



Picus[®] 2

Sustainability Fact Sheet

Simplifying Progress

SARTORIUS

Overview

Sartorius Picus® 2 electronic pipettes are the ideal tools to accompany you in daily, repetitive liquid handling in laboratories, where the accuracy of your results, speed and ergonomics are important. Their fully electronic control guarantees consistent, user-independent results, and their lightweight, ergonomic design gives you total convenience.

Life Cycle Thinking

Adopting life cycle thinking is key to enhancing sustainability and considering the environmental impacts from raw materials to end-of-life. We continuously refine our design processes and manufacturing methods to increase efficiency, reduce waste, and lower energy consumption. Furthermore, we consider the implications of shipping and strive to optimize logistics to minimize carbon emissions.



Manufacturing

Our pipette manufacturing plants in Helsinki and Kajaani, Finland operate with 100% renewable electricity. The waste heat from production is redirected to provide heating for the manufacturing or storage facilities, thereby reducing the consumption of district heat. Additionally, we have upgraded all the lighting to LED lights.

More than half of Picus® 2 components are manufactured in-house, which means we are controlling more stages of the manufacturing process and its impact to environment. Our plastic component manufacturing total waste recovery rate is 95%.

Sartorius's manufacturing and office facilities in Finland have partnered with a local Carbon Neutral Waste Management service¹. The carbon footprint produced by our waste management is measured and the CO₂ emissions are offset through Gold Standard certified afforestation projects. In 2023, we invested in permanent carbon sinks to neutralize our waste emissions. Alongside diligent waste sorting and reduction efforts, this service helps to counteract any remaining emissions, playing a vital role in our strategy to reduce our operational carbon footprint.

Our quality and environmental management system is certified by independent third party to meet ISO 9001, ISO 13485, and ISO 14001 standards.

Learn more about carbon neutral waste service (in Finnish): Hiilineutraali jätehuolto | L&T



Sartorius manufacturing plant in Kajaani, Finland



Product Design and Use

Long-lasting products reduce the need for frequent replacements, thereby conserving resources and minimizing waste. Pipettes are typically used for many years, commonly more than 10 years can easily be achieved with a proper maintenance program. Designing Picus® 2 pipettes to be easily maintainable and repairable extends their usable life and prevents unnecessary disposal. High quality materials and manufacturing technologies ensures stable performance of the product year after year. Sartorius also is committed to secure and long-term supply of spare parts for our pipettes.

A pipette and its corresponding tip constitute an integrated system, designed to function seamlessly together. When Sartorius pipettes are paired with either Optifit tips or Safetyspace® filter tips, they achieve their full intended performance. The compatibility between the pipette and tip and the overall quality of the manufacturing process is paramount for ensuring reliable, reproducible results. This in turn minimizes the necessity for unnecessary pipetting repetitions thereby reducing the overall usage of pipette tips, as well as the liquids pipetted.

The tips, a vital part of your daily work, generate significant portion of the total environment impact when using Picus® 2 pipette. For a deeper understanding of our approach to offer more sustainable pipette tips, we invite you to read more from the Sartorius pipette tip fact sheet. Make a choice that benefits both your research and the environment with Sartorius tips.

Packaging Design and Distribution

Packaging has a critical role in maintaining the integrity and quality of the products during transport. Packaging safeguards pipettes from potential damage and degradation while ensuring that they reach end users in pristine condition. By minimizing the packaging size, we can streamline logistics, improve storage efficiency, and reduce the material usage.

Picus® 2 packaging is made with 100% FSC certified cardboard.

After Sales Service

Longevity and reliability of pipette are pivotal to your success. That's why we are committed to providing repair, maintenance, and calibration services. Extend the life of your instruments with our expert technicians who are dedicated to keeping your instruments in peak condition. Regular maintenance can prevent unexpected breakdowns, extend the lifespan of your equipment, and save you from the cost and environmental impact of frequent replacements.

Regular calibration for optimal accuracy and repeatability is the cornerstone of quality. Our accredited calibration services guarantee that your instruments perform to their original specifications, ensuring consistent results and reducing the likelihood of errors that could lead to waste. By optimizing the functionality and extending the life of your instruments through spare part availability and our services, you're actively reducing waste and conserving resources.

In summary, product service plays a crucial role in enhancing product sustainability by promoting longevity, efficiency, and responsible end-of-life management, all of which contribute to reducing the overall environmental impact of our products.

End-of-Life

The product sales unit packaging that is made from cardboard, is fully recyclable. Disposing electronic devices, like electronic pipettes, should be done according to local guidelines. For least environmental harm you should look for a solution that separates components like plastics, metal, electronics, and battery, and that the components are not glued together. Picus pipettes are easy to disassemble for recycling. Sartorius service centers can help you with recycling electronic pipettes.

Waste from Electrical and Electronic Equipment (WEEE) must be sorted and disposed accordingly. This enables recovery and recycling of rare and valuable raw materials and prevents hazardous materials from ending up in the environment.

To minimize electrical waste, Picus® 2 pipettes are delivered with a universal USB charging cable. Other charging options like charging stands and adapters can be ordered separately as needed.

Sustainability at Sartorius

Sartorius is committed to contributing to a future where more people have access to better medicine. At the same time, we take on responsibility for the impact of our operations, wherever they occur.

Taking into account the concerns of its stakeholders, Sartorius has defined six strategic sustainability topics:



Climate Action



Resources and
Circularity



Water and Effluents



Supply Chains



Social Responsibility



Corporate
Governance

For more information, please visit:
www.sartorius.com/en/company/sustainability



Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

Singapore

Sartorius Singapore Pte. Ltd
Mapletree Business City
30 Pasir Panjang Road
#06-31A/32
Singapore 117440
Phone +65.6872.3966



For further information, visit,
www.sartorius.com