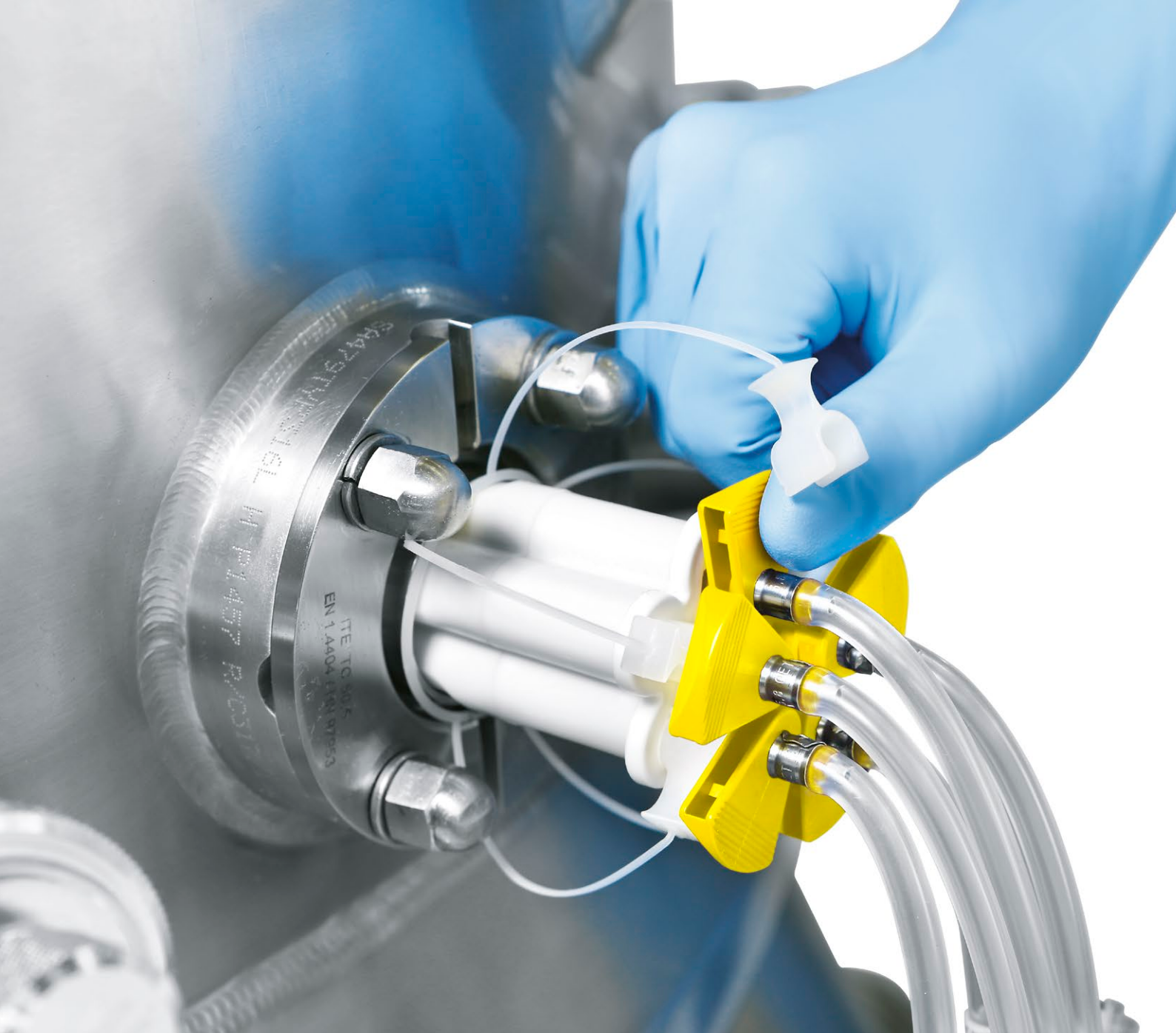




TakeOne® Simplified Sampling Solutions for Stainless Steel Equipment

Simplifying Progress



Sustainability Fact Sheet

SARTORIUS

Overview

TakeOne® aseptic sampling solutions have been designed to integrate seamlessly with stainless steel processes and collect perfectly representative samples with no risk of false-positives for each process step and scale. TakeOne® products are delivered fully assembled, in individual packs, gamma sterilized, integrity tested and ready for immediate use allowing you to bypass parts washing, equipment prep and assembly with confidence.

