

PMA.Evolution – The New PMA Standard Economic. Efficient. Effective.



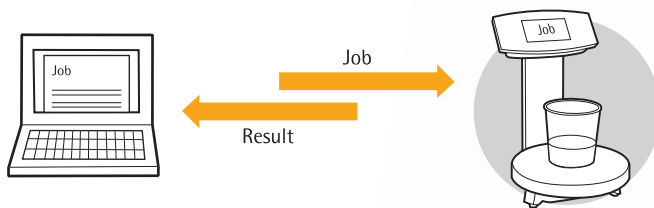
turning science **into solutions**

Efficient and Easy Paint Mixing

Increase your paint mixing efficiency while reducing your total cost of ownership: Benefit from the precise, reliable and fast results of the new PMA.Evolution. Its superior readability – both from an angle and from a distance – make it a practical and time-saving tool to work with. With a variety of accessories you can enhance the functionality and usability of this high-quality Made-in-Germany product.

Internal Mixing Module

- Integrated mixing module for PC-independent formulation processing



Integrated Applications

- Recalculation functionality in case of over-pouring a component
- Factor calculation to adapt a recipe to any desired can size
- The scale can be calibrated externally with 1 kg, 2 kg or 5 kg.
- The resolution of 0.1 g can be manually switched to 0.05 g for a range of up to 999.95 g. The fine range can be used up to the maximum capacity through taring.

Clean Work Environment



- Plane, solvent-resistant glass display and stainless steel weighing pan for easy cleaning



- Spillage-protected interface at the back of the head

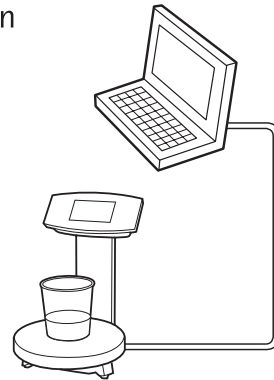


- In-use covers for display, column and weighing pan to prevent stains



Amazingly Fast Installation

- Unique one-cable solution for power and data via USB cable
- USB driver is Microsoft® certified for "plug & play" installation



New High-contrast Display



- Wide viewing angle and perfect readability, even from a distance



- Sensor touch with visual feedback



- Bargraph with zoom functionality as dosing assistance | capacity indicator

Enhanced Safety Mechanisms ▷

Compliance with International Regulations ▷

Exclusive PMA.Management Software ▷

Extended Features

Prevention of Unauthorized Usage



- Instant lock function: activated manually by pressing the lock symbol
- Communication lock function: automatic display lock if there is no ongoing communication between scale and PC



- Theft protection with fixation point on the back



Compliance with International Regulations

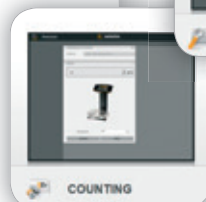
- Accordance with legal regulations
- Certified as explosion proof



Exclusive PMA.Management Software



- Scale settings easily adjustable with intuitive menu
- Intro text for individualized scale messages (e.g. branding)
- Basic counting application, e.g. for inventory control





sartorius

F

·0|T·

↔

Max 1500 g

d = 0.1 g

0.0 g



C



OK



PMA.Evolution

Functional Accessories

Keep your PMA.Evolution working like new for the longest time:
with covers to protect your scale from paint stains and weights to recalibrate it.
See what other useful accessories might be convenient for you.

In-Use covers

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



Cables

Link cable to converter, length 10 m (32.8 ft.)	YCC01-0052M10
Link cable to converter, length 20 m (65.6 ft.)	YCC01-0052M20
Link cable to converter, length 30 m (98.4 ft.)	YCC01-0052M30
Grounding cable, 2 m (6.5 ft.)	YCC01-X046M2
USB cable, length 5 m (16.4 ft.)	YCC01-0040M5



Power supply

USB power plug 5 VDC 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



Calibration weights

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



Technical Data

Metrological data

Capacity	7500 g
Readability	0.1 g
Tare range (subtractive)	-7500 g
Allowable operating temperature	5 ... 40 °C (41 ... 104 °F)
External calibration weight	1, 2, 5 kg (class F2 or better)

Dimensions | Weight

Weighing pan (Ø)	233 mm
Housing (W×D×H)	233 × 321 × 341 mm
Net weight	approx. 2.4 kg

Power supply

Power supply	via USB 2.0 or higher or power plug YPS06-USB
Input voltage power consumption	5 VDC 2 W
Ex-link converter (Ex zone 1 scale only)	Energy limiting unit included

Approval Ex zone 1 | Class I, Div. 1

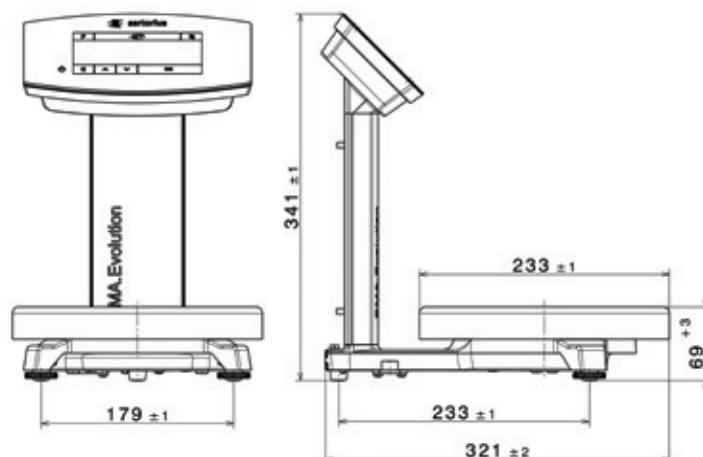
IECEX	Ex ib IIB T4 Gb
ATEX	II 2G Ex ib IIB T4 Gb
FM (USA)	IS CL I, DIV 1, GP C,D T4; CL I, Zone 1, AEx ib IIB T4
FM (CAN)	Ex ia CL I, DIV 1, GP C,D T4; Zone 1, Ex ib IIB T4

