

Göttingen, Germany, September 18, 2023

## The New Picus® 2 Sets a New Standard for Connected Electronic Pipettes

- Open connectivity enhances lab productivity and sets a new standard for electronic pipettes
- Fully electronic pipetting and companion app ensure reproducibility across workflows
- Unbeatable ergonomics and intuitive user interface transform the pipetting experience

The Life Science Group Sartorius introduces the Picus® 2 Electronic Pipette, the latest addition to its portfolio of electronic pipettes. With a combination of advanced connectivity and excellent performance, Picus® 2 sets a new standard, enabling efficient, repeatable dispensing across biopharmaceutical and QC applications.

Pipettes play an integral role in the connected lab by enabling real-time data recording and integration with other equipment, ultimately enhancing productivity and data quality. Picus® 2 electronic pipettes are equipped with Bluetooth connectivity, allowing seamless and flexible connection to any data management system. A unique offering with Picus® 2 is the Sartorius Pipetting app for mobile devices, thoughtfully designed as a comprehensive solution for pipetting workflows.

"With Picus® 2 we created a fully connected electronic pipette that transforms the pipetting experience for our users," said Joni Åke, Product Manager of Liquid Handling, Sartorius. "Thanks to its open connectivity, customers can either integrate Picus® 2 with their existing system, or pair it with the companion Sartorius Pipetting app for access to our expansive library of configurable protocols, step-by-step guides, software updates, and other pipette management tools that simplify daily progress."

Building on the success of its predecessors, Picus® 2 features its trademark unbeatable ergonomics, wide range of pipetting modes, and precise electronic pipetting. The top-to-bottom ergonomic construction, with soft-touch operation and electronic tip ejection, not only improves overall comfort and efficiency, but also reduces the risk of injury. Large digital display and intuitive menus enable easy access to everyday operations, while secure password protection and programmable calibration reminders support compliance needs.

"We designed Picus® 2 for maximum usability and comfort, two critical factors for getting reliable pipetting results," said Sandra Söderholm, Application Development Scientist for Lab Essentials, Sartorius. "Despite its ability to perform complex functions, Picus® 2 is as easy to operate as a mechanical pipette. Picus® pipettes are also synonymous with exceptional ergonomics, which promotes effortless, error-free pipetting. The ergonomic design, ease of use, and flexible connectivity work together to ensure accuracy and reproducibility in liquid handling workflows."

The launch of Picus® 2 further establishes Sartorius as a leading supplier of premier lab essentials that streamline processes and accelerate research in the modern, connected laboratory. Picus® 2 pipettes are available in single-channel models, covering a volume range of 0.2 µL to 10,000 µL, and in multichannel

models from 0.2 µL to 1,200 µL. Sartorius also supplies high-quality standard and specialized pipette tips that are validated for use with Picus® 2 in a wide range of applications.

#### Further information

- [Picus® 2 electronic pipette](#)
- [Pipettes and pipette tips by Sartorius](#)

#### A profile of Sartorius

The Sartorius Group is a leading international partner of 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 and vaccines safely, rapidly and economically. The company, based in Goettingen, Germany, has a strong global reach with around 60 production and sales sites worldwide. Sartorius delivers significant organic growth and regularly expands its portfolio through the acquisition of complementary technologies. In fiscal 2022, the company generated sales revenue of around 4.2 billion euros. At the end of 2022, around 16,000 employees were working for customers around the globe.

#### Contact

Lisa Dabney  
Head of Marketing Programs LPS  
+1 734-474-7117  
[Lisa.Dabney@Sartorius.com](mailto:Lisa.Dabney@Sartorius.com)

Follow Sartorius on X (Twitter) [@Sartorius\\_Group](#) and on [LinkedIn](#).