



sartorius

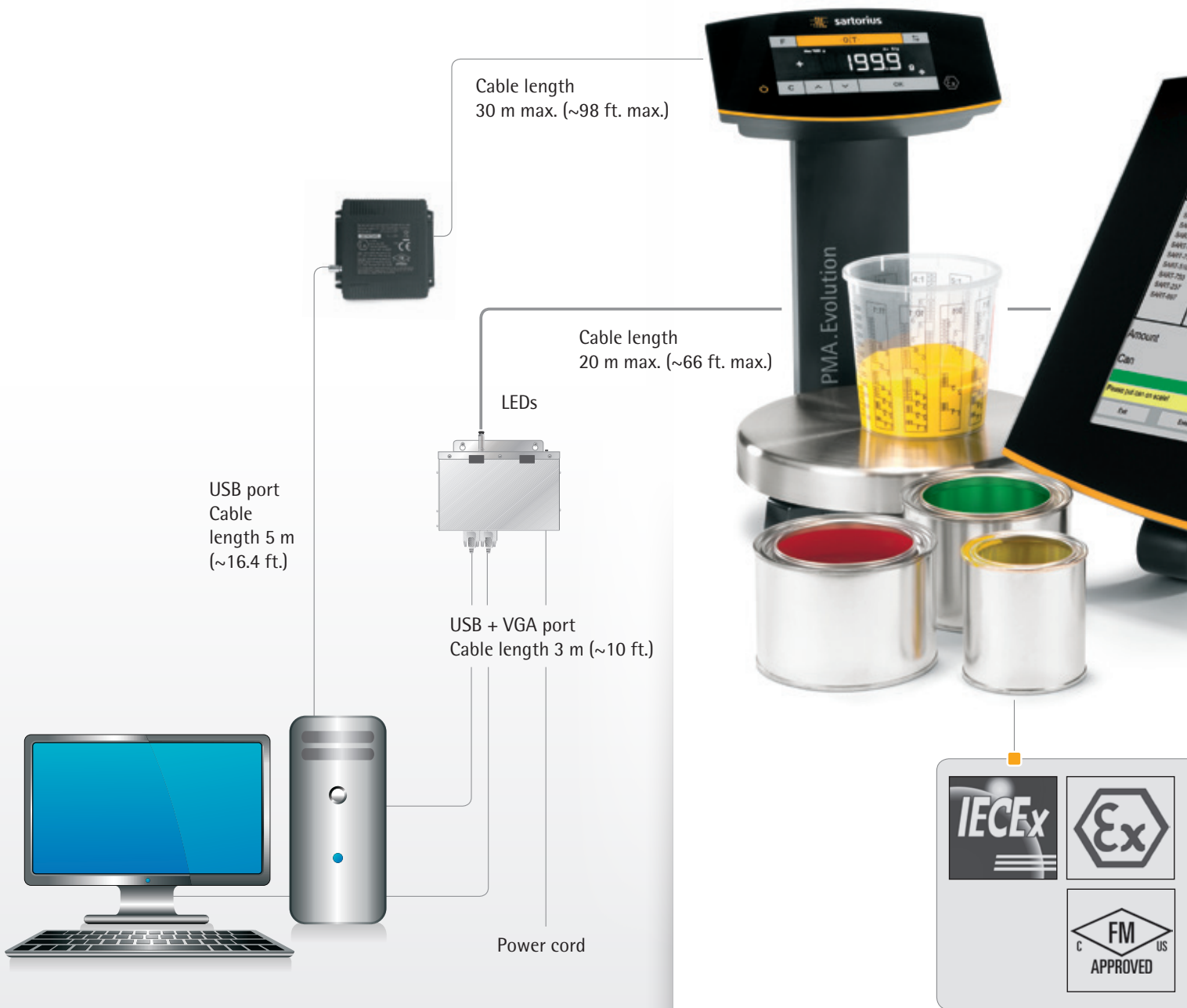
TopMix2.Touch The Perfect Touch for Your Color



turning science into solutions

TopMix2.Touch

No two reds are alike – finding the exact match requires solid craftsmanship as well as many years of experience and a feel for the right color mix. Just one look at the new Sartorius TopMix2.Touch and you'll see why its features define the new benchmarks in the paint mixing room: Its 15" multi-touch display along with a compact footprint offers a unique level of operating convenience in Zone 1 | Class I, Div. 1, hazardous areas.



Non-Ex area (safe area)



New Display Technology

The 15" color display with the Multi Touch feature can be used either with gloves or none.



Our Touch Pen with a holder will make it easy for you to input data. Just effortlessly attach it to the TopMix housing.



Easy to Clean

The new glass front display assures a comfy and fast cleaning to save time and money.



An in-use cover is also available to minimize your cleaning effort by preventing stubborn paint stains from building up.



Unique Connectivity

The new One Cable technology designed for use inside the Zone 1 | Class I, Div. 1 area enables more efficient installation, saving you valuable time.



reddot award 2014 winner

Safety Comes First

TopMix2.Touch has been tested and approved according to recognized international standards. This means your safety is ensured at all times in your paint-mixing area.

Sales and Service Contacts

For further contacts, visit www.sartorius.com

Europe

Germany
Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 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 RUS"
Uralskaya str. 4, Lit. B
199155 St. Petersburg
Phone +7.812.327.53.27
Fax +7.812.327.53.23

Spain & Portugal
Sartorius Spain, S.A.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)
Phone Spain +34.913.586.095
Phone Portugal +351.800.855.800
Fax Spain +34.913.589.623
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
LLS "Sartorius RUS"
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.
Libramiento Norte de Tepetzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepetzotlan,
Estado de México,
C.P. 54605
Phone +52.55.5562.1102
Fax +52.55.5562.2942
leadsmex@sartorius.com

Peru
Sartorius Peru S.A.C.
Av. Emilio Cavenecia 264 San Isidro
15073 Lima, Perú
Phone +51.1.441 0158
Fax +51.1.422 6100

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 (Shanghai) Trading Co., Ltd.
3rd Floor, North Wing, Tower 1
No. 4560 Jinke Road
Zhangjiang Hi-Tech Park
Pudong District
Shanghai 201210, P.R. China
Phone +86.21.6878.2300
Fax +86.21.6878.2882

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