



SafetySpace Filter Tips

Filter Tip to Protect from Contamination

Product Information

Filter tips are an efficient way to avoid cross-contamination of your sample and/or pipette and reduce the need for pipette maintenance. Sartorius SafetySpace Filter Tips have more space between the sample and the filter than conventional filter tips, which prevents the risk of the sample being absorbed into the filter. SafetySpace Filter Tips are ideal for use with Sartorius pipettes, enabling maximum accuracy and precision.



Description

SafetySpace Filter Tips are made of virgin polypropylene and the filter of polyethylene without any self-sealing additives to avoid contamination of the sample, which may affect the results. All the steps in the manufacturing process strictly follow ISO 9001 and ISO 13485 quality standards, and the tips are manufactured in ISO 8-classified cleanroom conditions to avoid any contamination. Every production batch of SafetySpace Filter Tips is tested for DNase, RNase, and endotoxins. The tips are released for sale once they have fulfilled the test requirements. Certificates are available at www.sartorius.com.

The additional space between the sample and the filter differentiates SafetySpace Filter Tips from other filter tips. This feature is particularly useful in reverse pipetting, pipetting foaming liquids (such as buffers and proteins), or when using Multi-Dispensing mode in electronic pipette.

Features

- Available in volumes from 10 to 1200 µl
- All production batches certified RNase, DNase, and endotoxin free
- Informative rack labels indicating tip volume, product number, lot number, date, and purity level for improved tip identification and traceability

- All SafetySpace Filter Tips are E-beam pre-sterilized
- Perfect fitting and sealing with all Sartorius pipettes
- Smooth and ergonomic tip attachment and ejection ensured by compatibility with the spring-loaded Optiload mechanism of Sartorius pipettes
- Color-coded tip trays make it easy to match tip with appropriate Sartorius pipette
- Available in Single Tray flip-top racks that are easy to open and close with one hand
- Materials used in tips, trays, and racks are 100% recyclable or can be incinerated as energy waste
- Universal design provides compatibility with most of the pipette brands

SafetySpace Filter Tips are also available as Low Retention tips. The Sartorius tip offering also includes Optifit non-filter tips and Optifit Low Retention non-filter tips (see catalog for details).

Applications:

- Molecular biology
- Microbiology
- Cell culture applications
- Radioactive work
- Any application where preventing cross-contamination is vital

Intended Use

The Sartorius SafetySpace Filter Tips are intended, designed and manufactured to be used in liquid dispensing in a variety of applications and to be used in combination with Sartorius pipettes. The Sartorius pipette and tip combination, fall within the scope of in-vitro diagnostics, and can be used as a diagnostic medical device in related applications. Thereby, Sartorius pipettes and Sartorius tips fulfil the relevant demands of the Directive 98/97/EC of the European Parliament.

It is recommended that SafetySpace Filter Tips are used with Sartorius pipettes to ensure optimum compatibility and performance.

Technical Data

Technical Specifications

Sterilization	Electron beam (beta irradiation) in accordance with EN 552 and ISO 11137 and validated by using 25 kGy as the minimum dose with SAL (sterility assurance level) of 10 ⁻⁶ .
Purity certificate	Every production batch (lot number) of tips is certified free of DNase, RNase, and endotoxins.
Endotoxin purity testing	Validated test result with LAL gel clot method for endotoxin free pipette tip is < 0,005 IU/ml (EU/ml).*
DNase purity testing	DNase-free pipette tips show no evidence of DNase activity in the fluorometric assay with the detection level of < 6,25*10 ⁻⁵ U/μl when DNase I is used as a standard.
RNase purity testing	RNase-free pipette tips show no evidence of RNase activity in the fluorometric assay with the detection level of < 3,125*10 ⁻⁹ U/μl when RNase A is used as a standard.
Standards	ISO 9001 Quality management standard ISO 14001 Environmental standard ISO 13485 Medical device quality management standard ISO 14644-1 Class 8 cleanroom manufacturing standard
Autoclavability	No
Transportation conditions	Temperature: -50°C to +50°C Humidity: 10%-90% Shocks: <10 G
Storage conditions	Temperature: -20°C to +40°C Humidity: 10%-60%
Filter pore size	20-40 μm
Expiry date	5 years from manufacturing date*

* Extended 10 μl tip: endotoxin purity testing result < 0,05 IU/ml (EU/ml), expiry date 3 years from manufacturing date.

Materials and Recyclability

Tip	Polypropylene (PP) Recyclable, can be incinerated as energy waste
Filter	Polyethylene (PE) Recyclable, can be incinerated as energy waste
Tray and Rack	Polypropylene (PP) Recyclable, can be incinerated as energy waste
Cardboard package	Cardboard Recyclable, renewable waste

Package Options



Single Tray racks
purity-certified tip racks

Ordering Information

SafetySpace Filter Tips

Order Code	Volume Range [µl]	Length [mm]	Package	Purity Level	Tips/Unit
				Certified free of RNase, Pre-sterilized DNase, endotoxins	
790011F	● 0,1-10	31,5	Single Tray	● ●	10 x 96
783201*	● 0,1-10 Extended	46	Single Tray	● ●	10 x 96
790021F	● 0,5-20	51	Single Tray	● ●	10 x 96
790101F	● 2-120	51	Single Tray	● ●	10 x 96
790201F	● 5-200	52,5	Single Tray	● ●	10 x 96
790301F	● 5-300	52,5	Single Tray	● ●	10 x 96
791001F	● 50-1000	78	Single Tray	● ●	10 x 96
791211F	● 50-1200	90	Single Tray	● ●	10 x 96

* Filter tip with regular air gap



Contact Details

Sartorius Biohit Liquid Handling Oy

Laippatie 1

00880 Helsinki, Finland

Phone +358.9.755.951

E-mail: lhinfo.finland@sartorius.com

Sartorius Lab Instruments GmbH & Co. KG

Weender Landstrasse 94-108

37075 Goettingen, Germany

Phone +49.551.308.0

Fax +49.551.308.3289

www.sartorius.com

The status of the information, specifications and illustrations in this catalogue is indicated by the date given below.

Sartorius reserves the right to make changes to the technology, features, specifications and design of the equipment without notice.

All trademarks are Sartorius property unless otherwise stated. Patents granted or pending.



◀▶ www.sartorius.com