

Calibration Certificate with ISO | IEC 17025 Accreditation

Pipettes and Dispenser



Benefits

- Gravimetric calibration according to ISO 8655
- Reassurance during quality audits
- Complete traceability to national standards
- International acceptance and recognition
- Calibration of pipettes and dispensers of all manufacturers
- Short calibration times

Product Information

The demands on the accuracy and reliability of laboratory data are constantly increasing. Pipettes must be calibrated regularly in order to achieve precise results and meet regulatory requirements. Our Sartorius Service has the right calibration for your needs. Calibration certificates with ISO | IEC 17025 accreditation are internationally recognized and provide the necessary security for your testing materials monitoring. The calibration ensures traceability to national standards and thus confidence in the accuracy of the measurement results. The uncertainty of measurement is also displayed in detail.

Specification

- Gravimetric calibration according to ISO 8655 (calibration of piston-operated volumetric instruments)
- Compliance with environmental conditions specified in ISO | IEC 17025
- 10 repeatability measurements per channel, each with three different volumes
- Calculation of the uncertainty of measurement
- Guaranteed traceability certificate to the national standard
- Affixation of a calibration mark
- Wide international recognition and acceptance of the calibration certificate
- Easy-to-read display of all measured values in the calibration

Customer Prerequisites

- Decontamination Declaration for devices
- For on-site service, provision of a suitable room (laboratory) for the installation of our calibration laboratories

Optional Services

- As-found calibration for documentation of the calibration results before maintenance work
- ISO 8655 calibration certificate for pipettes, dispenser, dilutors and burettes

You are interested?


Get in touch with us right away. We will be happy to provide you with a non-binding service offer. You can also find further information on our website: www.sartorius.com/en/services

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Straße 20
37079 Göttingen
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