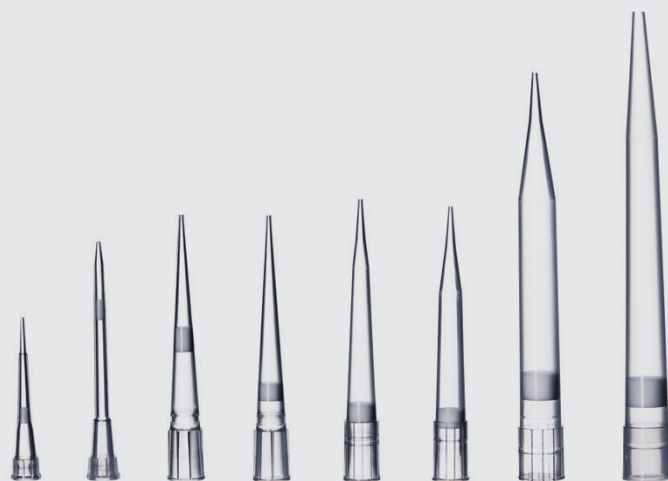


# Safetyspace<sup>®</sup> Filter Tips

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<sup>®</sup> 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<sup>®</sup> Filter Tips are ideal for use with Sartorius pipettes, enabling maximum accuracy and precision.

## Description

Safetyspace<sup>®</sup> 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. Safetyspace<sup>®</sup> Filter Tips are certified and tested for DNase, RNase, human DNA and endotoxins. The tips are released for sale once they have fulfilled the test requirements. Certificates are available at [www.sartorius.com](http://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 5,000 µL
- Tips are tested and certified to be free of DNase, RNase, human DNA and endotoxins
- 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

The Sartorius tip offering also includes Optifit non-filtered tips (see catalog or datasheet 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.

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	Tips are tested and certified to be free of DNase, RNase, human DNA and endotoxins
Endotoxin purity testing	Validated test result with Chromogenic kinetic method according to Ph.Eur. 2.6.14, Method D 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 $< 6.25 \cdot 10^{-5}$ U/ $\mu$ L 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 $< 1 \cdot 10^{-8}$ U/ $\mu$ L.
Human DNA purity testing	Human DNA free pipette tips show no evidence of human DNA in quantitative PCR (qPCR) using human DNA specific primers and with a detection level of $< 1$ pg/ $\mu$ L.
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: $-50^{\circ}\text{C}$ to $+50^{\circ}\text{C}$ Humidity: 10% - 90% Shocks: $< 10$ G
Storage conditions	Storage in room temperature is recommended Temperature: $-20^{\circ}\text{C}$ to $+40^{\circ}\text{C}$ Humidity: 10% - 60%
Filter pore size	20 - 40 $\mu\text{m}$
Expiry date	5 years from manufacturing date*

\* Extended filter tip sterility assurance level, DNase, RNase, human DNA and endotoxin testing detection limits may differ from values reported here. You can find batch specific certificates at [www.sartorius.com](http://www.sartorius.com)

## Materials and Recyclability

Tip	Polypropylene (PP) Recyclable, can be incinerated as energy waste
Polyethylene (PE)	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 DNase, RNase, human DNA, endotoxins	Pre-sterilized	
790011F	■ 10	31.5	Single Tray	■	■	10 x 96
783201*	■ 10 Extended	46	Single Tray	■	■	10 x 96
790021F	■ 20	51	Single Tray	■	■	10 x 96
790101F	■ 120	51	Single Tray	■	■	10 x 96
790201F	■ 200	52.5	Single Tray	■	■	10 x 96
LH-XF780201*	■ 200 Extended	77.5	Single Tray	■	■	10 x 96
790301F	■ 300	52.5	Single Tray	■	■	10 x 96
791001F	■ 1,000	78	Single Tray	■	■	10 x 96
LH-XF791001	■ 1,000 Extended	119	Single Tray	■	■	10 x 96
791211F	■ 1,200	90	Single Tray	■	■	10 x 96
LH-795001F	■ 5,000	150	Single Tray	■	■	1 x 50

\* Filter tip with regular air gap

### 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

### Finland & Baltics

Sartorius Liquid Handling Oy  
Tulppatie 1  
00880 Helsinki  
linfo.finland@sartorius.com

 For further information, visit  
[www.sartorius.com](http://www.sartorius.com)