

Preventive Maintenance and Service Contracts

Laboratory Balance



Benefits

- Consistent quality of measurement results
- Trouble-free operation, reduction of downtime
- Optimized device performance
- Extended device life
- Regulatory compliance

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 laboratory balance and ensure maximum productivity.

Specification

- Visual inspection for any equipment damage
- Workplace-related functional inspection of system components
- Filter settings and device leveling check
- Clean outside of device if possible
- Interior cleaning when function is impaired
- Measurement technology-related equipment inspection
- Internal calibration test (if applicable)
- Device adjustment (adjustment value, linearity, corner load) if necessary

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

Customer Prerequisites

- Device is at the maintenance site and has been sufficiently acclimatized
- The device is freely accessible

Optional Services

- Calibration certificate with ISO | IEC 17025 accreditation
- USP <41> certificate
- USP <1251> test report
- Balance test report
- Maintenance report

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