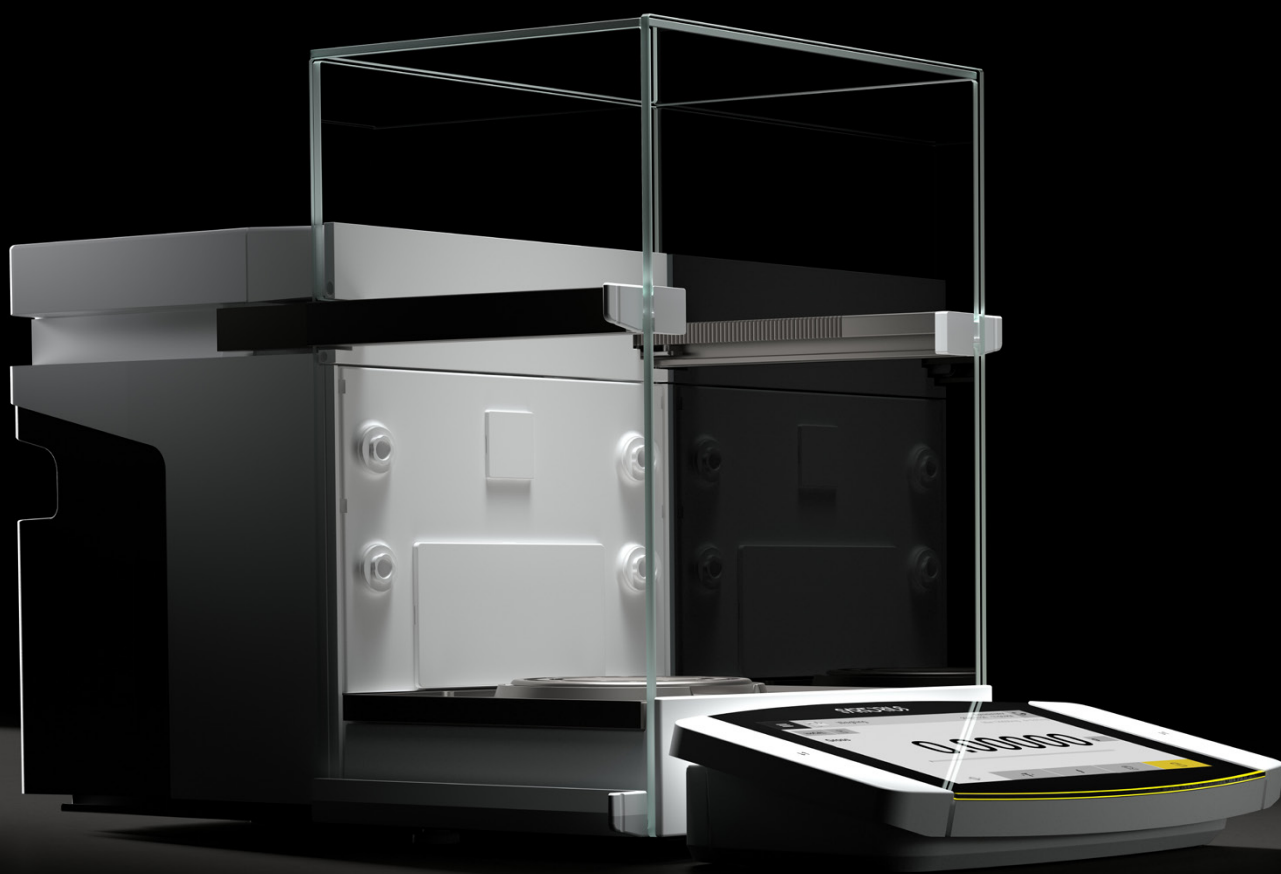


SARTORIUS

Cubis® II meets OMNIS

Seamless Integration
for Greater Efficiency,
Compliance, and Data
Integrity in the Lab





Cubis® II is part of OMNIS

In today's laboratories, precision, efficiency, and regulatory compliance are critical to success. The powerful combination of Metrohm OMNIS and the Sartorius Cubis® II premium laboratory balance delivers an integrated solution for automated sample preparation—from accurate weighing to titration.

This seamless workflow eliminates manual data transfer, saves valuable time, and supports full compliance with 21 CFR Part 11 and EU Annex 11 requirements for data integrity. Thanks to its modular system architecture, the solution adapts to your lab's evolving needs—now and in the future.

Cubis® II and OMNIS—The smart integration for maximum process security in regulated environments

Effortless, Secure Data Handling

More than numbers: Cubis® II sends encrypted results and metadata to OMNIS, ensuring data integrity and compliance.

