

Sartorius Guides Your Every Drop. Calibrate Your Pipettes with Confidence

Whether you're seeking a straightforward routine check or a thorough compliance control service, our adaptable approach ensures that your pipette metrology requirements are fulfilled with precision and care. At Sartorius, we offer a comprehensive range of solutions designed to meet your specific needs in piston-operated volumetric apparatus (POVA) verification, ensuring that the ISO 8655 standard is fully adhered to in every service we provide.

Explore our guide and discover the perfect solution tailored to your pipette metrology requirements. Our guide is designed to help you navigate through the various options we provide, enabling you to select the service that best fits your needs, being:

- Prevention of Experimental Deviations & Instant Pipette Verification
- Compliant Internalization of Pipette Calibration
- Robust Set-up for On-Site Mobility
- Compliant Externalization of Pipette Calibration
- Certified Training to Learn Mastery of Pipetting Practices

With Sartorius, your pipettes will be verified and maintained to the highest standards, ensuring accuracy and reliability in all your critical laboratory processes. Our commitment to quality and customer satisfaction means that we are here to support you every step of the way, from routine checks to full compliance services.



Contact Sartorius to Discuss Your Pipette Calibration Needs,
visit **ISO 8655 | Sartorius**



SARTORIUS



Prevention of Experimental Deviations. Instant Pipette Verification

Application

- Rapid pipette performance check through critical pipetting series or before dispensing precious samples (pass/fail)
- Pipette accuracy assurance after autoclaving or regular cleaning (as found/as left)
- Detection of deviations requiring maintenance, cleaning, or metrology check-up
- Confirmation of pipette & tips compatibility - compliance
- Ergonomic setup with minimal operation time
- Guided operating protocol to limit human errors
- Training on good pipetting techniques for new incomers and beginners

Instruments

- Semi-micro balance with precision 0.01 mg
- Lab Water System

Software & Accessories

- Simple calculation of errors in % and μL , accommodating environmental conditions (temperature, humidity, air pressure)
- Evaporation trap

Service & Training

- Installation
- Best pipetting practice



Shopping
Cart

Instruments

Cubis® II MCA-225P-3S00-DMDS for nominal volume $\geq 20 \mu\text{L}$

Arium® Advance EDI Benchtop Type 2 H2O-EDI-1-T

Arium® Bagtank H2O-AOV-50

Software & Accessories

Climate module YCM20MC

Evaporation Trap YCP07MC recommended for tested volumes $< 50 \mu\text{L}$

Ingenix software for pipette checking DSX5PC01LH-FAC01 (or)
Preinstalled application in the balance QAPP005

Service & Training

Balance Installation

Water System Installation

Basic Training on Best Pipetting Practices



Contact Sartorius
to Discuss Your
Pipette Calibration
Needs



Compliant Internalization of Pipette Calibration

Application

- Compliance to ISO 8655 standard for Piston-operated volumetric apparatus (POVA)
- Detection of deviations requiring maintenance, cleaning, or metrology check-up (as found)
- Confirmation of pipette & tips compatibility - compliance
- Ergonomic setup with minimal operation time
- Daily working sheet, automatically listing pipettes to be checked
- Flexible pipette fleet management with preloaded pipette types and ISO 8655 requirements
- Guided operating protocol to limit human errors
- Report: conformity assessment, POVA out-of-tolerance, full traceability, uncertainty calculations
- Calibration certificates with ISO IEC 17025 accreditation* (optional)
- Certified training on good pipetting techniques & Standards

Instruments

- Semi-micro balance with precision 0.01 mg
- Microbalance with precision 1 µg
- System to check several channels in parallel for multi-channel pipettes
- Lab Water System

Software & Accessories

- Climate module
- Software for assistance in complying with ISO 8655, ISO 17025 & 21CFR part 11
- Unambiguous reports for increased audit readiness

Service & Training

- Installation
- Calibration
- Training

Multi Channel Pipettes (MCP)

Speedcal Model 4, 8 or 12 channels -
Capacity 21 g - 0.01 mg / channel
Ingenix Pipette Advanced Calibration Software
DSX5PC01LWGADC01
Balance Installation & Calibration
Arium® Advance EDI Benchtop Type 2 H2O-EDI-1-T
Arium® Bagtank H2O-AOV-50
Water System Installation
Training on Pipetting Standards & Maintenance

Single Channel Pipettes (SCP)

Cubis® MPS105S for calibration of pipettes from 20 µL to 5 mL
(or)
Cubis® MPS6.6S for calibration of pipettes down to 0.5 µL
Ingenix Pipette Advanced Calibration Software
DSX5PC01LWGADC01
Balance Installation & Calibration
Arium® Advance EDI Benchtop Type 2 H2O-EDI-1-T
Arium® Bagtank H2O-AOV-50
Water System Installation
Training on Pipetting Standards & Maintenance



