

SARTORIUS

Pipette Essentials

Make Your Choice



Check These Features to Choose Your Perfect Pipette

- **Reliable Accuracy and Precision**

Check that the pipette manufacturer complies with international standards and regulations (e.g., ISO).

- **Ergonomic Pipette Handle**

Test that the pipette feels comfortable in your hand and is convenient to operate.

Check that the finger hook suits your hand without any need to squeeze or strain your muscles.

- **Smooth and Light Tip Ejection**

Test that the tip ejection function is light and ergonomic to use.

- **Light Weight**

Test the weight of the pipette. Light weight is essential in long pipetting series to prevent muscle strain.

- **Easy to Maintain**

Check that the pipette is easy to disassemble without tools and has only a few parts to clean.

Mechanical Pipettes

- **Light Operation of the Pipette**

Check that the pipette's plunger is light to press at both maximum and minimum volumes.

- **Easy-to-Read Display**

Check that the volume is displayed in numbers, down to the last decimal.

- **Display Visible While Working**

Check that the volume is easy to read while working.

- **Volume-Lock**

Check that the pipette has a volume-lock to avoid accidental volume changes.

- **Fully Autoclavable**

Check that the pipette is fully autoclavable. Autoclaving is the most efficient way to decontaminate a pipette.

Electronic Pipettes

- **Electronic Tip Ejection**

Check that the pipette has electronic tip ejection.

- **Plug & Play Operation**

Check that the pipette has the basic pipetting functions available for use intuitively (pipetting, volume and speed adjustment).

- **Advanced Features**

Check that the electronic pipette supports the performing of complex workflows, (e.g., plate tracker function for microwell plate assays).

If used in regulated laboratories, check that the electronic pipette has the following features: calibration reminders, possibility to prevent unauthorized changes in important settings.

Also investigate whether you could benefit from connectivity features, like performing guided workflows with the help of an app.



Test the Pipette in Your Lab —Ask for a Demo Pipette from Sartorius!

www.sartorius.com

Sartorius Liquid Handling Oy
Tulppatie 1
00880 Helsinki, Finland
lhinfo.finland1@sartorius.com

Specifications subject to change without notice. Copyright Sartorius Lab Instruments GmbH & Co. KG. All Rights Reserved. Printed in the EU or US on paper bleached without chlorine. Version 2 | 02 | 2025