



Biosart® 250 Funnel

Sustainability Fact Sheet

Simplifying Progress

SARTORIUS

Biosart® 250 Funnel:

Sustainable and Efficient Filtration Solution

Reuse of Biosart® 250 Funnel for a Precise and Eco-Friendly Filtration

The Biosart® 250 Funnel is designed for efficient microbiological testing of pharmaceuticals, food, beverages, and water. It allows for quick performance of the filtration steps required in routine processes and ensures reliable results. These sterile funnels can be used with Sartorius gridded membrane filters placed onto the stainless-steel bases of a vacuum manifold. Biosart® 250 Funnels offer the additional benefits of reduced waste and resource savings through the option for reuse.

At a Glance

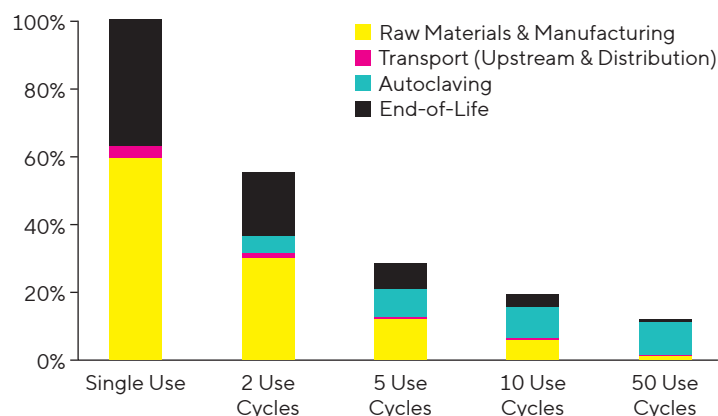
- Sterile, ready-to-use 250 mL plastic funnels, compatible with Sartorius vacuum manifolds and membrane filters
- Autoclavable up to 50 times
- Possesses elasticity due to its polypropylene (PP) construction, ensuring optimal sealing with a bayonet-type closure
- Marked graduations at 50, 100, 150, 200, and 250 mL for a convenient volume measurement
- Large inner diameter ensures a high flow rate. No liquid retaining due to its conical shape



Relative Reduction of Carbon Footprint Single vs. Reuse of Biosart® 250 Funnel

Reusing Biosart® 250 Funnels can significantly reduce the resource demand and carbon footprint compared to a single-use application. Biosart® 250 Funnels are made of polypropylene. The analysis below compares different scenarios - single use versus multiple use cycles including autoclaving after each use cycle. The carbon footprint of the Biosart® considers the impact of raw materials, injection molding in Goettingen using renewable electricity, transport of materials to Sartorius and transport of the product to an average customer in Europe. For the end-of-life, the incineration processes including energy recovery were modelled. The electricity demand for autoclaving was measured and considered in the analysis assuming an average European electricity mix.

Relative Reduction of CO₂ Footprint Single vs. Reuse Biosart® 250 Funnel



Sustainability Benefits

- **Reusable:** Autoclavable up to 50 times, reducing resource use and waste by up to 95%
- **Lower Carbon Footprint:** Up to 88% less CO₂ emissions compared to the single-use option
- **Renewable Energy:** Utilizing renewable electricity for cleaning and autoclaving can further reduce environmental impacts

Order Information

Description	Order No.
50 Biosart® 250 funnel packed in bags	16407--25----ALK
50 Biosart® 250 funnel individually packed	16407--25----ACK

For further information and orders, visit [sartorius.com](https://www.sartorius.com)



Sustainability at Sartorius

Sartorius is dedicated to shaping a future where improved medicine is more accessible to many. Concurrently, we acknowledge and address the impacts of our operations globally. 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 Göttingen
Phone +49 551 308 0

USA

Sartorius Corporation
3874 Research Park Drive
Ann Arbor, MI 48108
Phone +1 734 769 1600

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**
[sartorius.com](https://www.sartorius.com)