

Equipment Qualification (IQ | OQ)

Sterisart® Universal Pumps



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 Sterisart® pump 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

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

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