

# Equipment Qualification (IQ | OQ)

## Arium® Water Purification Systems



### Benefits

- Compliance with regulatory requirements (GLP | GMP)
- Risk-based verification of device performance
- Proof of suitability for your devices
- Extensive documentation and certificates
- Audit security

### Product Information

You must demonstrate proof of suitability for the equipment in your GLP | GMP regulated laboratory. The Sartorius equipment qualification is documented proof that your water purification system has been installed correctly, works reliably and is suitable for use. Our IQ | OQ qualification follows the "risk-based approach", as recommended for equipment inspections in regulated industries. Specially trained Sartorius service engineers perform all the necessary tests directly on-site at your premises. The comprehensive and customer-specific documentation can easily be integrated into existing quality management systems.

## Specification

- On-site qualification (IQ | OQ) by a specially trained service technician
- Performance of device-specific tests for the Installation Qualification (IQ). This includes, among others:
  - Checking the equipment delivery for completeness and any transport damage
  - Documentation of installation conditions
  - Verification of correct installation of equipment components
- Performance of device-specific tests for the Operational Qualification (OQ). This includes, among others:
  - Function tests
  - Initial metrological equipment check
  - Training of the operating personnel
- Complete documentation in special IQ | OQ templates
- Consultation on device maintenance and maintenance cycles

## Customer Prerequisites

- Device is at the installation site and has been sufficiently acclimatized
- The device is freely accessible
- To wall-mount the device, the included wall bracket must be pre-installed by the customer
- Supply lines available at the installation site, max. 1 m away from the location of the system
- Feed water pressure according to device specifications
- Atmospherically vented drain with odor trap (max. 1 m away from the location of the system)
- Splash-proof wall outlet (240 V, 50 Hz), protected with 16 A fuse, max. 1 m away from the location of the system

## 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)