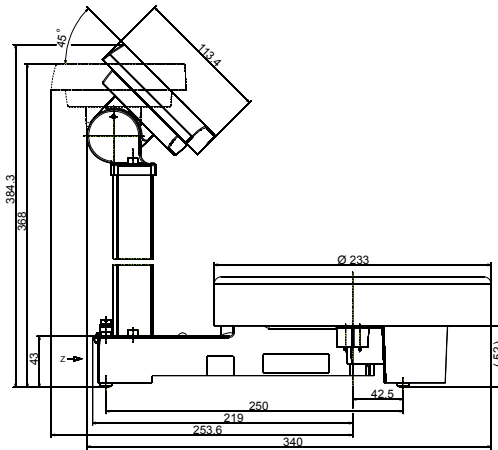
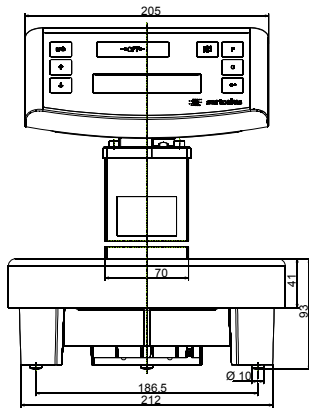




Sartorius PMA.Quality (Model PMA7501)



Product features:

- Large 14-segment display
- Display lock function
- Communication symbol
- Scrolling text function
- Correction of formulation over-pours
- Calculation by factor function
- Housing resistant to solvents
- Round stainless steel pan, wrap around design
- Protected against overload
- Raised display adjustable (removable)
- Standard serial interface port
- Power supply provided with rating for specific countries from 90-264V, 50-60 Hz.
- Can be used in flammable environments



Optional equipment:

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

Other models:

- Approved for Ex zone 1 | Ex zone 2

General technical data

Weighing capacity	999.95 g 7500 g
Readability	0.05 g 0.1 g
Tara range (subtractiv)	-999.95 g -7500 g
Allowable ambient operating temperature	0 up to + 40 °C
Pan size (∅)	233 mm
Scale housing (WxDxH)	233 x 340 x 385 mm
Net weight	Approx. 3.2 kg
Interface	RS232
Interface port	9-pin
Calculation factors	0.25 0.5 0.75 1.0 1.5 2.0 2.5 3.0 3.5 4.0 4.5 5.0 liter