This fact sheet presents the current sustainability status of TakeOne® aseptic sampling devices, using the 1.5" TC version with Tuflux® TPE tubing. It highlights our commitment to enhancing the sustainability of our products, with continuous improvements being made over time.

Collection containers are considered in separate Sustainability Fact Sheets, for example; Flexsafe® bags for storage and shipping.

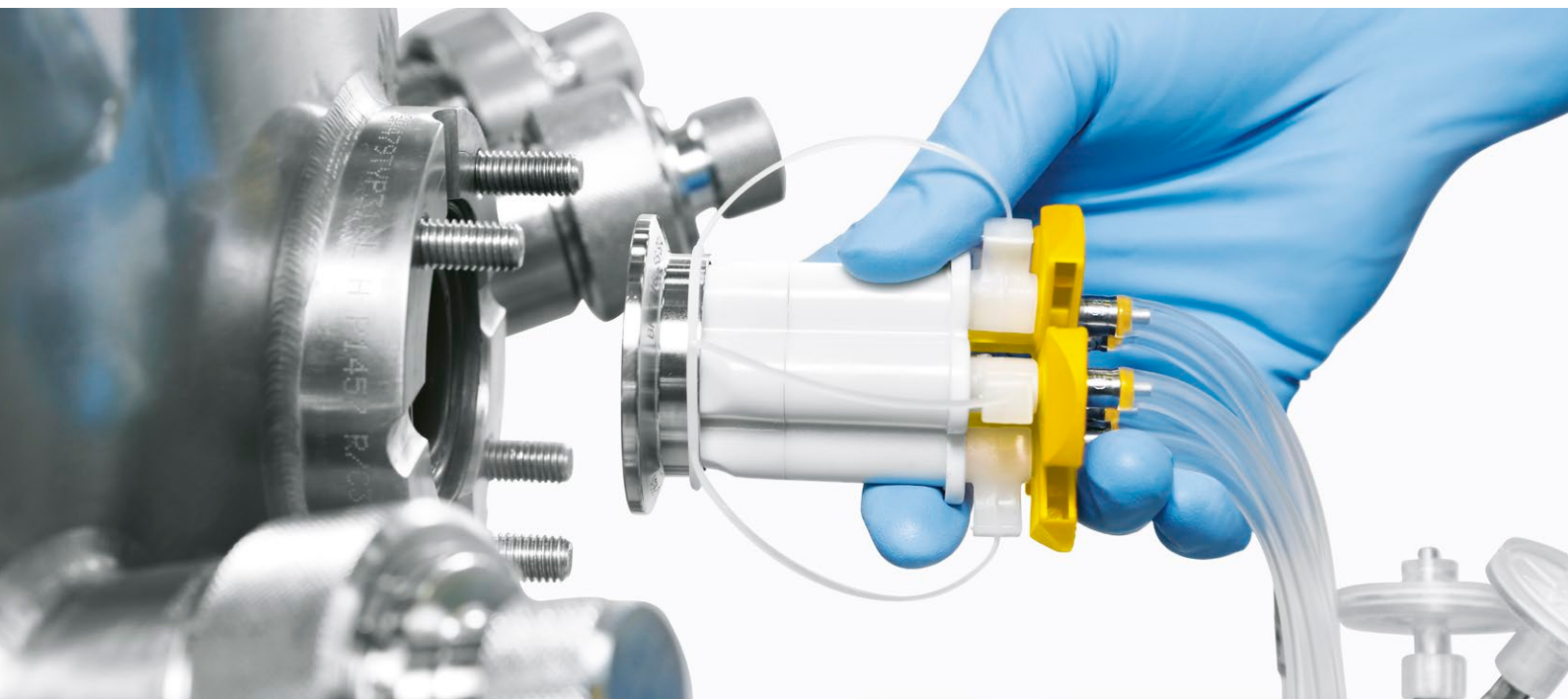
Life Cycle Thinking

At Sartorius, we are committed to sustainability and are actively seeking innovative ways to reduce the ecological footprint of our products.

