

Göttingen, February 27th, 2024

The Fastest Stabilization Time Meets Everyday Analytical Weighing

- Fastest stabilization time in its class for analytical weighing; results in under 2 seconds
- Low minimum sample weight starting at 4.1 mg sets new benchmark for semi-microbalances
- Innovative Cleaning QApp and tool-free assembly simplify compliance and routine maintenance
- Advanced ionizing technology fully eliminates static electricity, optimizing accuracy

The Life Science Group Sartorius introduces the Cubis® II Ultra-High Resolution Semi-Micro balance, a perfect addition to the Cubis® II portfolio of premium laboratory balances. Designed for a wide range of weighing applications, this balance is the quintessential instrument for everyday use, offering impressive speed, precision, and ease-of-use.

Equipped with five-digit readability and a 220g capacity, the balance integrates cutting-edge weighing technology and a low minimum sample weight of 4.1 mg (or 8.2 mg for some models), surpassing typical requirements. Additionally, its industry-leading stabilization time of less than 2 seconds significantly enhances efficiency in the field.

Automatic compensation for environmental changes, such as temperature, humidity, air drafts, and air pressure, ensures consistent performance under real-world conditions. Crucial to this achievement is an innovative ionization technology that effectively neutralizes static electricity.

Recognizing the importance of cleanability, the new Cleaning QApp dramatically simplifies routine maintenance with clear, step-by-step instructions. Tool-free disassembly and reassembly of balance components further facilitate the cleaning processes, enabling compliance with standard operating procedures.

"The Cubis® II Semi-Micro balance is a game-changer for labs looking for a balance that excels in everyday weighing tasks," says Eszter Paldy, Manager of Product Management Premium Weighing Solutions at Sartorius. "With its fast stabilization time, low minimum sample weight, and ease of cleaning, this balance is a real solution that adapts to the dynamic needs of modern labs."

The new balance is fully equipped with the Cubis® II family's standard features, including modularity, compliance, connectivity, and digital data management.

For more information, please visit:

[Cubis® II Ultra-High Resolution Semi-Micro balance](#)

[Premium Balances from Sartorius](#)

About Sartorius

Sartorius is a leading international partner of life sciences research and biopharmaceutical manufacturing. 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 safely, rapidly, and economically. The company, based in Göttingen, Germany, 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 2023, the company generated sales revenue of around 3.4 billion euros. Currently, around 14,600 employees are working for customers around the globe.

Contact

Lisa Dabney

Head of Marketing Programs LPS

+1 734-474-7117

Lisa.Dabney@Sartorius.com

Follow Sartorius on [LinkedIn](#).