

## SafetySpace Filter Tips Available as 5,000 $\mu$ l



- Bring 5,000  $\mu$ l filter tips to your workspace in a pre-sterilized purity certified single tray rack.
- No more need to compromise sample purity with bulk tips.
- Sartorius Certified: DNase, RNase, Endotoxin free.
- Pair the tip with Picus<sup>®</sup> NXT, Picus<sup>®</sup> or Tacta<sup>®</sup> 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 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



	Product Code	Volume Range	Tip Length	Packaging	Certified		Tips Per Unit
<b>Filter Tips</b>							
SafetySpace filter tip 5,000 µl	LH-795001F	100-5,000 µl	150 mm	Single tray rack	DNase, RNase, endotoxin free	Pre-sterilized	1 x 50
<b>Standard Tips</b>							
Optifit tip 5,000 µl	780304	100-5,000 µl	150 mm	Single tray rack	DNase, RNase, endotoxin free	Non-sterile	1 x 50
	780305	100-5,000 µl	150 mm	Single tray rack	DNase, RNase, endotoxin free	Pre-sterilized	1 x 50
	780300	100-5,000 µl	150 mm	Bulk in bag		Non-sterile	1 x 100
	780308	100-5,000 µl	150 mm	Bulk in bag		Non-sterile	5 x 200

For further information contact your Sartorius local representative.

Sartorius Lab Instruments GmbH & Co. KG  
 Otto-Brenner-Strasse 2037079  
 Goettingen, Germany  
 Phone +49.551.308.0  
 www.sartorius.com

USA Toll-free +1.800.635.2906  
 UK +44.1372.737159  
 France +33.1.70.62.50.00  
 Italy +39.0362.5557.11  
 Spain +34.913.586.095  
 Russian Federation +7.812.327.53.27  
 Japan +81.3.6478.5200

Sartorius Biohit Liquid Handling Oy  
 Laippatie 1  
 00880 Helsinki  
 Finland  
 Phone +358.9.755.951  
 linfo.finland1@sartorius.com