

Questions to Ask When Choosing Your Pipette Tips

August 2023 | Content supplied by Maria Voutilainen, PhD. and Sandra Söderholm, PhD. Published by Lab Manager

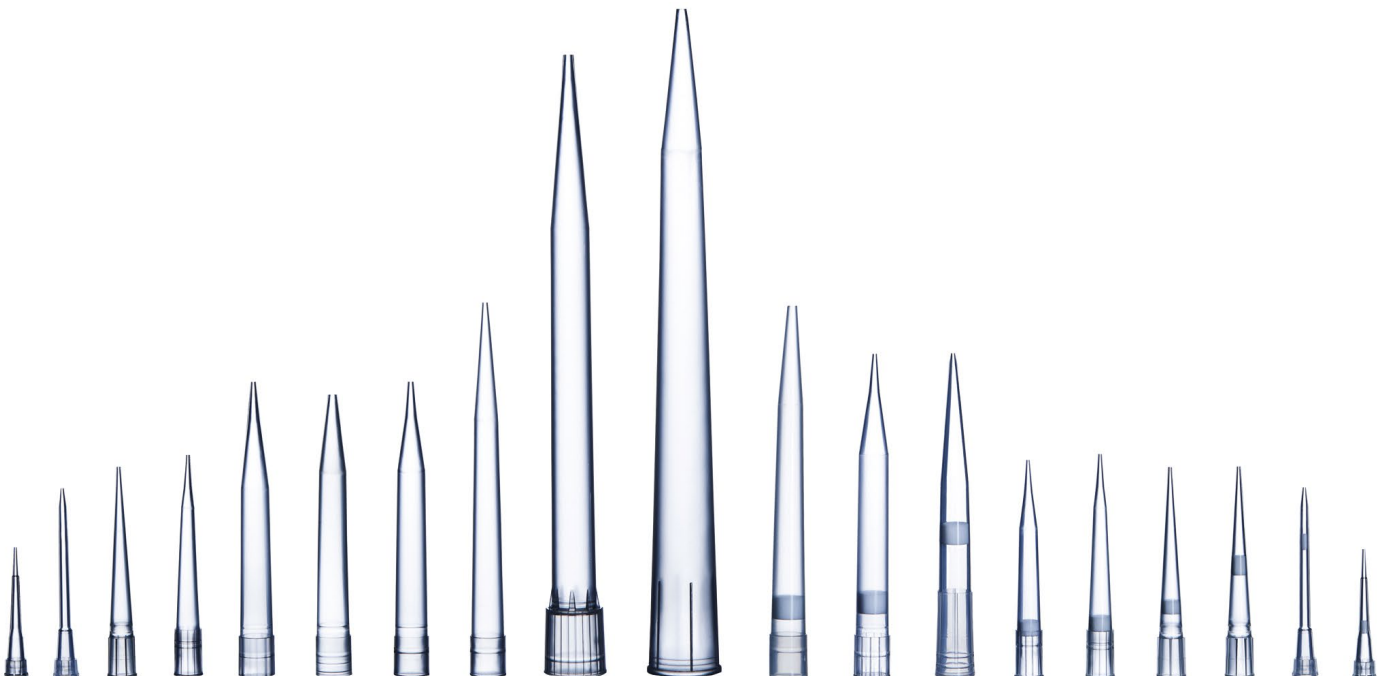


Careful selection of tips ensures accurate and reliable pipetting results and successful experiments.



Pipettes and pipette tips are used in laboratories for various different liquid transfer tasks. Among the many common laboratory consumables, high-quality pipette tips are crucial for successful experiments in the laboratory. The best results are achieved when the pipette and tip match perfectly with each other and the tip variant is optimal for the liquid being transferred and for the demands of the application. Though the cost of the tips is important, the price is not the only factor that should be taken into consideration when choosing your tips.

Proper evaluation and consideration of pipette tips on criteria such as compatibility, accuracy, and precision with your pipette, purity, packaging, ergonomics and ease of use, and the overall quality of the tips, can go a long way to streamlining and optimizing laboratory workflows and experiments. The use of high-quality tips that match your pipette may end up being a more cost-effective solution as the pipetting results are more reliable and fewer repetitions – and eventually fewer tips – might be required.





About Maria Voutilainen

Maria Voutilainen is a Product Manager at Sartorius. In her current position, she is responsible for managing liquid handling consumables in Sartorius' Laboratory Products and Services portfolio. Maria has an extensive experience in working with various laboratory instruments and consumables across different phases of the product life cycle. Maria holds a Ph.D. in genetics from the University of Helsinki, Finland.

Maria Voutilainen
Product Manager

6 Questions to Ask When Purchasing Pipette Tips

1. Are the pipette tips compatible with my pipette model?
2. How ergonomic are my pipette and tips as a system?
3. What are the certificates of tip purity I need for my workflow (e.g., DNase, RNase, and endotoxin-free)?
4. Does my application require the use of pre-sterilized and/or filtered tips?
5. Does my application require the use of specialty tips such as extended or wide bore tips?
6. What packaging options are available for the tips?

Sartorius Standard Optifit and Safetyspace® Filter Tips

Sartorius pipettes and tips are designed to be a seamlessly compatible system that ensures reliable and repeatable pipetting results. The high-quality standard Optifit tips are a preferred choice for a wide range of laboratory applications and pipetting tasks, while the Safetyspace® Filter tips ensure contamination-free work by protecting both your samples and the pipette. The unique Safetyspace® feature leaves extra room between the filter and your sample allowing the use of the full volume of the tip with any pipetting mode. The broad selection of packaging and purity options for the tips offers suitable alternatives for the specific demands of different laboratory settings. For optimal accuracy and precision, it is important to calibrate your pipette using the pipette tip variants (e.g., standard, pre-sterilized, filter tip) utilized in your work.

Bulk and Refill Tips – An Affordable and Less Waste Generating Option

Laboratories with high consumption rates of pipette tips often lean toward purchasing bulk and/or refill tips as a more affordable alternative. Bulk and refill tips reduce waste generation by allowing empty tip boxes to be reused rather than being discarded. In selecting your tips, it is important to acknowledge the differences that can exist between manufacturers. Things to consider include:

1. Ensuring that the tips can withstand autoclaving without deformation;
2. Checking the compatibility and perfect fitting of the tip to the pipette;
3. Confirming that the tips are made from high-quality virgin polypropylene with molding technology that ensures a smooth and consistent tip surface;
4. Choosing tips packed in durable containers and are easy to load onto racks.

To learn more, visit:


www.sartorius.com/en/products/pipetting/pipette-tips

Germany

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

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 650 322 1360

 For further information, visit
www.sartorius.com