

# TopMix2.Touch Package (Model TMP020-X00002)

For use in Zone 1 hazardous areas.

Paint Mixing Terminals



- Up-to-date display technology: prepared for today's and tomorrow's paint mixing applications (size, resolution, multi-touch)
- Glass front display: easy to clean, save's time and money
- Unique – one cable technology inside the Class I Div 1 area: allows a more efficient time saving installation
- Compact design (weight & dimensions): easy handling, save on storage space
- VESA physical mounting interface (VESA MIS-E, 200 x 100 mm)

## Technical Specification

### Terminal TM02-X

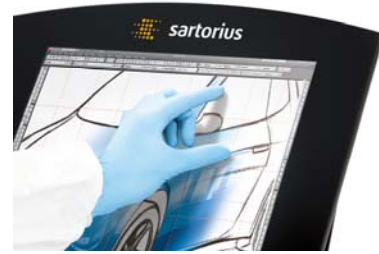
Display	15" XGA (1024 x 768)
Housing (WxDxH)	420 x 315 x 322 mm (incl. mount bracket)
Weight	approx. 8.5 kg (incl. mount bracket)
Data interface	RS232, integrated Power/VGA in
Allowable operating temperature	0 ... +40 °C   +32 ... 104 °F
Approval	Ex zone 1   Class I, Div 1
IECEX	Ex ib mb IIB T4 Gb (IECEX DEK 12.0072 X)
ATEX	II 2G Ex ib mb IIB T4 Gb (DEKRA 12ATEX0246 X)
FM (US)	IS CL I, DIV 1, GP C,D T4   CL I, Zone 1, AEx ib mb IIB T4
FM (CA)	Ex ia   CL I, DIV 1, GP C,D T4   Zone 1, Ex ib mb IIB T4
Electromagnetic Compatibility	In accordance with EN 61326-1:2006

### Ex-link Converter YC012-Z

Housing (WxDxH)	375 x 175 x 73 mm
Weight	approx. 1.7 kg
Power voltage	100 – 240 V   50 – 60 Hz   25 VA
Data interface	Integrated Power/VGA out, USB in (Typ B), DVI in, DVI out, Power in
Allowable operating temperature	0 ... +40 °C   +32 ... 104 °F
Approval	Ex zone 1   Class I, Div I
IECEX	[Ex ib Gb] IIB (IECEX DEK 12.0072 X)
ATEX	II (2)G [Ex ib Gb] IIB DEKRA 12ATEX0246 X)
FM (US)	AIS CL I, DIV 1, GP C,D   CL I, Zone 1, [AEx ib] IIB
FM (CA)	Ex ia   CL I, DIV 1, GP C,D   Zone 1, [Ex ib] IIB

### Standard equipment

TopMix2.Touch terminal/control panel for Zone 1 potentially explosive areas	TM02-X
Ex-link converter	YC012-Z
Ex-link cable (20 m)   (65.6 ft)	YCC01-0049M20
Video cable DVI to VGA (3 m)   (9.8 ft)	YCC01-0051M3
USB cable (3 m)   (9.8 ft)	YCC01-0040M3
D-SUB 9-pin cable (3 m)   (9.8 ft)	YCC01-0041M3
Bracket for wall and table installation	YWB06
3x Grounding cable (2 m)	YCC01-X046M2



### Optional equipment

Video cable DVI to DVI (3 m)   (9.8 ft)	YCC01-0050M3
---	--------------



YC012-Z, Ex-link converter

### Technical Drawing