Adopting life cycle thinking is key to enhancing sustainability and considering the environmental impacts from raw materials to end-of-life disposal. We are dedicated to refining our production methods, boosting efficiency, minimizing waste. We also consider the environmental toll of shipping practices and are committed to optimizing logistics to reduce carbon emissions.

Our ongoing research into materials and designs aims to lessen environmental impact and enhance the recyclability of our products. Guided by Product Carbon Footprint (PCF) screenings, we gain valuable insights that drive the development of more sustainable products and deepen our understanding of their ecological footprints.

Integrating life cycle thinking into our operations not only supports the sustainability of our products but also enables our customers to make environmentally conscious decisions with confidence.



Environmental Overview

Material selection

All materials of construction are selected with care. The polymers used in plastic components and support materials, along with packaging materials like cardboard, are of high quality and largely recyclable.

Raw material acquisition

Raw materials and components are supplied from sources close to the manufacturing sites where feasible. Subassembly manufacturing and final assembly are performed in the same facility in order to reduce the environmental impact of internal transport.

Material processing

Components are manufactured on modern equipment in efficient processes that take material, energy and water consumption into account.

Certification

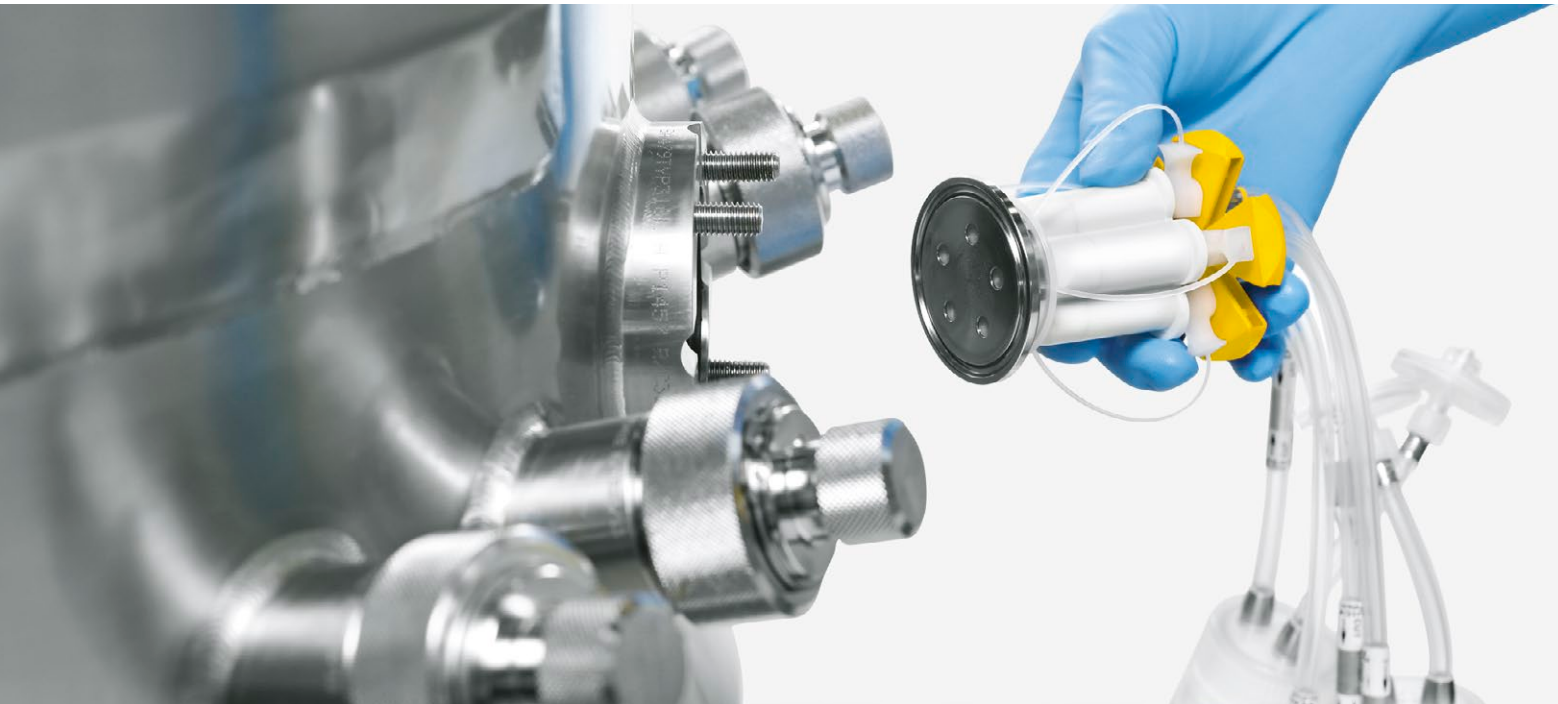
To enhance supply resilience and improve proximity to customers, production is expanding to our ISO 14001 certified sites in M'Hamdia, Tunisia, and Yauco, Puerto Rico.