Shopping
Cart



Contact Sartorius
to Discuss Your
Pipette Calibration
Needs

Robust Set-up for On-Site Mobility



Application

- Compliance to ISO 8655 standard for Piston-operated volumetric apparatus (POVA)
- Ergonomic setup with minimal operation time
- Daily working sheet, automatically listing pipettes to be checked
- Flexible pipette fleet management with preloaded pipette types and ISO 8655 requirements
- Guided operating protocol to limit human errors
- Certified training on good pipetting techniques & standards
- Service report: conformity assessment, POVA out-of-tolerance, full traceability, uncertainty calculations
- Transport trolley or case to secure the weighing systems
- Calibration weights in cases



Instruments

- Semi-micro balance with precision 0.01 mg
- Microbalance with precision 1 µg
- System to check several channels in parallel for multi-channel pipettes

Software & Accessories

- Climate module
- Software for assistance in complying with ISO 8655, ISO 17025 & 21CFR part 11
- Unambiguous reports for increased audit readiness

Service & Training

- Training



Shopping
Cart

Multi Channel Pipettes (MCP)

Speedcal Model 4, 8 or 12 channels -
Capacity 21 g – 0.01 mg / channel

Ingenix Pipette Advanced Calibration Software
DSX5PC01LWGADC01

Calibration weight set – several options

Single Channel Pipettes (SCP)

Cubis® II MCA225P-3S00-D MDS for nominal volume $\geq 20 \mu\text{L}$
(or)

Cubis® II MCA66P-3S00-D MDS for nominal volume $<20 \mu\text{L}$

Evaporation Trap YCP07MC

Climate module YCM20MC

Ingenix Pipette Advanced Calibration Software

DSX5PC01LWGADC01

Transport box for Cubis® II Balances YDB01MC-3

Calibration weight set – several options



Contact Sartorius
to Discuss Your
Pipette Calibration
Needs



Compliant Externalization of Pipette Calibration

Application

- Delegation of calibration, maintenance and/or repair of Piston-operated volumetric apparatus (POVA)
- Compliance to ISO 8655 standard for Piston-operated volumetric apparatus (POVA) and calibration certificates with ISO IEC 17025 accreditation*
- Flexibility to address internal customer requirements
- Service report compliant to ISO 8655-6:2022: conformity assessment, POVA out-of-tolerance, full traceability, uncertainty calculations

Sartorius Service Solution

- Onsite service* or at Sartorius metrology labs
- Full traceability of calibrations with undisputable POVA identification and unambiguous reports for an increased audit readiness
- Original spare parts from main other manufacturers
- Functional test, careful cleaning, lubrication, replacement of defective spare parts, check for damage and leak tightness, adjustment, calibration according to ISO 8655
- Verification of any volume from 0.1 μ L to several mL
- Measurement uncertainty calculation (on request)

*Not applicable to all countries. Please contact us.



Questions to Clarify

Fleet size
Number of single channel pipettes (SCP)
Number of multi channel pipettes (MCP)
Number and type of other POVA
Number of tip references per POVA
As found + as left certificates: Y/N
Maintenance: Y/N
Volumes to be checked

Number of repetitions / volume
On site or at Sartorius labs
On site: characteristics of the calibration lab
ISO 17025 compliance: Y/N
Frequency
Decontamination
ISO 8655 tolerances: Y/N
Max. pipette immobilization time
Budget



Contact Sartorius
to Discuss Your
Pipette Calibration
Needs



Certified Training to Learn Mastery of Pipetting Practices

Application

- Certified training on good pipetting techniques for new incomers and beginners
- Update and alignment on pipetting practices among experienced lab professionals
- Application-specific best pipetting handling to limit analytical variances
- Ergonomics improvement for long series of pipetting
- Avoid typical mistakes made when calibrating a weighing instrument

Trainings

- ISO 9001 certified Pipetting Academy
- Comprehensive combination of theory and hands-on practices
- Several modules developed for in-person or Virtual training
- Preliminary interview with your training manager to adapt the content and align the message with your lab challenges
- Training certificate sent to each participant after successfully completing the course
- Certified trainers



Shopping
Cart



Contact Sartorius
to Discuss Your
Pipette Calibration
Needs

Pipetting Academy

8620B6 to be selected or combined

Basics of pipetting
Ergonomics
Pipetting techniques
Pipetting in cell culture applications
Pipetting standards & maintenance
Best practices for working with proteins
Best pipetting practices for nucleic acid work
Best practices for cell-based sample preparation
Working with electronic pipettes

Lab Weighing Academy

8620C6 to be selected or combined

General Weighing Facts
Balance Calibration and Qualification
Quality Management and
Regulations

Germany

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



For further contacts, visit
www.sartorius.com

USA

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