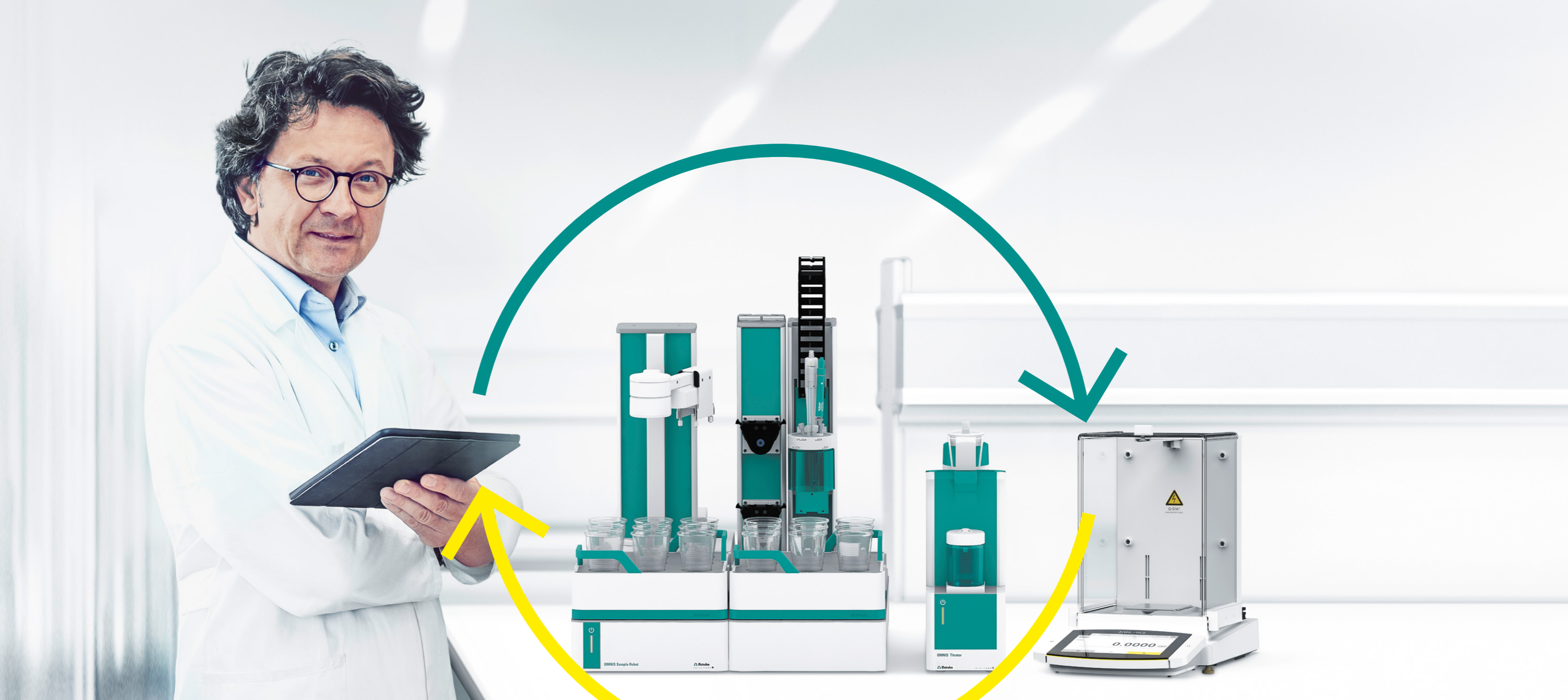


SARTORIUS

Cubis[®] II meets OMNIS

Seamless Integration into
OMNIS, the Analytical
Platform from Metrohm





Cubis® II is now part of OMNIS

With the new software release (available in May 2020) Cubis® II now is a full member of the Metrohm OMNIS analytical platform. OMNIS is a user- and application friendly analytical software platform, focusing on what matters (sample-batch-approach), reducing errors (self-monitoring, alerts and system checks) and ensuring a high level of regulatory compliance (21 CFR Part 11; EudraLex, Volume 4, Annex 11).

Data security and integrity

Cubis® II not only sends the measurement results to OMNIS, but also the complete metadata for every weighing-in or weighing-out performed on the balance. When integrated in a local network, encryption of any data exchanged makes sure the highest standards of data integrity and security are met at all times.

Central user management

The user administration for the Cubis® II balance is managed and verified centrally in OMNIS. The complete Audit Trail of all events on the balance is recorded in OMNIS. All requirements regarding regulatory compliance (FDA and EudraLex) are fully met.

Maximum efficiency

The weighing-in or weighing-out of a sample or even of a complete sample series can be started right from the balance – without the need to even start OMNIS. Tolerance limits regarding sample weights are graphically represented on the display of the balance. When part of a local network, several OMNIS Clients can «share» the same Cubis® II balance.



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 from Sartorius 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**
 The Cubis® II balance offers best-in-class performance, including high-capacity and high readability, regardless of sample type
- 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.
- Error-free Operations**
 Ensure a high level of trust and transparency by using Cubis® II's functionality to assess relevant factors during data acquisition.
- Reliable Support**
 Our lab weighing team provides expert support to help ensure your success, even when making very specific and/or unique requests.

Sales and Service Contacts

For further contacts, visit sartorius.com

Germany

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

USA

Sartorius Corporation
 5 Orville Drive, Suite 200
 Bohemia, NY 11716
 Phone +1 631 254 4249
 Toll-free +1 800 635 2906