



PMA.Quality
For easy interactive
communication

The new PMA.Quality series

Rugged, reliable and full-featured



The new PMA.Quality paint-mixing scale from Sartorius sets new standards in operation, features and applications. Whether used as a stand-alone scale or connected to a PC or mixing system, the PMA.Quality has the winning features that only a genuine Sartorius paint-mixing scale can offer.



Its enhanced, large 14-segment display provides outstanding viewing quality for showing weights, color components and other information important to the body shop specialist. Even relatively long texts are easy and convenient to read thanks to the scale's scrolling text function.



The display unit is backlit and tiltable, allowing easy adjustment to your viewing angle and thus ensuring excellent readouts even under poor lighting conditions. In addition, you can set up the display unit separately as dictated by space requirements.



The keypad overlay features a completely new design, and the operator is guided intuitively by icons familiar from computer technology. Large keys with positive click action let the user know when a key is pressed.

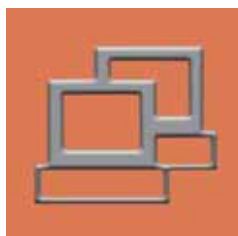
Did you under- or over-pour a component? No problem! With its recalculation feature, the PMA.Quality will quickly and accurately calculate the exact quantities that you need to add so that your formula will be accurate – to put you on the safe side. You can just as easily use this built-in formula calculator to calculate quantities for paint can sizes that differ from the standard 1-liter size and conveniently pour in the individual components.

On the PMA.Quality, the Dual-Range function enables you to simply increase the reading accuracy to 0.05 g – over a weighing range of 1,000 g – which is especially convenient when working with the smallest paint amounts.

The overload and side impact protection makes the weigh cell exceptionally resistant to extremely tough environmental conditions. Of course, all parts of the housing are resistant to solvents and are therefore easy to clean. The wrap-around design of the stainless steel pan protects the weigh cell from dirt build-up and paint spillage.

Thanks to its intrinsically safe design, the explosion-protected version of the new PMA.Quality can be safely operated in zone 1 hazardous areas and Class I Division 1 locations without any added and difficult-to-install safety systems. Just like many of the current Sartorius paint-mixing products, the PMA.Quality can also be safely used in flammable environments (according to HD 384.4.482 S1, TÜV NORD CERT).





If the scale is in the interactive communication mode, this is conveniently displayed for the user by the communication icon.



With the display lock function developed by Sartorius, the PMA.Quality scale can be locked to prevent unauthorized use.



Uncompromisingly excellent

- : Large 14-segment display
- : New keypad overlay design
- : Scrolling text function
- : Programmable information displayed during power-up
- : Recalculation function
- : Built-in formula calculator
- : Dual-range scale
- : Display lock function
- : Communication symbol
- : Detachable display unit
- : Housing parts resistant to solvents
- : Serial RS-232C interface port
- : Multi-language setup menu
- : Service available worldwide
- : May be used in flammable environments
- : USB adapter cable (optional)

Product	Order Number
PMA.Quality for zone 1, Class I Division 1	PMA7501-X

Accessory:

In-use cover	YDC01PMA
--------------	----------

Connection to TopMix|EcoMix:

RS-232C data cable (3 m)	YCC01-0041M3
--------------------------	--------------

Technical Specifications:

Weighing capacity ranges	999.95 g/7,500 g
Readability	0.05 g/0.1 g
Weighing pan diameter	233 mm
Net weight	3.2 kg

The PMA7501-X... can be installed in zone 1 hazardous areas.

Approval Information

The PMA7501-X has been tested and approved for use in zone 1 hazardous areas. It was tested by KEMA Quality B.V. according to the harmonized European Standards EN 50014 and EN 50020.

The European Standards previously mentioned are equivalent to Factory Mutual (FM) Approval Standards Class 3600 and Class 3610 as well as the CSA Approval Standards C22.2 No. 0, C22.2 No. 142; and C22.2 No. 157, E60079-0 and E60079-11, according to which this equipment has been tested and approved for the North American market.

You can request these certificates from Sartorius AG, Goettingen, Germany.

Sartorius AG
Weender Landstrasse 94-108
37075 Goettingen, Germany

Phone +49.551.308.3234
Fax +49.551.308.3612

www.sartorius.com