These certifications demonstrate our commitment to environmental responsibility, ensuring that our environmental management system drives continuous improvement, compliance with legal requirements, and achievement of sustainability objectives.

Distribution

The production and distribution of TakeOne® products is managed with logistics solutions designed to minimize environmental impacts linked to transport to final customers.

By expanding production to additional sites in M'Hamdia, Tunisia, and Yauco, Puerto Rico — closer to key markets — we reduce transportation distances and improve delivery speed. This approach ensures reliable availability while lowering the carbon footprint associated with long-haul transport.



Product

Recyclability	44%
Recycled content	20%

The recyclability of our product stands at 44% and contains 20% of recycled materials, reflecting our commitment to sustainability and environmental responsibility.

Recyclability is defined as the characteristic of products that retain useful physical or chemical properties after serving their original purpose. Once these products are separated and sent to recycling, they can be reintroduced into manufacturing as raw materials¹. In this context, we refer to the technical recyclability of a material or component, meaning that technological solutions exist to recycle them, either mechanically or through advanced recycling methods.

The TakeOne® sampling hubs are designed to facilitate and encourage the recycling and responsible disposal of the used stainless steel sampling hub face plates. For complete guidance, please refer to the Instructions for Use TakeOne® Aseptic Sampling Solutions.



Elements of the product: Options at the end-of-life

Component	Material	Recyclable
Device face plate and faceplate components such as septa	SS, Sil (Pt)	Yes
Faceplate components such as holder	PBT with GF	No
Topcap components such as springs, washer, ring and other	SS, PA	Yes
Topcap components such as top cap and retainer	PBT with GF	No
Cannula and cannula components such as diaphragm	SS, Sil (Pt)	Yes
Cannula components such as thumb tab	PBT with GF	No
Finished article components such as tubings	TPE	No
Finished article components such as plugs, cable ties and clamps	PP, ETFE, SS	Yes

Note. For abbreviations see page 7.

Definitions: ¹Based on European Environmental Agency GEMET – Environmental thesaurus
²Only pre-consumer and post-consumer materials shall be considered as recycled content, in alignment with ISO 14021. The numbers provided herein are the best available approximations.

