

# Installation

## Moisture Analyzers



### Benefits

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

### Product Information

The reliability and service life of your moisture analyzers 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 moisture analyzer 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
- Connection of peripheral devices (such as printers)
- 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

- Installation report
- Calibration certificate with ISO | IEC 17025 accreditation
- Temperature adjustment with protocol
- Reproducibility test with protocol
- 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](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)