

Preventive Maintenance and Service Contracts

Manual Mass Comparator



Benefits

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

Product Information

Mass Comparators 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 manual mass comparator 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
- 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

- Comparator Test Protocol

Are you interested?

We would be pleased to hear from you! We are happy to provide you with a non-binding service quote. For more information about our service portfolio, please visit:
www.sartorius.com/en/services/instrument-service

Learn More:



Germany

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

USA

Sartorius Corporation
3874 Research Park Dr.
Ann Arbor, MI 48108
Phone +1 734 769 1600

 For further contacts, visit
www.sartorius.com