



Revolutionize Pipette
Calibration —
From 2 μ L to 5 mL with
the Cubis® II MPS116S

Simplifying Progress

SARTORIUS

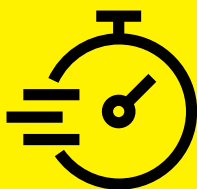
Cubis® II MPS116S: Redefining Pipette Calibration— Efficient, Mobile, and Space-Saving

In modern calibration labs, every centimeter of bench space, every minute saved, and every error avoided matters. The **Cubis® II MPS116S** pipette calibration solution meets these demands with an all-in-one approach that replaces multiple instruments, streamlines workflows, and simplifies handling—whether you're in-house or on the move. Capable of calibrating pipettes from **2 µL up to 5 mL**, it covers nearly all pipette volumes in a single device



Optimizes Use of Lab Space

By integrating three balances into a single device, the MPS116S frees up valuable bench space — ideal for labs where every inch counts.



Faster Throughput, Less Downtime

Minimal setup time, faster stabilization time and fewer balance changes mean technicians can calibrate more pipettes per day. And if another balance fails, the Cubis® II MPS116S seamlessly steps in — ensuring no workflow interruptions.



One Device. Total Mobility

Optimized for field use, the Cubis® II MPS116S requires only a single instrument to cover a broad volume range (2 µL to 5 mL). It adapts quickly to new environments and is ready for accurate measurements within one hour of setup — saving time and simplifying logistics for mobile service teams.



Ergonomics That Work for You

Designed for long calibration routines, the Cubis® II MPS116S features intuitive, ergonomic operation. The automatic lid reduces evaporation, while the integrated moisture trap ensures stable results — pipette after pipette.



Cost Savings

With only one balance to maintain and calibrate instead of two or more, labs benefit from fewer service contracts, less downtime, and lower long-term costs.



Environmental Monitoring Built In

The Cubis® II MPS116S connects directly with the Cubis® II climate module and supports integration of third-party sensors into the water reservoir. This gives you full control over critical environmental parameters like water temperature — fully traceable and documented.

Testimonial

"The combination of micro, semi-micro and analytical balances makes the device **extremely versatile**.

The **very short stabilization time** enables **efficient work**.

The clever geometry of the weighing vessel **minimizes evaporation and thus enables precise work** down to the lower volume range."

— Jan Klinkner, GM, Klinkner & Partner



Dr. Roman Klinkner and successor Jan Klinkner

About Klinkner & Partner:

Klinkner & Partner GmbH operates a DAkkS-accredited calibration laboratory for pipettes, dispensers, and other liquid handling instruments. Established in 2014 in Lahr (Black Forest), Germany, the laboratory provides maintenance, repair, and calibration for all piston-operated devices on the market – regardless of manufacturer or model – to customers in Germany and abroad.

Accredited according to DIN EN ISO/IEC 17025 and DIN EN ISO 8655, the laboratory employs 15 staff members and is known for delivering reliable, precise, and high-quality services.

With a comprehensive full-service approach – including accredited calibration, routine maintenance, equipment repairs, rental services, sales consulting, and user training – Klinkner & Partner meets the highest standards of quality and service "Made in Germany."

In addition to its laboratory services, this family-owned company has been advising and training laboratories across all industries since 1994. Through seminars, courses, and expert forums, Klinkner & Partner supports customers in the areas of laboratory management, quality assurance, and analytics.





Upgrade Your Existing Cubis® II Balance

Already working with a Cubis® II balance? The YCP01MPS accessory gives your existing equipment a powerful new role – transforming it into a fully functional pipette calibration system without the need for a separate device. It's the ideal way to protect your investment and expand your lab's capabilities at minimal cost.

Universal Compatibility

The YCP01MPS is designed to work seamlessly with all Cubis® II balances of the 3rd generation, including the following models:

- MCA36P | MCA66S | MCA116S
- MCA226S | MCA125S | MCA225S|P

Recommended Combinations for Best Performance

To ensure optimal results and user experience, we recommend pairing:

- MCA36P + YCP01MPS – Ideal for calibrating low-volume pipettes (starting from 2 µL) with high precision.
- MCA125S + YCP01MPS – Perfect for covering a wide pipette volume range up to 5 mL.

Maximize Efficiency with Minimal Effort

Installation is quick and intuitive. Once connected, the accessory unlocks all dedicated pipette calibration functions directly within your Cubis® II balance—giving you access to robust software features and automated processes tailored to ISO 8655 calibration workflows.

Whether you're expanding your service offerings or standardizing your internal lab equipment, the YCP01MPS gives you the flexibility to do more—without additional hardware investments.

Pipette Volume	2 µL	10 µL	20 µL	100 µL	200 µL	1000 µL	5000 µL
ISO 8655 Requirements	0.001 mg			0.01 mg			0.1 mg
MPS116S	2 µL to 5,000 µL						
MCA36P* S + YCP01MPS	2 µL to 200 µL						
MCA66S P + YCP01MPS	2 µL to 200 µL						
MCA116S 226S + YCP01MPS			20 µL to 200 µL				
MCA125S* 225S P + YCOP01MPS			20 µL to 5,000 µL				

* recommended Cubis® II balances to use with YCP01MPS



Technical Specifications

Maximum Capacity	g	111
Readability	mg	0.001 (option to turn off up to three 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 ISO 8655)	µL	2 – 5,000
Typical Stabilization Time at 0.001 mg Repeatability	s	3
Typical Stabilization Time at 0.01 mg Repeatability	s	1.5
Typical Stabilization Time at 0.1 mg Repeatability	s	1.2
Norms Supported		ISO 8655
Dimensions (L x W x H)	mm	510 x 240 x 301

* This specification is within ISO 8655 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 (maximum)	W	64.5
Ambient Temperature	°C	+5 to +40
Relative Humidity	%	80% for temperatures up to 31 °C, decreasing linearity to 50% relative humidity for 40 °C
IP Level		IP30

Accessories

Order Number

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	YCM20MCDAKKS
Wheeled Transportation Case	YDB02MPS

Weighing Tables

Order Number

Made From Synthetic Stone With Vibration Dampening	YCP505-SP01
Made From Wood With Synthetic Stone	YCM20MC
Wall Console	YWT04



Germany

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

USA

Sartorius Corporation
3874 Research Park Drive
Ann Arbor, MI 48108
Phone +1 734 769 1600