



Pipette Calibration
Solutions

Simplifying Progress

SARTORIUS

Taking care of your pipettes is important to ensure the quality of work performed in the laboratory. In addition to cleaning and performing routine maintenance, pipette accuracy and precision should be checked at regular intervals. Smart solutions for pipette calibration can help laboratories achieve:



Compliance according to ISO 8655:2022



Fewer Errors



Increased Efficiency



Lower Calibration Costs

Overview



Product	MPS6.6S MPS105S	Speedcal Mobile	YCP04MS VF988 YCP07MC
Description	Single Channel Pipette Calibration System	Multichannel Pipette Calibration System with 4,8 or 12 channels	Moisture Trap for Cubis® II Balances
Maximum Capacity	6.1 100 g	Up to 12 x 21 g	6.1 to 520 g*
Readability	0.001 0.01 mg	0.01 mg	0.001 0.01 0.1 mg*
Volume of pipetting vessel	6 mL 100 mL	21 mL per channel	21 mL
Pipette calibration range (acc. to ISO 8655:2022)**	From 0.5 µL to 5,000 µL	From 0.5 µL to 1,000 µL	From 0.5 µL to 1,000 µL
Ingenix compatibility	Yes	Yes	Yes
Automatically opening lid	Yes	No	No

*Depending on the Cubis® II balance being used

**without emptying the vessel during one 3 x10 calibration cycle

Cubis® MPS

The Solution for Calibration of Single Channel Pipettes

Cubis® MPS is a portfolio of specialized systems designed to perform fast, ergonomic and precise pipette calibrations. These highly specialized instruments are for customers who use their balances exclusively for pipette calibrations and need fast stabilization times and optimal ergonomic handling.

Automatic Operation

Cubis® MPS comes with a touch-free automatic draft shield supporting error-free and ergonomic workflows. This motion sensor technology recognizes the user approaching the moisture trap from a wide-angle range and ensures proper closing after the dispense. It closes again only if the user removes the pipette.

A Complete Calibration Lab in One

Cubis® MPS is a standalone solution for climate data management and water temperature control. It can be equipped with a climate tower and temperature-controlled liquid reservoir. Workflows are integrated to guide users through the pipette calibration routine (acc. to ISO 8655:2022), calculate weight values into volumes (incorporating environmental data as temperature and barometric pressure) and evaluate according to ISO 8655:2022 tolerances.

Two models of Cubis® MPS are offered:

- MPS6.6S for higher resolution and calibration of pipettes down to 0.5 μL
- MPS105S for higher maximum capacity and calibration of pipettes from 20 μL to 5 mL



Speedcal Mobile

The Solution for Calibration of Multi Channel Pipettes

Speedcal Mobile is a patented, unique multi-channel pipette calibration system which allows for parallel measurement of up to 12 volumes. Speedcal Mobile can be also used to calibrate single channel pipettes. With its integrated semimicro balances (readability of 0.01mg), it is possible to calibrate pipettes from 0,5 μ l according to ISO 8655:2022.

Speedcal Mobile benefits:

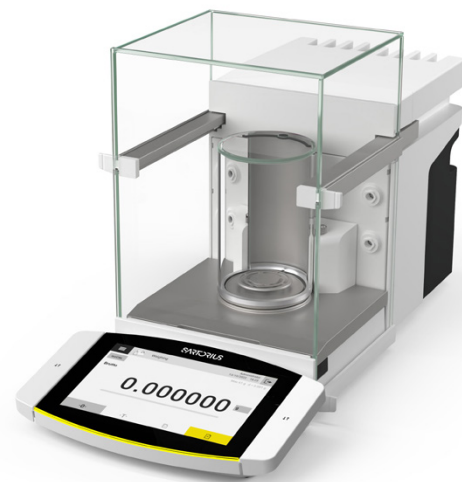
- Fast calibration (3 x 10) of a multichannel pipette thanks to up to 12 weighing cells in one instrument
- Parallel measurement ensuring ISO 8655:2022 compliance
- Speedcal Mobile is available in 4, 8 or 12-channel versions (later upgrade possible).
- High degree of mobility for on-site calibrations due to ergonomic, wheeled transport trolley



Moisture Traps

for Universal Cubis® II Balances

Your Sartorius Cubis® II Balance can easily be adapted for checking and calibrating your pipette by equipping it with an evaporation trap. Together with the Pipette Check QApp (QAPP005), the user is guided through the complete process for checking pipettes according to ISO 8655:2022 tolerances.



Ingenix Suite Advanced Pipette Calibration Module

ISO Compliant Software for Calibration and Pipette Management

The Advanced Pipette Calibration Module is an integral part of the Ingenix Suite software family when it comes to simplifying workflow for your pipette calibration.

The convenient interface offers a guided workflow for your pipette calibration tasks. It allows easy scheduling of calibration intervals and management of the respective pipette fleet.

As every other module, it can be installed and run as a stand-alone software solution for your pipette calibration tasks, or in combination within the Ingenix Suite family of modules.

It is completely compliant with ISO 8655:2022, ISO 17025 and 21 CFR Pt. 11. Network capabilities with automatic transfer to a SQL database provide central data management and the highest level of security standards.

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 contacts, visit

www.sartorius.com