

Göttingen, Germany | June 3, 2026

Cubis® III: Compliant weighing, direct integration, and scalable connectivity, without extra software

- Weighing data sent directly and securely into existing lab systems, including LIMS, ELN, and CDS, without extra software
- Built-in audit trails, electronic signatures, user access controls, and guided weighing workflows help labs meet regulatory requirements
- Simplified validation, maintenance, and fleet management help reduce IT effort and control long-term costs

Sartorius launches Cubis® III, the latest innovation in its premium line of lab balances, delivering compliant, connected weighing without relying on extra software typically required to bridge instruments and lab systems. With built-in connectivity, native compliance features, and configurable on-instrument workflows, Cubis® III integrates directly into any digital ecosystem, simplifying validation, IT effort, and long-term maintenance.

“For compliance-driven labs, a balance is a critical data source. But in many setups, meeting regulations depends on an additional middleware layer,” said Eszter Paldy, Manager of Product Management Premium Weighing Solutions at Sartorius. “With Cubis® III, compliance and connectivity are handled on the instrument, so labs can choose how they want to integrate and connect directly from the device.”

Cubis® III delivers direct, standards-based connectivity so weighing data can flow securely into existing informatics environments. An embedded OPC UA server and support for enterprise protocols such as HTTPS, FTPS, and SMB enable direct integration into LIMS, ELN, and CDS without custom, point-to-point solutions. Role-based access control and LDAP integration bring enterprise IT practices to the device, supporting secure user management across balance fleets.

Compliance features are fully embedded in the instrument. On-instrument audit trails and electronic signatures support 21 CFR Part 11 requirements, along with guided, pharmacopeia-aligned weighing workflows that enforce consistency across sites and users. Digitized cleaning workflows (Cleaning QApp) capture cleanings with authenticated sign-off, improving documentation and traceability without additional paperwork.

By removing the dependency on standalone middleware, Cubis® III reduces the scope of systems that must be installed, qualified, and maintained, streamlining lifecycle management and making it easier to scale balance fleets over time. High accuracy with small sample quantities also supports efficient use of valuable materials.

With two packages, Cubis® III and Cubis® III Max, Sartorius enables labs to choose the level of data handling and workflow sophistication that best fits their needs today, while retaining the flexibility to expand capabilities over time through additional workflows and integrations. For labs facing growing demands around compliance, data governance and cost control, Cubis® III lab balances provide a scalable foundation for a more connected, efficient and future-ready lab.

More information about Cubis® III:

www.sartorius.com/cubis-III

A profile of Sartorius

Sartorius is a leading international partner for life sciences research and the biopharmaceutical industry. With innovative laboratory instruments and consumables, the Group's Lab Products & Services Division focuses on laboratories performing research and quality control at pharmaceutical and biopharmaceutical companies as well as academic research institutes. The Bioprocess Solutions Division, with its broad product portfolio focusing on single-use solutions, helps customers manufacture biotech medications, vaccines, and cell and gene therapies more safely, rapidly, and sustainably. Based in Göttingen, Germany, Sartorius has a strong global reach with around 60 production and sales sites worldwide. Sartorius regularly expands its portfolio through the acquisition of complementary technologies. In 2025, a sales revenue of around 3.4 billion Euros was generated with more than 14,000 employees working for customers around the globe.

For more information visit our [Newsroom](#) or follow us on [LinkedIn](#).

Contact

Attila Aranyos

Head of Marketing

Lab Products & Services Division

+33 673 26 59 51

Attila.Aranyos@Sartorius.com