

# Installation

## MD8 Air Sampler



### Benefits

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

### Product Information

The reliability and service life of your MD8 Air Samplers 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 MD8 Air Sampler is immediately ready for operation and gives reliable and reproducible results right from the start.

## Specification

- Checking the scope of delivery for completeness and any transport damage
- Installation of the device at the place of use
- Equipment installation
- 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

- GLP | GMP compliant equipment qualification
- Calibration certificate (airflow rate and sampling time)

## 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)