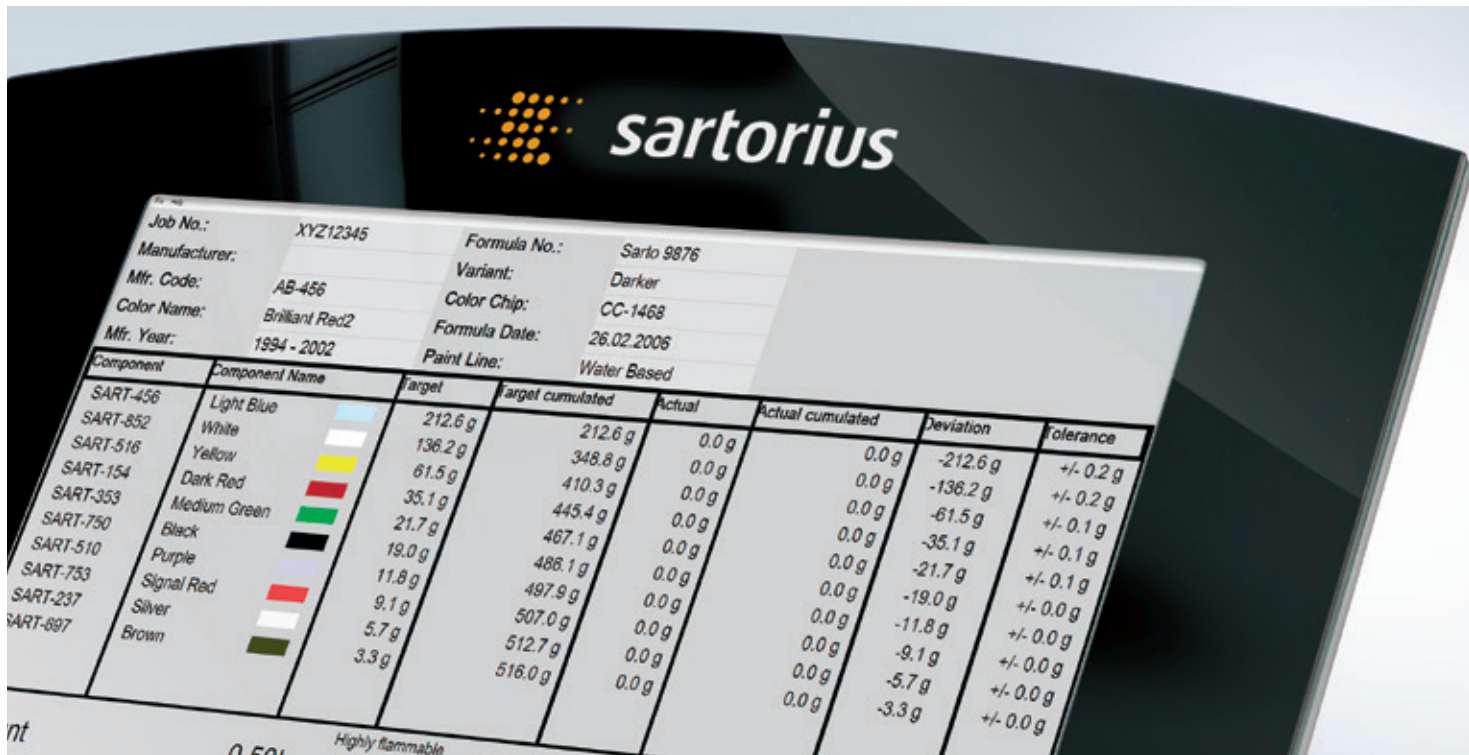


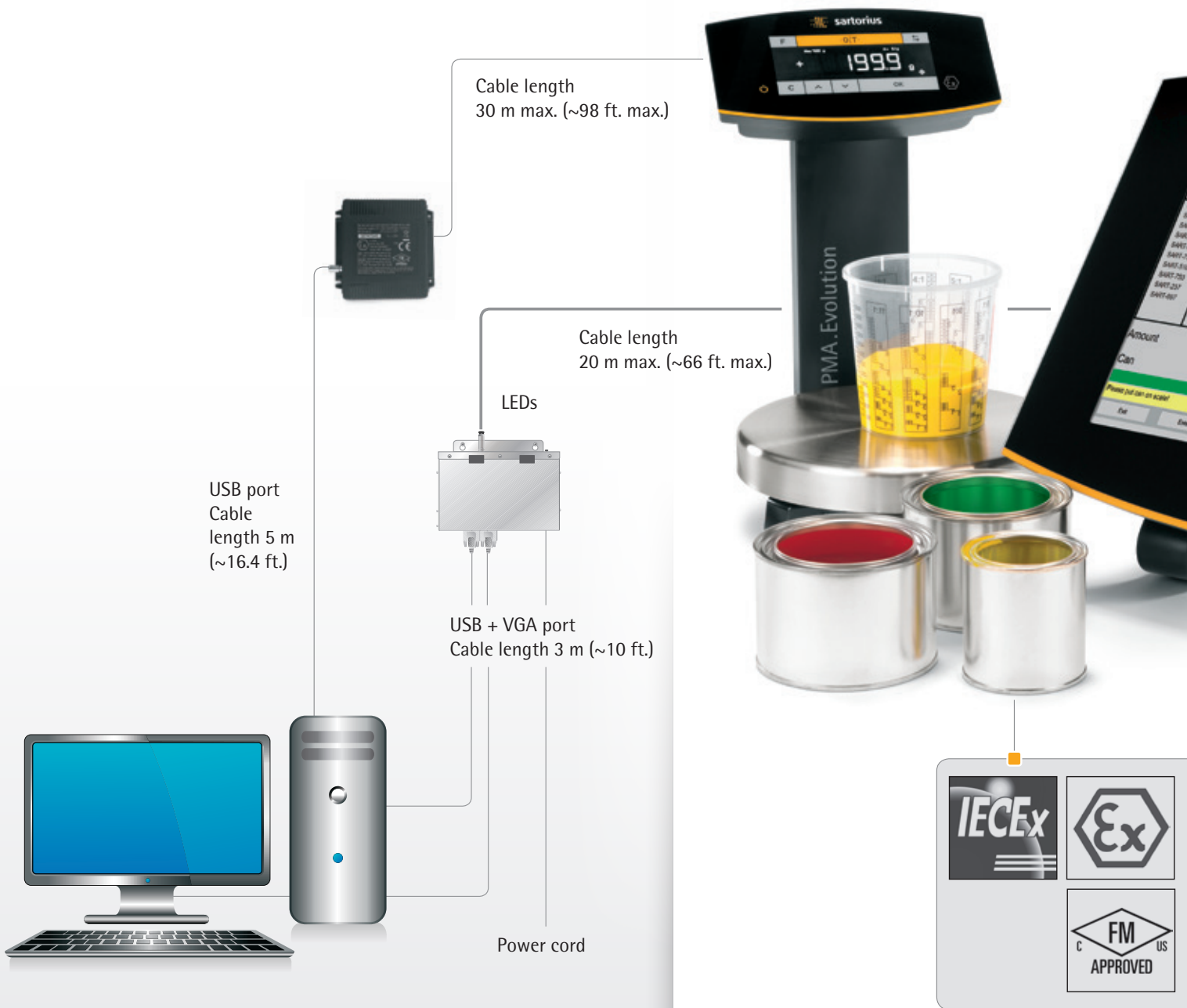


TopMix2.Touch The Perfect Touch for Your Color



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.

Ex area (hazardous area)

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

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



◀ www.sartorius.us