

Göttingen, Germany | September 30, 2025

Introducing the Cubis® II Pipette Calibration System: a complete game-changer for fast, ISO-compliant calibration of single channel pipettes

- One instrument for fast, reliable, and ISO-compliant calibration of pipettes ranging from 2 to 5 mL
- Immediate efficiency gains with reduced instrument fleet costs and increased throughput
- Fast and accurate measurements, and enhanced usability due to larger receiving vessel and durable plastic beaker

Sartorius announces the launch of the Cubis® II Single Channel Pipette Calibration System, integrating the capabilities of multiple systems into a single instrument for fast, reliable, and ISO-compliant calibration for pipettes ranging from 2 to 5,000 µL. Designed to transform workflows, this streamlined solution delivers immediate cost and efficiency gains for on-site calibration service providers.

“Pipette calibration is crucial for research labs, yet providers of this essential service are constrained by inefficient setups that require multiple systems to calibrate one set of pipettes,” says Lucas Foerster, Product Manager Lab Weighing at Sartorius. “The Cubis® II calibration system provides ISO-compliant calibration for 95 percent of all single-channel pipettes, eliminating the need for users to switch between multiple instruments. This also means fewer instruments to purchase, manage and transport to the customer site. It’s a complete game changer.”

Powered by Sartorius’ industry-trusted Cubis® II weighing technology, the calibration system ensures accurate and rapid measurements in any lab environment. Its larger receiving vessel reduces emptying frequency, while the durable plastic beaker prevents breakage, streamlining the calibration workflow and minimizing interruptions. Advanced motion sensorics and an automated lid further enhance the user experience during repetitive calibration tasks.

Sartorius continues to foster innovation in the life sciences sector, supporting researchers with state-of-the-art solutions that drive efficiencies. Supporting the pipette calibration workflow are the integrated QApp and the Ingenix Suite software to simplify compliance with the latest regulatory requirements. Laboratories can also upgrade their existing Cubis® II balance into a Single Channel Pipette Calibration System using the new YCP01MPS accessory, maintaining the versatility of a universal, premium lab balance.

More information about this product

<https://www.sartorius.com/en/products/weighing/pipette-calibration>

A profile of Sartorius

Sartorius is a leading international partner to 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, the company 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 2024, the company generated sales revenue of around 3.4 billion euros. More than 13,500 employees work for customers around the globe.

Visit our [Newsroom](#) or follow us on [LinkedIn](#).

Contact

Lisa Dabney

Head of Strategic Marketing

Lab Products & Services Division

+1 734-474-7117

lisa.dabney@sartorius.com