

Installation

Manual Mass Comparator



Benefits

- Perfectly functioning devices
- Reliable, reproducible results
- Professionally trained operating
- Extended device life

Product Information

The reliability and service life of your manual mass comparator depends crucially on the correct installation and configuration. Our Sartorius Service installs and configures your devices properly, taking into account the regulatory requirements and quality standards. Your manual mass comparator is immediately ready for operation and gives reliable and reproducible results right from the start.

Specification

- Checking the equipment delivery for completeness and any transport damage
- Professional installation of the equipment components
- Setup of default parameters
- Adaptation to ambient conditions at the place of use
- Initial startup with functional tests
- Initial metrological equipment check
- Instruction of operating personnel (no application training)
- Consultation on device maintenance and maintenance cycles

Customer Prerequisites

- Device is at the installation site and has been sufficiently acclimatized
- The installation place 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