Approval Ex zone 2

IECEX	Ex ic nA IIB T4 Gc
ATEX	II 3G Ex ic nA IIB T4 Gc

IP protection

Dust and water protection rating in accordance to EN 60529 IEC 60529:	IP40
---	------



Communication

Data output	USB Type B
Protocol	Sartorius SBI XBPI
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
----------------------	--

User interface	High contrast display 14 segment Bargraph as dosing assistance or capacity indicator Text menu in 12 languages
----------------	--

Sales and Service Contacts

For further contacts, visit www.sartorius.com

Europe

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Weender Landstrasse 94-108
37075 Goettingen

Phone +49.551.308.0
Fax +49.551.308.3289

France & Suisse Romande

Sartorius France
2, rue Antoine Laurent de Lavoisier
ZA de la Gaudrée
91410 Dourdan

Phone +33.1.70.62.50.00
Fax +33.1.64.59.76.39

Austria

Sartorius Austria GmbH
Modecenterstrasse 22
1030 Vienna

Phone +43.1.7965760.0
Fax +43.1.7965760.24

Belgium

Sartorius Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles

Phone +32.2.756.06.90
Fax +32.2.481.84.11

Finland & Baltics

Sartorius Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki

Phone +358.9.755.951
Fax +358.9.755.95.200

Hungary

Sartorius Hungária Kft.
Kagyó u. 5.
2092 Budakeszi

Phone +3623.457.227
Fax +3623.457.147

Ireland

Sartorius Ireland Ltd.
Unit 41, The Business Centre
Stadium Business Park
Ballycoolin Road
Dublin 11

Phone +353.1.8089050
Fax +353.1.8089388

Italy

Sartorius Italy S.r.l.
Viale A. Casati, 4
20835 Muggiò (MB)

Phone +39.039.4659.1
Fax +39.039.4659.88

Netherlands

Sartorius Netherlands B.V.

Phone +31.30.60.53.001
Fax +31.30.60.52.917
info.netherlands@sartorius.com

Poland

Sartorius Poland sp.z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn

Phone +48.61.6473830
Fax +48.61.6473839

Russian Federation

LLC "Sartorius ICR" and LLC "Biohit"
Uralskaya str. 4, Lit. B
199155, Saint-Petersburg

Phone +7.812.327.5.327
Fax +7.812.327.5.323

Spain & Portugal

Sartorius Spain, S.A.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)

Phone Spain +34.902.123.367
Phone Portugal +351.800.855.800
Fax Spain +34.91.358.96.23
Fax Portugal +351.800.855.799

Switzerland

Sartorius Mechatronics Switzerland AG
Ringstrasse 24a
8317 Tagelswangen (ZH)

Phone +41.44.746.50.00
Fax +41.44.746.50.50

U.K.

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9QQ

Phone +44.1372.737159
Fax +44.1372.726171

Ukraine

LLC "Biohit"
Post Box 440 "B"
01001 Kiev, Ukraine
Phone +380.44.411.4918
Fax +380.50.623.3162

Americas

USA

Sartorius Corporation
5 Orville Drive, Suite 200
Bohemia, NY 11716

Phone +1.631.254.4249
Toll-free +1.800.635.2906
Fax +1.631.254.4253

Argentina

Sartorius Argentina S.A.
Int. A. Ávalos 4251
B1605ECS Munro
Buenos Aires

Phone +54.11.4721.0505
Fax +54.11.4762.2333

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900
Fax +55.11.4362.8901

Canada

Sartorius Canada Inc.
2179 Dunwin Drive #4
Mississauga, ON L5L 1X2

Phone +1.905.569.7977
Toll-Free +1.800.668.4234
Fax +1.905.569.7021

Mexico

Sartorius de México S.A. de C.V.
Circuito Circunvalación Poniente
No. 149
Ciudad Satélite
53100, Estado de México
México

Phone +52.5555.62.1102
Fax +52.5555.62.2942

Asia | Pacific

Australia

Sartorius Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175

Phone +61.3.8762.1800
Fax +61.3.8762.1828

China

Sartorius Scientific
Instruments (Beijing) Co., Ltd.
33 Yu An Road, Airport Industrial Park
Zone B,
Shunyi District, Beijing 101300,
P.R.China

Phone +86.10.8042.6300
Fax +86.10.8042.6486

Hong Kong

Sartorius Hong Kong Ltd.
Unit 1012, Lu Plaza
2 Wing Yip Street
Kwun Tong
Kowloon, Hong Kong

Phone +852.2774.2678
Fax +852.2766.3526

India

Sartorius Weighing India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra,
Nelamangala Tq
562 123 Bangalore, India

Phone +91.80.4350.5250
Fax +91.80.4350.5253

Japan

Sartorius Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.3740.5408
Fax +81.3.3740.5406

Malaysia

Sartorius Malaysia Sdn. Bhd
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622
Fax +60.3.8996.0755

Singapore

Sartorius Singapore Pte. Ltd
1 Science Park Road,
The Capricorn, #05-08A,
Singapore Science Park II
Singapore 117528

Phone +65.6872.3966
Fax +65.6778.2494

South Korea

Sartorius Korea Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, Bundang-Gu
SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700
Fax +82.31.622.5799

Thailand

Sartorius (Thailand) Co. Ltd.
129 Rama 9 Road,
Huaykwang
Bangkok 10310

Phone +66.2643.8361-6
Fax +66.2643.8367



◀ www.sartorius.com/paintmixing