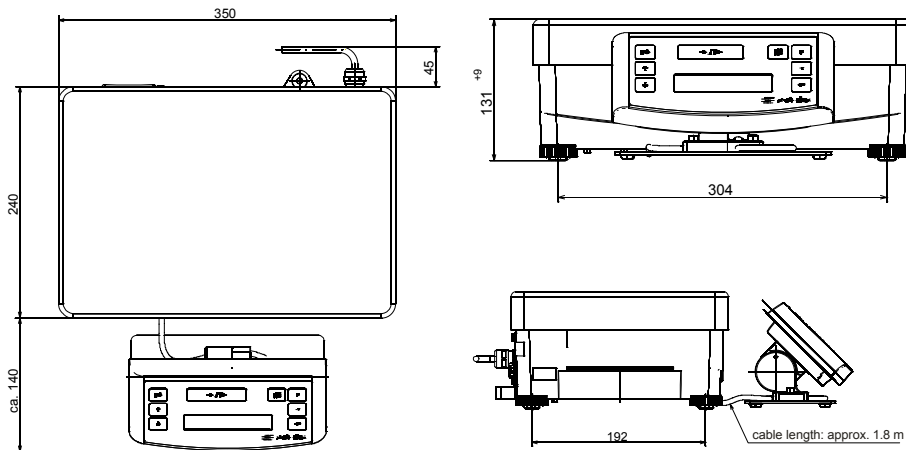




## Sartorius PMA.Power (Model PMA35001)



### 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 2

### General technical data

Weighing capacity	35000 g
Readability	0.1 g
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   1.0   1.5   2.0   2.5   3.0   3.5   4.0   4.5   5.0   10.0   15.0   20.0   25.0 litres