SARTURIUS

Safetyspace® Filter Tips Available as 5,000 µL

- Bring 5,000 μL filter tips to your workspace in a pre-sterilized purity certified single tray rack.
- DNase, RNase, human DNA and endotoxin free.
- Pair the tip with Picus® Nxt, Picus® or Tacta® for perfect fit, superior ergonomics and maximum accuracy and precision.



The Safetyspace® feature prevents liquid contacting the filter even in reverse pipetting. The filter helps prevent aerosol and liquid contamination and cross-contamination, and reduces the need for pipette maintenance.

Ideal for:

- Cell culture
- Environmental laboratories
- Microbiology

Applications:

- Pipetting foaming liquids such as buffers and proteins
- Pipetting contaminating liquids like microbial cultures
- Pipetting solvents
- Multiple dispensing functions of electronic pipettes
- Reverse pipetting

Features:

- Packed in color-coded single tray racks
- e-beam pre-sterilized
- Certified DNase, RNase, human DNA and endotoxin free
- Expiry 5 years from date of manufacturing
- Manufactured according to ISO 9001, ISO 14001 and ISO 13485
- Manufactured in an ISO Class 8 cleanroom environment



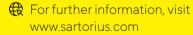




Filter Tips	Product Code	Volume Range	Tip Length	Packaging	Certified		Tips Per Unit
Safetyspace® filter tip 5,000 μL	LH-795001F	100 - 5,000 μL	150 mm	Single tray rack	DNase, RNase, human DNA and endotoxin free	Pre-sterilized	1x50
Standard Tips							
Optifit tip 5,000 μL	780304	100 - 5,000 μL	150 mm	Single tray rack	DNase, RNase, human DNA and endotoxin free	Non-sterile	1x50
	780305	100 - 5,000 μL	150 mm	Single tray rack	DNase, RNase, human DNA and endotoxin free	Pre-sterilized	1x50
	780300	100 - 5,000 μL	150 mm	Bulk in bag		Non-sterile	1x100
	780308	100 - 5,000 μL	150 mm	Bulk in bag		Non-sterile	5x200

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
Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki
Phone +358.9.755.951
Fax +358.9.755.95.292
Ihinfo.finland@sartorius.com