

## Toward Less Waste, Less Plastic, and a More Sustainable Future

### Our Commitment

At Sartorius, our commitment and responsibility to sustainability is non-negotiable. From design and manufacturing to use and disposal, our innovative laboratory solutions are thoughtfully designed to minimize environmental impact. Because we believe even small changes make a big difference.

### What We're Doing

We are actively reducing waste and plastic use through several initiatives. These include rethinking product design, using renewable materials, and implementing reuse strategies. As a result, Sartorius is delivering measurable progress in waste and plastic reduction across its Lab Essentials portfolio.

### Our Focus



#### Circular Materials:

Using renewable and recycled materials, including ISCC PLUS-certified plastics and aluminum with recycled content.



#### Minimized Packaging Waste

Reducing plastic content, weight and complexity, and using mechanically recycled materials.



#### Innovative Product Design

Developing lab tools that last longer, can be maintained easily, and support reuse where possible



#### Optimized Distribution

Improving packaging and pallet design to reduce shipping volume and lower transport-related emissions

### Sustainability at a Glance

Goal	Progress
Plastic Reduction	Up to 70% less in pipette tip refill packaging
Product Reuse	95% less waste with Biosart® 250 Funnels
Renewable Electricity	100% in key facilities (Finland, Germany)
Waste Recovery in Production	Up to 95% for plastics (Picus® 2)
Water & Energy Efficiency	50% fewer cleaning cycles (Arium® iJust)
Recyclability & Materials	Recycled content in packaging and parts

# Products



## Vivaflow® SU Filtration Cassettes

- 30.5% less plastic than previous models
- ISCC PLUS-certified renewable polycarbonate
- Replaced plastic bubble wrap with cardboard inserts
- Tubing kits tailored to reduce unused parts



## Pipette Tips

- Up to 70% plastic reduction in Refill Tower and FlexiBulk tip products
- Bulk sourcing and storage of raw material



## Quintix® Pro Standard Lab Balance

- Baseplate made with 20% recycled aluminium
- 70% of the packaging is made from recycled materials
- Power-saving mode reduces energy use by over 50% during standby



## Sartolab® RF|BT Vacuum Filtration

- 22% reduction in plastic film for 150–250 mL products
- Optimized product and pallet design reduces total plastic per shipment
- 40% thinner pallet stretch wrap used in shipping



## Biosart® 250 Funnels

- Reuse reduces total plastic waste by up to 95%
- Built from durable polypropylene
- Up to 50 autoclave cycles per funnel



## Picus® 2 Electronic Pipette

- Built for long-term use and repairability
- 95% plastic waste recovery in production
- FSC-certified recyclable packaging
- 100% renewable electricity used in production



## Arium® Lab Water System

- Up to 50% fewer cleaning cycles per year with iJust software
- Built for longevity, with replaceable and repairable parts
- Take-back program for UV lamps in certain regions

## What's Next

We are continuously rethinking and redesigning our solutions while developing new innovations that support both laboratory performance and environmental responsibility.

Our focus is on helping labs reduce their environmental footprint by promoting resource efficiency and enabling circular practices. This includes minimizing operational waste, enhancing product recyclability, and reducing reliance on fossil-based materials.

Sartorius is aiming for climate neutrality by 2045 and transitioning to 100% renewable electricity by 2030.