SARTURIUS

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 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: locking of protocols, user profiles, calibration notifications.



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

www.sartorius.com

Sartorius Biohit Liquid Handling Oy Laippatie 1 00880 Helsinki, Finland Ihinfo.finland@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 1 | 06 | 2021