

Test Protocol Temperature Test Protocol Reproducibility Moisture Analyzers



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

- Test protocol temperature:
 - Test according to Sartorius testing instructions
 - Check with calibrated temperature adjustment set
 - Documentation of test results in an easy-to-read test protocol
- Test protocol reproducibility:
 - Test according to Sartorius testing instructions
 - Reproducibility check with saline solution
 - 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

Optional Services

- Calibration certificate with ISO | IEC 17025 accreditation for the weighing system of the moisture analyzer

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