



TopMix2.Touch El Toque Perfecto para Su Color

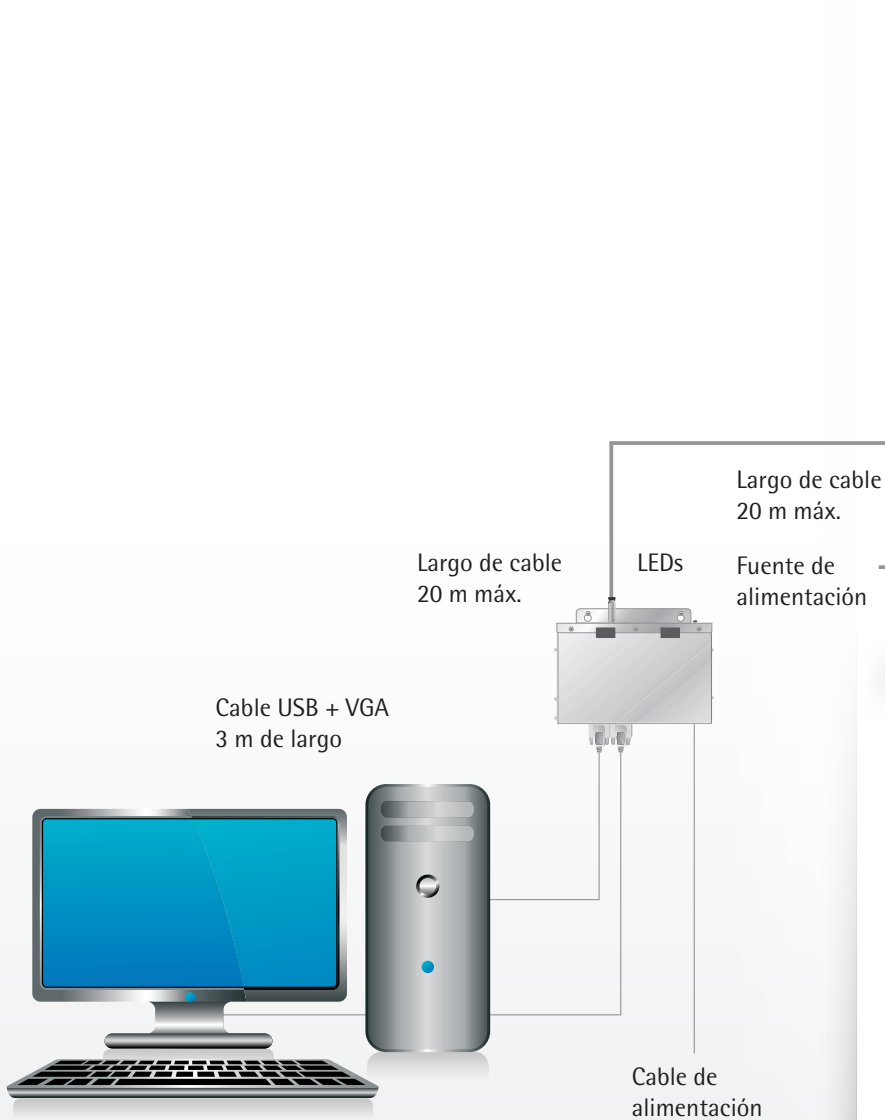
Job No.: XYZ12345 Formula No.: Sarto 9876
 Manufacturer: Variant: Darker
 Mfr. Code: AB-456 Color Chip: CC-1468
 Color Name: Brilliant Red2 Formula Date: 26.02.2006
 Mfr. Year: 1994 - 2002 Paint Line: Water Based

Component	Component Name	target	target cumulated	Actual	Actual cumulated	Deviation	tolerance
SART-456	Light Blue	212.6 g	212.6 g	0.0 g	0.0 g	-212.6 g	+/- 0.2 g
SART-852	White	136.2 g	348.8 g	0.0 g	0.0 g	-136.2 g	+/- 0.2 g
SART-516	Yellow	61.5 g	410.3 g	0.0 g	0.0 g	-61.5 g	+/- 0.1 g
SART-154	Dark Red	35.1 g	445.4 g	0.0 g	0.0 g	-35.1 g	+/- 0.1 g
SART-353	Medium Green	21.7 g	467.1 g	0.0 g	0.0 g	-21.7 g	+/- 0.1 g
SART-750	Black	19.0 g	486.1 g	0.0 g	0.0 g	-19.0 g	+/- 0.0 g
SART-510	Purple	11.8 g	497.9 g	0.0 g	0.0 g	-11.8 g	+/- 0.0 g
SART-753	Signal Red	9.1 g	507.0 g	0.0 g	0.0 g	-9.1 g	+/- 0.0 g
SART-237	Silver	5.7 g	512.7 g	0.0 g	0.0 g	-5.7 g	+/- 0.0 g
SART-697	Brown	3.3 g	516.0 g	0.0 g	0.0 g	-3.3 g	+/- 0.0 g

Highly flammable

TopMix2.Touch

No hay dos rojos parecidos - encontrar la mezcla exacta requiere un conocimiento sólido, al igual que muchos años de experiencia y tacto para la mezcla de color indicada. Solo vea la nueva Sartorius TopMix2.Touch y verá por que sus funciones definen nuevos puntos de referencia en la sala de mezclado de pintura: Su pantalla multi-touch de 15" y tamaño compacto ofrece un nivel único de conveniencia de operación en Zona 1 | Clase 1, Div. 1 áreas peligrosas.



Zona No Peligrosa (Área No Ex)



Nueva Tecnología de Display

La pantalla de 15" a color con función multi-touch puede ser usada con o sin guantes.



Fácil de Limpiar

La nueva pantalla de vidrio asegura una limpieza rápida y confortable para salvar tiempo y dinero.



Conectividad Única

La nueva tecnología de un solo cable dentro de la Class | Div Area lleva a una instalación más eficiente y rápida.



reddot award 2014
winner

Seguridad Primero

La TopMix2.Touch ha sido probada y aprobada de acuerdo a reconocidos estándares internacionales. Esto significa que su seguridad siempre está asegurada en todo momento en su área de mezclado de pintura.

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 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 +34.902.123.367
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

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/es