

Cubis[®] MPS6.6S

Pipette Calibration Balance

Benefits

- Motion Control, quick opening of the draft shield: convenient and efficient pipetting
- Integrated humidity sensor: controls humidity inside the weighing chamber
- White LED light: illumination of the full weighing area
- Red LED light: instant alert if climate parameters are out of tolerance

Product Information

For professional pipette calibration, you need the right balance — one that combines accuracy, precision, speed and convenience.

Accuracy and precision to comply with international regulations. Speed to meet requirements for cost- and time-efficient performance. And convenience for working in a comfortable posture. Does all this sound impossible? Not if you find the right balance. Meet the Sartorius Cubis[®] MPS: The perfectly balanced system designed to reconcile all your needs for accuracy, precision, speed and convenience in pipette calibration.



Technical Specifications


General		Specific	
Power supply:	100–240 V~, -15% +10%, 50–60 Hz, 1.0 A	Readability [mg]	0.001
Power consumption:	<35 VA	Weighing capacity [g]	6.1
Ambient temperature:	Operation +10 °C to +30 °C	Typical response time [s]	5
Highest relative humidity:	80% for temperatures up to 31 °C, decreasing linearly to 50% relative humidity for 40 °C	Repeatability [\pm mg]	0.001 Standard deviation
Safety of electrical equipment:	According to EN 61010-1:2010: Safety requirements for electrical equipment for measurement, control, and laboratory use – Part 1: General requirements	Linearity [\pm mg]	0.004
Electromagnetic compatibility:	According to EN 61326-1:2013: Electrical equipment for measurement, control, and laboratory use – EMC requirements – Part 1: General requirements	Eccentric load [mg] (Test load [g])	0.007 (5 g)
Defined immunity to interference:	Suitable for use in industrial areas	Number of possible pipette calibration at 10.3 ISO Standard	33 (10 μ L pipette)
Interference emission	Class B (suitable for use in residential areas and areas that are connected to a low voltage network that also supplies residential buildings). The device can therefore be used in both areas.		

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further information, visit
www.sartorius.com