Disclaimer
The recyclability of the product may be influenced by its use, such as the presence or absence of agents defined as hazardous, as well as local regulations and the capabilities of local companies to manage those materials. Data refers to one TakeOne® aseptic sampling devices, using the 1.5" TC version with Tuflex® TPE tubing.

Packaging

Recyclability	100%
Renewable content	49%
Recycled content	46%

Recyclability¹

Characteristic of products that retain useful physical or chemical properties after serving their original purpose. Once these products are separated and sent to recycling, they can be reintroduced into manufacturing as raw materials.

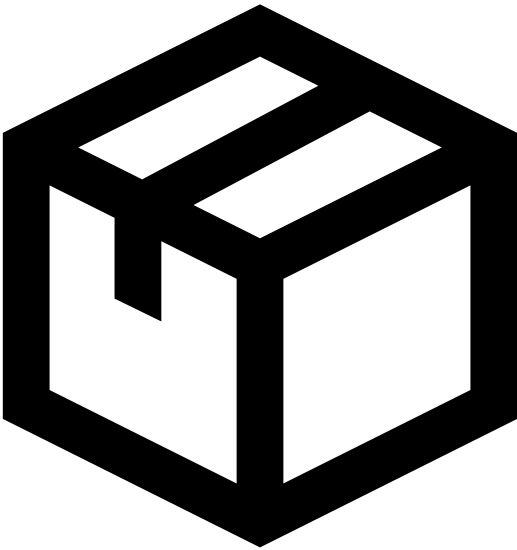
In this context, we refer to the technical recyclability of packaging materials, meaning that technological solutions exist to recycle them, either mechanically or through advanced recycling methods.

Renewable content²

Materials that are derived from resources that are quickly replenished by ecological cycles or agricultural processes, so that the services provided by these and other linked resources are not endangered and remain available for the next generation. In our product packaging, we refer specifically to cardboard and paper.

Recycled content³

Proportion of the mass of recycled material on the total mass of the packaging.



Elements of the primary and secondary packaging: Options at the end-of-life

Category	Packaging element	Material	Recyclable
Plastics	Polybags	LDPE	Yes
Paper and cardboard	Shipping case	Cardboard	Yes

Note. For abbreviations see page 7.

Definitions: ¹ Based on European Environmental Agency GEMET – Environmental thesaurus
² Corporate Sustainability Reporting Directive (CSRD)
³ Only pre-consumer and post-consumer materials shall be considered as recycled content, in alignment with ISO 14021. The numbers provided herein are the best available approximations.

Disclaimer
The recyclability of the packaging may be influenced by local regulations as well as the capabilities of local companies to manage those materials. Data refers to one TakeOne® aseptic sampling devices, using the 1.5" TC version with Tuflux® TPE tubing. As the Packaging and Packaging Waste Regulation (PPWR) finalizes its method for calculating recyclability, Sartorius is committed to adjusting its practices accordingly to ensure compliance and alignment with regulatory standards.

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 further information on sustainability, visit
www.sartorius.com/en/company/sustainability

List of Abbreviations

ABS	Acrylonitrile butadiene styrene
AL	Aluminum
CS	Carbon Steel
EPDM	Ethylene propylene diene elastomer
EVA	Ethylene vinyl acetate
ETFE	Ethylenetetrafluoroethylene
EVOH	Ethylene vinyl alcohol
GF	Glass fiber
HDPE	High density polyethylene
LDPE	Low density polyethylene
LLDPE	Linear low density polyethylene
PA	Polyamide
PBT	Polybutylene terephthalate
PC	Polycarbonate
PE	Polyethylene
PEEK	Polyether ether ketone
PES	Polyethersulfone
PET	Polyethylene terephthalate
PP	Polypropylene
PS	Polystyrene
PSU	Polysulfone
PU/PUR	Polyurethane
PVC	Polyvinylchloride
Sil	Silicone
Sil (Pt)	Platinum cured silicone
SS	Stainless steel
TPE	Thermoplastic elastomer

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Göttingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178



For more information, visit
[sartorius.com](https://www.sartorius.com)