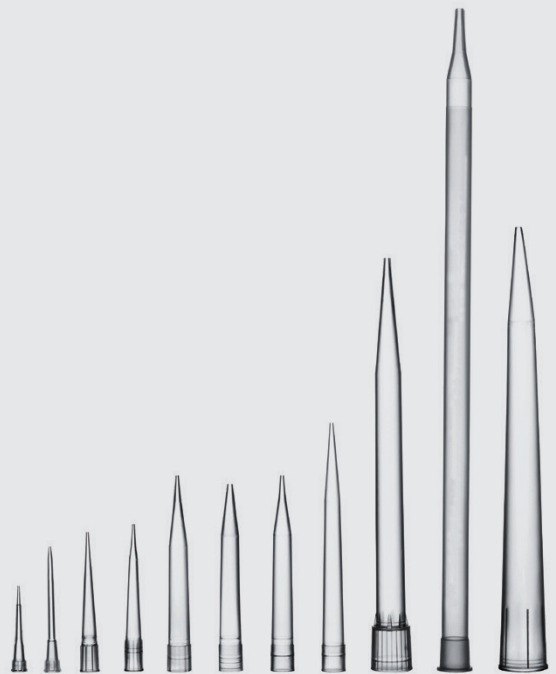


Optifit Tips

Standard Multipurpose
Tips for Various Applications



Product Information

Sartorius Optifit Tips are ideal for use with for Sartorius pipettes, enabling maximum accuracy and precision, as well as comfortable ergonomics. Sartorius pipettes feature a unique spring-loaded Optiload mechanism that makes it easy to attach and eject tips while still ensuring a perfect seal. Non-filter Optifit Tips are available in various packaging types, purity certified and pre-sterilized.

Description

Optifit Tips are manufactured from virgin polypropylene. Strict quality standards (ISO 9001 and ISO 13485) are followed at every step of the process, from R&D to production and delivery. The tips are manufactured in ISO 8 classified cleanroom conditions to avoid any contamination and are certified and tested for DNase, RNase, human DNA and endotoxins. Presterilized tips are also available for laboratory work where purity requirements are of critical importance. The tips are released for sale once they have fulfilled the test requirements.

Certificates are available at www.sartorius.com.

Features

- Available in volumes from 10 µl to 10 ml including extended and wide-bore tips
- Products in single tray, refill pack and FlexiBulk® packages are certified as DNase, RNase, human DNA and endotoxin free
- Informative rack labels indicating tip volume, product number, lot number, expiration date, and purity level for improved tip identification and traceability
- E-beam pre-sterilized tips available for sterile work
- Perfect fitting and tip sealing with all Sartorius pipettes
- Smooth and ergonomic tip attachment and ejection ensured by compatibility with the Optiload feature of Sartorius pipettes
- Color-coded tip trays make it easy to match tip with appropriate Sartorius pipette
- 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

- Tip racks can be refilled with tips from Bulk or Refill packages. The racks and tips are fully autoclavable (121°C, 20 min, 1 bar).

The Sartorius tip offering also includes Safetyspace™ Filter Tips (see catalog or datasheet for details).

Intended Use

The Sartorius Optifit 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 Optifit 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*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 < 1*10 ⁻⁸ U/μL.
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/μL.
Standards	ISO 9001 Quality management standard ISO 14001 Environmental standard ISO 13485 Medical device quality management standard ISO 14644-1 Class 8 cleanroom
Autoclavability	Tips, trays, and racks are autoclavable (121°C, 20 min, 1 bar)
Transportation conditions	Temperature: -50°C to +50°C Humidity: 10% - 90% Shocks: < 10 G
Storage conditions	Storage in room temperature is recommended Temperature: -20°C to +40°C Humidity: 10% - 60%
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

Materials and Recyclability

Tip	Polypropylene (PP) Recyclable, can be incinerated as energy waste
Tray and Rack	Polypropylene (PP) Recyclable, can be incinerated as energy waste
FlexiBulk® pack	Polyethylene terephthalate (APET) / Polyethylene (PE) Recyclable, can be incinerated as energy waste
Cardboard package	Cardboard Recyclable, renewable waste

Package Options



FlexiBulk® Pack



Single Tray



Refill Pack



Bulk packaging



Refill Tower

Ordering Information

Optifit Tips

Order Code	Volume Range	Length	Packaging	Purity Level		Tips/Unit
				Free of DNase, RNase, human DNA & endotoxins	Pre-sterilized	
790010	■ 10 µL	31.5 mm	Single Tray	■		10 × 96
790011			Single Tray	■	■	10 × 96
790012			Refill Tower			10 × 96
790013			Refill Pack	■	■	20 × 96
790014			Bulk in Bag			1,000
790200	■ 200 µL	51 mm	Single Tray	■		10 × 96
790201			Single Tray	■	■	10 × 96
790202			Refill Tower			10 × 96
790203			Refill Pack	■	■	15 × 96
LH-B790204			FlexiBulk®	■		960
790350	■ 350 µL	54 mm	Single Tray	■		10 × 96
790351			Single Tray	■	■	10 × 96
790352			Refill Tower			10 × 96
790353			Refill Pack	■	■	15 × 96
LH-B790354			FlexiBulk®	■		960
791000	■ 1,000 µL	71.5 mm	Single Tray	■		10 × 96
791001			Single Tray	■	■	10 × 96
791002			Refill Pack	■		10 × 96
791003			Refill Pack	■	■	10 × 96
LH-B791004			FlexiBulk®	■		480
791020	■ 1,000 µL Wide bore tip	68.5 mm	Single Tray	■		10 × 96
791021			Single Tray	■	■	10 × 96
LH-B791024			FlexiBulk®	■		480
791200	■ 1,200 µL	71.5 mm	Single Tray	■		10 × 96
791201			Single Tray	■	■	10 × 96
791202			Refill Pack	■		10 × 96
791203			Refill Pack	■	■	10 × 96
LH-B791204			FlexiBulk®	■		480
780304	■ 5,000 µL	150 mm	Single Tray	■		50
780305			Single Tray	■	■	50
780300			Bulk in Bag			100
780308			Bulk in Bag			1,000
LH-780314	■ 10,000 µL	155 mm	Single Tray			35
LH-780316			Bulk in Bag			250

Extended Standard Tips

Order Code	Volume Range	Length	Packaging	Purity Level		Tips/Unit
				Free of DNase, RNase, human DNA & endotoxins	Pre-sterilized	
783210	■ 10 µL	46 mm	Single Tray	■		10×96
783211			Single Tray	■	■	10×96
LH-X780200	■ 200 µL	77.5 mm	Single Tray	■		10×96
LH-X780201			Single Tray	■	■	10×96
LH-X791000	■ 1,000 µL	119 mm	Single Tray	■		10×96
LH-X791001			Single Tray	■	■	10×96
791210	■ 1,200 µL	90 mm	Single Tray	■		10×96
791211			Single Tray	■	■	10×96
791212			Refill Pack	■		10×96
791213			Refill Pack	■	■	10×96

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