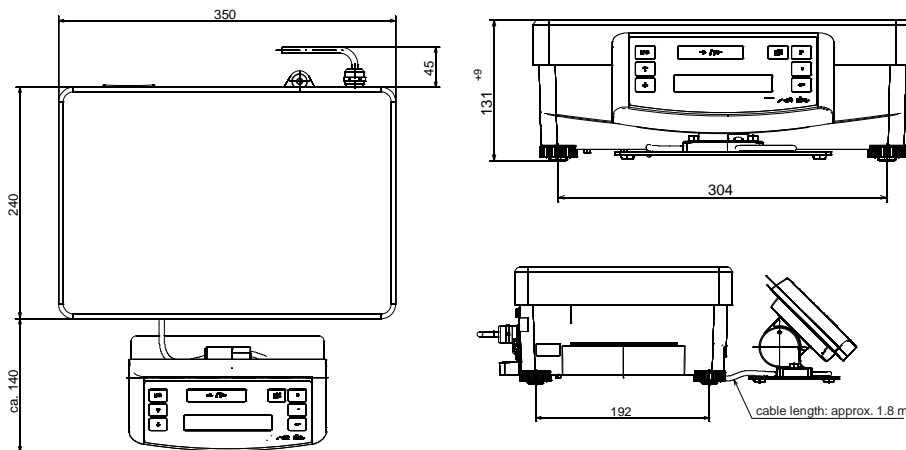


Sartorius PMA.Power for Ex zone 2 (Model PMA35001-Y)



Product features:

- Large 14-segment display can be arranged in any way on the table or at the wall (cable length 1.8m)
- Display lock function
- Communication symbol
- Scrolling text function
- Correction of formulation over-pours
- Calculation by factor function
- Stainless steel weighing pan, removable
- Housing resistant to solvents
- Protected against overload
- Standard serial interface port
- Internal power supply 100...240V
- Can be used in flammable environments



Standard equipment:

- Bracket for wall mounting
- Adapter 25 pin. to 9 pin.

Optional equipment:

- In-use cover for display
- Data cable to connect to a PC

Other models:

- Approved for Ex zone 1
- Standard non Ex approved

General technical data

Weighing capacity	35000 g
Readability	0.1 g
Manufacturer declaration of conformity according "Directive 94/9EC":	II 3 G Ex nA nL IIB T4 IP4x
Tara range (subtractiv)	-35000 g
Allowable ambient operating temperature	+10 up to + 30 °C
Load plate	350 x 240 mm
Scale housing (WxDxH)	350 x 240 x 132,5 mm
Net weight	Approx. 11 kg
Interface	RS232
Interface port	25-pin
Calculation factors	0.5 - litres