One system. Full control

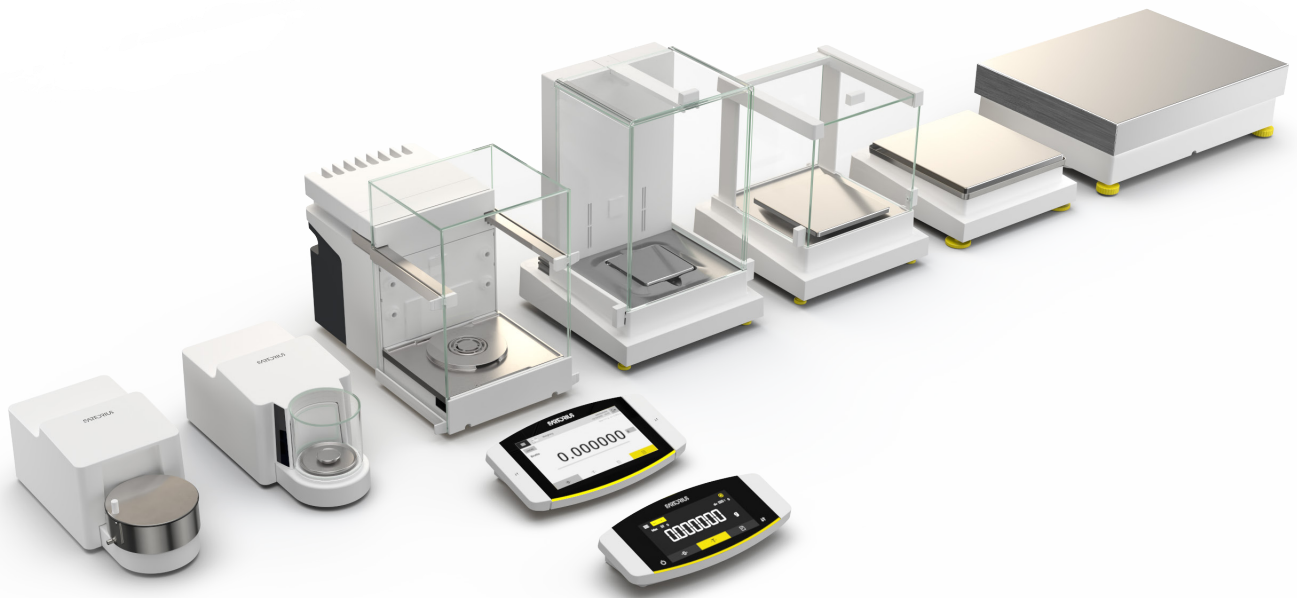
OMNIS centrally manages user access and balance activity with a real-time Audit Trail—fully compliant with FDA 21 CFR Part 11 and Annex 11, effortlessly.

Smart weighing at the balance

Start sample series from Cubis® II without opening OMNIS. Visual guides support weighing, while multiple OMNIS clients share one balance for efficient workflows. Cubis® II balance.

Discover the benefits from the integration of Cubis® II into the OMNIS world

	Cubis® II	Non-Cubis® Balances
Transfer of Weighing Values	Yes	Yes
Transfer of Metadata and Audit Trail	Yes	No
Plug-and-Play Integration	Yes	No
Balance Sharing with Multiple OMNIS Systems	Yes	No
Seamless Workflow Support	Yes	Limited
Operation Directly on the Balance's Display	Yes	No
Centralize User Management	Yes	No
Full Compliance	Yes	No



Cubis® II High-End Laboratory Balance

The New Generation of Modular Precision, Analytical and Microbalances

For scientists in R&D and analytical laboratories needing the most reliable lab weighing results, the Cubis® II is a completely configurable, high performance portfolio of balance hardware and software that will align with your unique demands and compliance requirements to maximize operational efficiency and experimental outcomes.

All modules of the Cubis® II balance were designed for an intuitive operation, further aided by intelligent diagnostic systems. These design elements guarantee a higher degree of repeatability for different workflows while lowering the probability of human error during measurement.

Trusted Results

The Cubis® II balance offers several advantages over other lab balances:

Leading Performance

Leading Performance The Cubis® II balance offers best-in-class performance, including high- capacity and high readability, regardless of sample type

Error-Free Operations

Ensure a high level of trust and transparency by using Cubis® II's functionality to assess relevant factors during data acquisition.

Full End-to-End Data Integrity

Ensure your data is reliable and cannot be manipulated with Cubis® II's user management function and LIMS connectivity.

Reliable Support

Our lab weighing team provides expert support to help ensure your success, even when making very specific and/or unique requests.

Discover More and Request your Demo



USA

Sartorius Corporation
3874 Research Park Dr.
Ann Arbor, MI 48108
Phone +1 734 769 16006

Germany

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



For further contacts, visit
www.sartorius.com