

MPS116S

Cubis® II Single Channel Pipette Calibration System



Benefits

- Cover about 95 % of all single channel pipettes with only one instrument thanks to 111 g weighing capacity combined with 120 mL weighing vessel
- Get the best performance in terms of resolution and speed with adaptable readability between 1 and 100 μg
- Easy and ergonomic handling due to an automated lid avoiding evaporation combined with a large moisture trap
- Keep all relevant data monitored: Together with the Cubis® II climate module and the option to implement third party sensors for water temperature into the instruments water reservoir

Technical Specifications

Max. Capacity	g	111
Readability	mg	0.001 (option to turn off up to 3 digits)
Repeatability at 5 %	mg	0.006
Typical repeatability at 5 %*	mg	0.004
Pipetting weighing vessel volume	mL	120
Nominal volume pipette calibration range** (in accordance to ISO8655)	µL	2 - 5,000
Typical stabilization time at 0.001 mg readability	s	3
Typical stabilization time at 0.01 mg readability	s	1.5
Typical stabilization time at 0.1 mg readability	s	1.2
Norms supported		ISO 8655
Dimensions (L x W x H)	mm	510 x 240 x 301

*This specification is within ISO8655 limit values for pipettes below 20 µL.

**Higher volume range is possible by emptying the vessel in between.

General Specifications

Power supply	V Hz A	100 - 240 ± 10 % 50 - 60 0.8
Power consumption (max.)	W	64.5
Ambient temperature	°C	+ 5 to + 40
Relative humidity	%	80% for temperatures up to 31 °C, decreasing linearly to 50% relative humidity for 40 °C
IP level		IP30

Accessories

	Order No.
Suction pump for emptying the weighing vessel	YCP05-SP01
Climate module, uncalibrated	YCM20MC
Climate module with DAkkS certificate	YCM20MC-DAKKS
Climate module with tower, uncalibrated	YCM20MC-Tower
DAkkS certificate for climate module	YCM20DAKKS
Wheeled transportation case	YDB02MPS

Weighing tables


	Order No.
Made from synthetic stone with vibration dampening	YWT03
Made from wood with synthetic stone	YWT09
Wall console	YWT04

Germany

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

USA

Sartorius Corporation
3874 Research Park Dr.
Ann Arbor, MI 48108
Telephone +1 734 769 16006

 For further contacts, visit
www.sartorius.com