-segment display enables the operator to clearly read both weights and color components as well as other important information.

In addition, the PMA Power display is equipped with a bracket that can be installed on a wall or set up on a bench. This gives you optimal flexibility in customizing your work area to suit your needs. It also ensures that you will have all important information available at a glance!

The special display lock function designed by Sartorius enables you to secure your PMA Power against unintentional use and hide weights on its display very easily to prevent unauthorized paint mixing.

Description	Order Number
PMA Power for use in safe areas	PMA35001
PMA Power for use in zone 2 hazardous areas	PMA35001-Y
PMA Power for use in zone 1 hazardous areas (Class I, Div 1) (main adapter not included with the equipment supplied)	PMA35001-X
Accessories for connection to a PC	
RS232/USB cable 25pin to USB	YCC01-USBM2
Adapter cable 25pin to 9pin	69PM0057
RS-232 data cable (2 m)	YCC01-0047M2
USB/RS-232 converter cable (1.8 m)	YCO12
Technical Specifications	
Weighing capacity	35,000 g
Readability	0.1 g
Load plate	350 x 240 mm
Net weight	11.4 kg
Accessories	
Dust cover for display unit	YDC01PMA
External calibration weight 10 kg (class F2)	YCW7148-00
External calibration weight 10 kg (class F2)with DAKKS certificate	YCW7148-02

For further technical specifications and information on approvals and certifications, please refer to our data sheets.

Germany

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

USA

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

 For further contacts, visit
www.sartorius.com