

1,000 μ L Extended Plus Tips

Reach Further Into Long and Narrow Vessels to Collect Your Sample
With the New 1,000 μ L Extended Plus Tips

Highlights

- Longest 1,000 μ L tip on the market
- Narrow and extended tip form to reach easier to the bottom of deep, narrow vessels
- Comes as standard or pre-sterilized either with or without Safetyspace[®] filter

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



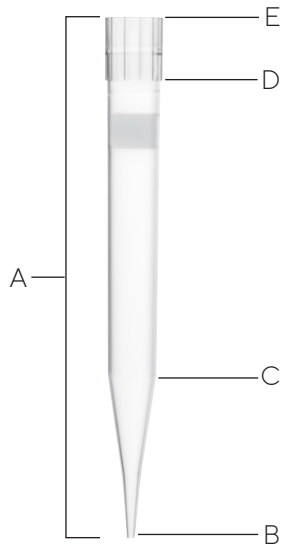
Meet the Newest Addition to the Sartorius Tip Portfolio, the New and Improved 1,000 μ L Extended Plus Tips

The Extended Plus tips has even further extended reach compared to other 1,000 μ L Extended tips available on the market, which allows you to reach those last drops in even narrower, deeper vessels.

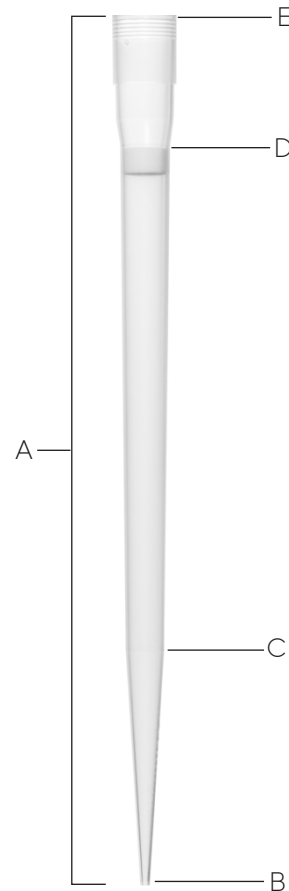
The 1,000 μ L Extended Plus tips are available as non-sterile and pre-sterilized with or without a Safetyspace[®] filter.

The Extended Plus tips have a universal design and optimized to be used in combination with Sartorius pipettes. With the extra length, the tip minimizes the risk of contamination, as the pipette does not need to touch the vessel wall. The long and narrow form also allows easier sample collection from the bottom of the vessel without tilting.

Dimensions	Optifit Safetyspace [®] 1,000 μ L	Optifit Safetyspace [®] Extended Plus 1,000 μ L
Length		
A	71.4 78.2 mm	119 mm
External Diameter		
B	1.3 mm; Internal diameter: 0.8 mm	1.3 mm; Internal diameter: 0.8 mm
C	6.9 7.4 mm	5.5 mm
D	8.0 8.1 mm	7.0 mm
E	8.7 8.9 mm	8.8 mm



Safetyspace[®] 1,000 μ L tip



Safetyspace[®] Extended Plus 1,000 μ L tip

Application

With the 1,000 µL Extended Plus tip, it is possible to reach to the bottom of long and narrow vessels such as 15 mL conical tubes, without contaminating the pipette.

Extended Plus tips are excellent when working with deep well plates (adding reagents and transferring samples), collecting samples from long (up to 12 cm) and narrow vials or tubes, collecting retention samples from Vivaspin® tubes (ultrafiltration devices), working with blood and serum samples (contamination risk of the pipette is minimized), needing to reach deep into tubes to collect supernatant without disturbing cell pellets or other pelleted samples, inoculating bacterial or yeast cultures into deep culture tubes or flasks.



Purity

The new Extended Plus tip comes in different purity variants, i.e., RNase, DNase, human DNA and Endotoxin free or pre-sterilized either with or without SafetySpace® filter.

Material	RNase, DNase, human DNA & Endotoxin Free	Pre-sterilized	SafetySpace® Filter
LH-X791000	X		
LH-X791001	X	X	
LH-XF791001	X	X	X

Definitions	Description	Purity Level
Human DNA	Sartorius purity certified tips are analyzed for the presence of DNA using quantitative PCR and human DNA specific primers.	< 1 pg/µL
DNase	Deoxyribonuclease (DNase) is any enzyme that catalyzes the degradation of DNA. The absence of DNase is tested by using a fluorometric assay.	< 6.25 × 10 ⁻⁵ U/µL
RNase	Ribonuclease (RNase) is an enzyme that catalyzes the degradation of RNA into smaller components and can be generally found from organisms. The absence of RNase is tested by using fluorometric assay.	< 1 × 10 ⁻⁶ U/µL
Endotoxin	Endotoxins are lipopolysaccharides found in Gram-negative bacteria, which can cause serious health effects in humans and animals. Limulus Amebocyte Lysate (LAL)* is used to detect the presence of endotoxins on the pipette tips.	< 0.005 IU/mL (EU/mL)
Sterility	We use e-beam irradiation to pre-sterilize the pipette tips.	—

*Our tips are currently being tested with a LAL assay, applying the EP. 2.6.14 Chromogenic kinetic method D

Germany

Sartorius Lab Instruments GmbH & Co. KG
 Otto-Brenner-Strasse 20
 37079 Göttingen
 Phone +49 551 308 0

USA

Sartorius Corporation
 3874 Research Park Drive
 Ann Arbor, MI 48108
 Phone +1 734 769 1600

 For further information, visit
sartorius.com