

Installation

Laboratory Centrifuge



Benefits

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

Product Information

The reliability and service life of your laboratory centrifuges 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 laboratory centrifuge 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
- Initial startup with functional tests
- Device test run
- 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

- Installation report
- Calibration and test certificates
- GLP | GMP compliant equipment qualification

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