

# Preventive Maintenance and Service Contracts

## Laboratory Centrifuge



### 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 centrifuge and ensure maximum productivity.

## Specification

- Visual inspection for any equipment damage
- Workplace-related functional inspection of the system components
- Careful cleaning
- Device test run
- Creation of a maintenance report

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 and test certificates

## 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](http://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](http://www.sartorius.com)