

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

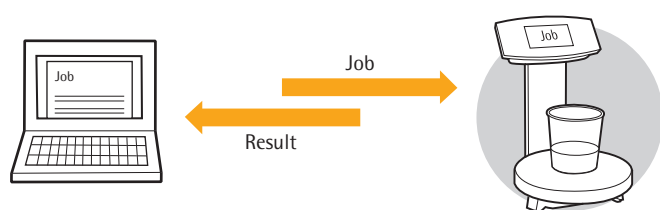


Efficient and Easy Paint Mixing

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



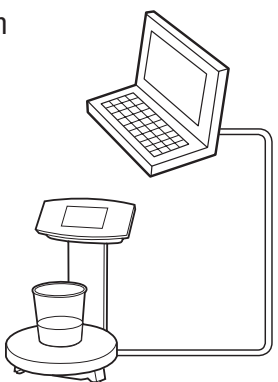
- | Spillage-protected interface at the back of the head



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

Amazingly Fast Installation

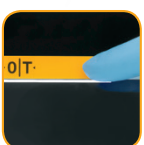
- | Unique one-cable solution for power and data via USB cable
- | USB driver is Microsoft® certified for "plug & play" installation



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

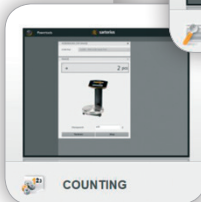
- | Accordance with legal regulations
- | Certified as explosion proof



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



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 5.2 VDC 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



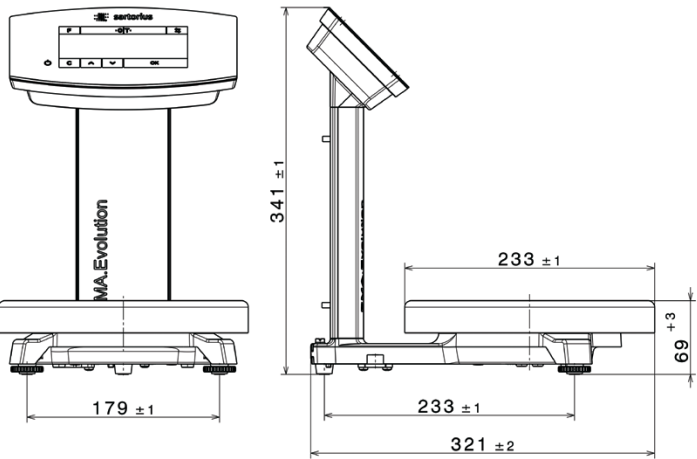
Calibration weights	
External calibration weight 1 kg (class F2)	YCW614-AC-00
External calibration weight 1 kg (class F2) with DAKKS certificate	YCW614-AC-02
External calibration weight 2 kg (class F2)	YCW624-AC-00
External calibration weight 2 kg (class F2) with DAKKS certificate	YCW624-AC-02
External 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 (∅)	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
PCEC (CN)	EX ib IIB T4 Gb

Approval Ex zone 2	
IECEx	Ex ic nA IIB T4 Gc
ATEX	II 3G Ex ic nA IIB T4 Gc
PCEC (CN)	Ex ic nA IIB T4 GC
DEKRA (JP)	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.
Kagyló 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. Avalos 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