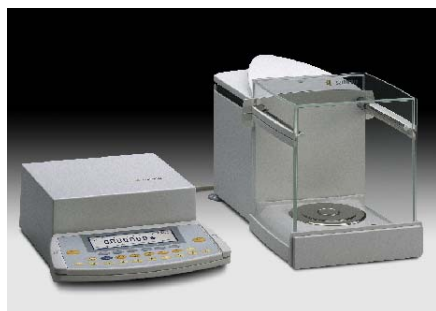




ME36S Microbalance



Premium Microbalance with outstanding weighing capacity.

- Outstanding weighing technology featuring exceptionally fast response times
- Fully automatic draft shield system for fast adaptation to any work routine
- Draft shield operation with learning function activated by user-friendly draft shield keys
- Ideal for use as measuring and test equipment in quality management systems
- SQmin function for displaying the minimum sample quantity according to the United States Pharmacopeia
- New S.U.R.E. function for dynamic display of the uncertainty of measurement according to the DKD Calibration Certificate
- Fully automatic calibration|adjustment function, isoCAL
- Backlit graphical display with plain-text prompts and info on all settings
- Highly resistant membrane keypad overlay with "bubble-shaped" keys
- Complete package of applications is standard
- RS-232C port plus additional Sartorius printer port

Model	ME36S
Readability	1 µg
Weighing capacity	31 g
Repeatability	± ≤ 2 µg
Max. linearity	± ≤ 10 µg
Avg. response time	14 - 18 s
Pan size	Ø 30 mm

Key accessories	Order No.
Data printer with date, time and statistics functions	YDP03-OCE
Foot switch for draft shield, tare and print functions	YPE01RC
Data transfer software SartoConnect with RS-232C standard cable (1.5 m);	YSC01L
Ionizing blower	YIB01-ODR
Balance table	YWT01