

# Test Protocol Rotation Speed, Timer, Temperature

## Laboratory Centrifuge



### Benefits

- Consistent quality of measurement results
- Documented device performance
- Trouble-free operation
- Easy-to-read certificates

### Product Information

It is imperative in ISO quality management systems that a test equipment management system be in place, and therefore that laboratory devices are calibrated regularly. Our Sartorius Service has the right calibration for your needs. The test protocol of rotation speed, timer and temperature documents the measurement results in an easily readable certificate.

## Specification

- Calibration of rotation speed:
  - Calibration at two measuring points according to Sartorius testing instructions
- Calibration of timer:
  - Calibration at two measuring points according to Sartorius testing instructions
- Calibration of temperature
  - Test according to Sartorius testing instructions
- Documentation of test results in an easy-to-read test protocol

## Customer Prerequisites

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

## 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)