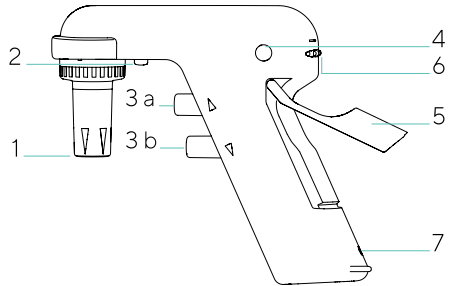


Midi Plus Electronic
Pipetting Controller



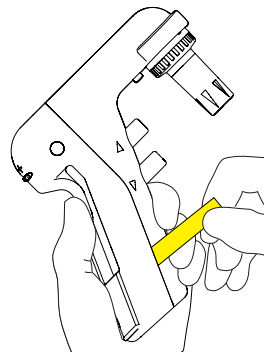
Features

1. Nose-cone, silicone adapter and hydrophobic filter
2. Locking button
3. Operating buttons
 - a. Aspiration
 - b. Dispensing
4. Handle cover release buttons
5. Fold-out stand
6. Speed adjustment wheel
7. LED battery indicator

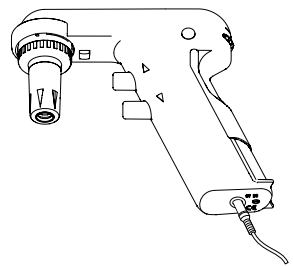


Getting Started

1. Midi Plus is delivered with the battery inserted. The battery poles are covered by a plastic protective film. Before first use, remove the plastic film by pulling it gently from below the battery cover.
2. Charge Midi Plus for at least 4 hours before first use.
 - To charge with AC-adaptor: connect Midi Plus to the charger cable and plug the charger into the mains power outlet.
3. The battery status is indicated by a LED. When the battery is fully charged the LED emits a green light, and the Midi Plus is ready for use.



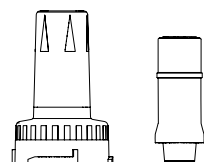
Midi Plus is equipped with an adapter set (nose-cone and silicone adapter) that fits all commonly used serological or volumetric pipettes.



Pipetting is possible while connected to the AC-adaptor

Operation

1. Make sure the hydrophobic filter is inserted
2. Inserting a pipette:
 - Gently push a pipette into the silicone adapter holding the pipette at the top. Make sure the pipette is attached firmly and is air-tight.
3. Adjust the speed by turning the speed adjustment wheel



Standard nose-cone and silicone adapter set

4. Aspiration:

- Immerse the pipette into the liquid
- Press the aspirating button until the pipette is filled just above the desired volume
- Lift the pipette from the liquid
- Press the dispensing button to slowly dispense the liquid until the level is exactly at the desired volume

5. Dispensing:

- Place the pipette against the receiving vessel wall
- Press the dispensing button until the pipette is empty
- To activate the blow-out, press the dispensing button completely

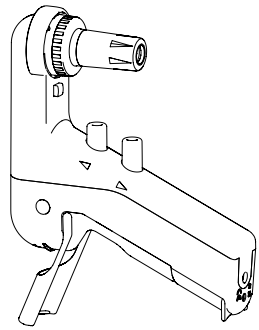
Note: Press the dispensing button gently to let the liquid flow out by the force of gravity. Fine-tune the speed by adjusting finger pressure while dispensing.

Note: To ease the aspiration of certain liquids, such as volatile or viscous ones, increase the speed, especially when using a large volumetric pipette.

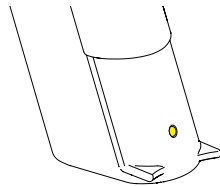
Other Recommendations

1. When not in use, Midi Plus should always be placed on a table to rest on the fold-out stand.
2. The LED indicator starts to blink when the battery charge level is low: connect the Midi Plus to the power supply.
 - Pipetting is possible while connected to the AC-adaptor.
3. Midi Plus can be used with Sartorius Optifit 5,000 μl and 10,000 μl tips. Adapters for the 5,000 μl and 10,000 μl Optifit Tips are available as accessories.

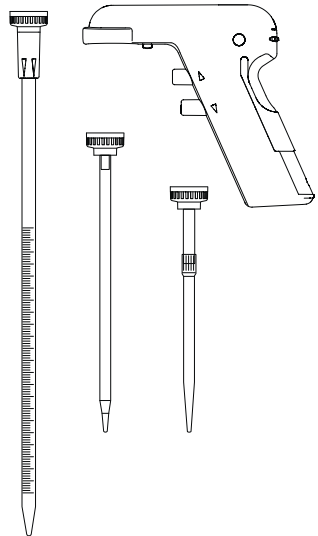
Midi Plus, with a variety of optional nose-cones and silicone adapters. Optional extras include the Optifit 5,000 μl and 10,000 μl tips, corresponding nose-cones and silicone adapters, and are available separately.



Fold-out stand



LED indicator starts to blink when the battery charge level is low



Midi Plus and Accessories

Order Code	Item	Qty Unit
710931	Midi Plus	1
LH-7129120	Replacement filter, 0,45 µm, non-sterile, autoclavable	5
LH-7129131	Replacement filter, 0,2 µm, sterile	5
LH-711019	Adapter set (standard) including nose-cone and the silicone adapter	1
LH-711017	Adapter set for 5,000 µl pipette tip, autoclavable, including nose-cone and the silicone adapter	1
LH-711018	Adapter set for 10,000 µl pipette tip, autoclavable, including nose-cone and the silicone adapter	1
LH-712146	Nose-cone (standard)	1
LH-712144	Nose-cone for 5,000 µl Optifit Tip	1
LH-712145	Nose-cone for 10,000 µl Optifit Tip	1
LH-712153	Silicone adapter (standard)	1
LH-712151	Silicone adapter for 5,000 µl Optifit Tip	1
LH-712152	Silicone adapter for 10,000 µl Optifit Tip	1
780300	Optifit Tip 5,000 µl	100
780308	Optifit Tip 5,000 µl	1,000
780310	Optifit Tip 10,000 µl (Extra long tip for Midi Plus)	100

Note: Only use the power supply provided with the Midi Plus.

This Quick Start Guide is for the basic operation of the Midi Plus pipette controller. For further information and more detailed instructions, please see the user manual at www.sartorius.com.

A paper print-out is available on request from lhinfo.finland@sartorius.com.

Sales and Service Contacts

For further contacts, visit www.sartorius.com



Finland

Sartorius
Biohit Liquid Handling
Laippatie 1
00880 Helsinki
Phone +358 9 755 951

Germany

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

USA

Sartorius Corporation
5 Orville Drive, Suite 200
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

UK importer

Sartorius UK Ltd.
Longmead Business Centre,
Blenheim Road
KT19 9 QQ Epsom, Surrey