

Göttingen, April 3, 2024

Sartorius Unveils Quintix® Pro: Redefining the Standard for Laboratory Balances

- A high-resolution, 7-inch touchscreen enhances the user experience and simplifies workflows
- Customizable user roles and an extensive application portfolio enable seamless integration into diverse lab environments
- Automatic, internal motorized leveling ensures the highest level of weighing accuracy
- Energy-efficient operation modes and a digital user manual significantly reduce environmental impact

The life science group Sartorius announces the launch of the new Quintix® Pro laboratory balance, a groundbreaking innovation that sets a new benchmark in standard laboratory weighing. Designed to meet the evolving requirements of modern laboratories, it excels in usability, flexibility, and sustainability.

Merging sophisticated engineering with premium features, the Quintix® Pro delivers an unmatched user experience. Its intuitive 7-inch, high-resolution, full-touch graphic display streamlines menu navigation and enhances productivity. The series is further differentiated by its automated motorized leveling system, designed to simplify daily operations, and deliver consistently accurate results.

"Our goal was to transform the user experience with a balance that embodies smart, progressive upgrades," explains Christian Weidner, Manager of Product Management Universal Weighing Solutions at Sartorius. "This balance is not just aesthetically pleasing; it also addresses everything that standard balances have been missing. The vivid touch display greatly enhances ease of use, while the motorized leveling provides both accuracy and peace of mind at the touch of a button."

Flexibility is at the heart of the Quintix® Pro. It incorporates 15 application programs and an extensive language selection, making it a true universal weighing instrument for a wide array of workflows. The advanced user management tools offer access controls and customization for tailoring to individual needs. High-quality, durable materials are used in the housing to facilitate adherence to stringent cleaning protocols.

This launch showcases Sartorius' dedication to environmental responsibility, featuring energy-saving modes, automatic dimming, and a user-defined automatic start timer that significantly lessen the ecological footprint. The digital user manual, accessible directly from the display, reduces paper usage and provides immediate access to essential information.

The Quintix® Pro replaces the successful Secura® and Quintix® product lines, as an eco-friendly, cost-effective balance that integrates advanced capabilities traditionally reserved for premium models. It is the quintessential choice for laboratories committed to environmental stewardship without compromising on quality or functionality.

For more information, please visit:

<https://www.sartorius.com/en/products/weighing/laboratory-balances/quintix-pro>

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

Lab Products & Services Division

+1 734-474-7117

Lisa.Dabney@Sartorius.com

Follow Sartorius on [LinkedIn](#).