



Simplify your Laboratory  
Processes with Sartorius  
Digital IQ | OQ  
Documentation

Go Digital - Stay Personal

Simplifying Progress

**SARTORIUS**

Whether you are already digital or looking to make the transition – we have the perfect solution for your needs. Discover the innovative digital IQ | OQ documentation as a sustainable alternative to the conventional paper-based method and simplify your laboratory processes.

Thanks to the well-known personal support from your service engineer throughout the entire process, you can switch over effortlessly and at no additional cost. Experience the benefits of innovation with Sartorius.



## Efficient Workflow

Our digital IQ | OQ solution enhances productivity by simplifying document approval and storage, reducing time and effort compared to traditional methods.



## Modern & Sustainable Solutions

Integrate a digital solution that reduces paper usage and optimizes resource management, aligning with your sustainability goals without compromising on quality.



## Secure & Compliant Documentation

Our platform adheres to stringent security standards, including electronic signatures compliant with 21 CFR Part 11, ensuring your documentation is both secure and reliable.



### Ready to simplify your equipment qualification documentation?

Visit our website or contact us directly to check the availability in your region and discover how Sartorius can enhance your laboratory operations.

<https://www.sartorius.com/en/services/instrument-service/lab-balances-weighing-instrument-services/digital-iqoq>

### Germany

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

### USA

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

 **For further information, visit**  
[sartorius.com](https://www.sartorius.com)