

Digital Equipment Qualification Documentation (IQ/OQ)

Sartorius Digital IQ/OQ offers a paperless solution for equipment qualification, enabling customers to securely approve and sign documents digitally via a web-based platform, making the IQ/OQ process more efficient and sustainable.

1. What is a digital equipment qualification (IQ/OQ)?

It is the digital version of the paper-based equipment qualification documentation (IQ/OQ).

2. How are the digital IQ/OQ documents provided to Sartorius customers?

Sartorius uses an electronic process via a web-based platform to provide the digital IQ/OQ documents. Customers can use this platform to digitally approve, sign, and download the documents.

3. Do I need to purchase or download any software or apps?

You do not need any additional software or apps; standard web browsers can be used.

4. How can I get a digital IQ/OQ?

When ordering an IQ/OQ service, you can decide whether you want the IQ/OQ documentation in paper or digital form at no extra cost.

5. Why is the digital IQ/OQ process a sustainable and efficient solution for laboratory equipment qualification documentation compared to a paper-based documentation?

The digital IQ/OQ process can optimize equipment qualification documentation by providing, approving, signing, and downloading via a single platform, saving paper and resources, which is more sustainable than the paper-based process.

6. Is Sartorius' digital IQ/OQ process secure?

Yes, because it uses electronic signatures in accordance with 21 CFR Part 11, offers secure storage and access

controls, applies the four-eyes principle, and ensures time-limited access to documents.

7. How is a digital IQ/OQ performed and who is involved?

The service technician documents the IQ/OQ activities digitally and makes the documentation available to the customer online via a web-based platform so that they can review, approve, and sign it.

8. How do I access the digital IQ/OQ documents?

You can access the documents via the secure links that will be sent to you by email.

9. How do I sign and approve the documents?

You approve and sign the documents electronically via the web-based platform.

10. How do I save the documents?

After final approval, you should download and save the IQ/OQ documents from the web-based platform, as the links are time-limited.

11. How can I get further information about digital IQ/OQ?

For questions, such as the availability of a digital IQ/OQ in your region, please use the [contact form](#) on our



Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
3874 Research Park Drive
Ann Arbor, MI 48108
Phone +1 734 769 1600