

ISO 8655 Calibration Certificate (Manufacturer Certificates)

Pipettes, Dispenser, Dilutors
and Burettes



Benefits

- Gravimetric calibration according to ISO 8655
- Ensured, reliable calibration results
- Competent service for pipettes, dispenser, dilutors and burettes from other manufacturers
- Easy-to-read certificates
- Short calibration times

Product Information

It is imperative in ISO quality management systems that a test equipment management system be in place, and therefore that devices are calibrated regularly. Our Sartorius Service has the right calibration for your needs. With an ISO calibration certificate (manufacturer certificates) from Sartorius, your pipettes, dispenser, dilutors and burettes are calibrated according to DIN EN ISO 8655. The accuracy and quality of the measurement results are documented in an easily readable certificate.

Specification

- Gravimetric calibration according to DIN EN ISO 8655 (calibration of piston-operated volumetric instruments)
- Proof of compliance with the required accuracy tolerances
- Affixation of a test label | calibration sticker
- Easy-to-read display of all measured values in the calibration
- High acceptance of the calibration certificate

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
- Calibration certificate with ISO | IEC 17025 accreditation for pipettes and dispenser

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