

On-Site Maintenance and Calibration

Electronical and mechanical
Pipettes, Dispenser, Dilutors
and Burettes



Benefits

- Consistently reliable pipetting results
- Low logistic effort, short service times
- Optimized device performance
- Extended device life
- Regulatory compliance
- Available also for devices from other manufacturers

Product Information

Precision devices require regular professional maintenance and calibration to deliver consistently accurate and reliable results and fulfill all regulatory requirements. We recommend that you take out a maintenance contract so that you can fully rely on your devices and ensure maximum productivity.

Specification

- Functional test
- Careful cleaning
- Lubrication of seals, O-rings and pistons
- Replacement of defective seals and O-rings
- Check for damage and leaktightness
- Adjustment
- Calibration according to ISO 8655

Repairs, original spare parts, travel costs and other expenses are not included in the scope of services and will be charged separately.

Customer Prerequisites

- Decontamination declaration for devices
- Provision of a suitable room (laboratory) for the installation of our calibration laboratory
- Sartorius checks the location before the calibration to ensure that the required ambient conditions are met
- Pipettes are delivered to the calibration place and are sufficiently acclimatized

Optional Services

- As-found (initial) calibration for documentation of the calibration results before maintenance work
- Calibration certificate with ISO | IEC 17025 accreditation for pipettes and dispenser
- ISO 8655 calibration certificate for pipettes, dispenser, burettes and dilutors

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