

Midi Plus

Ergonomic Large-Volume Pipetting

Simplifying Progress

SVISCISVS

Easy and Ergonomic Pipetting of Large Volumes from Bottles or Tubes

Ergonomic design protects you from RSI

Midi Plus is a light-weight electronic pipette controller, which allows easy aspiration from bottles and tubes. Its ergonomic design helps to minimize arm and hand raises that are often required while using serological or volumetric pipettes. In addition the electronic operation of Midi Plus reduces the needed finger force compared to a pipette filler with a squeezable bulb. These qualities ease the workload in long periods of pipetting and protect you from Repetitive Strain Injury (RSI).

Compatible with a wide range of pipettes and tips

- The Midi Plus is compatible with commonly used glass or plastic pipettes in the 1 100 ml volume range.
- It is also compatible with Sartorius 5 ml and 10 ml Optifit Tips with adapter sets, available separately.

Hydrophobic filters prevent over-aspiration

- The hydrophobic Sartorius filter prevents over-aspiration, which saves the pipetting controller from contamination.
- Filters are available as sterile and non-sterile, to match your purity requirements.

Save bench space by using the fold-out stand

- The integrated fold-out stand saves bench space.
- While Midi Plus is resting on the stand, the glass or plastic pipette is kept detached from the table surface, which prevents contamination.





Stepless speed control for accurate dispensing

- Aspiration and dispensing speeds can easily be controlled by using the adjustment wheel or by adjusting finger pressure on the operating buttons.
- The speed can be selected according to your application and pipette volume.
- Possibility to dispense by the force of gravity for slow, drop-by-drop dispensing. Just by applying a very small force on the dispensing button.

Ordering Information 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 μ m, pre-sterilized	5
LH-711019	Adapter set (standard) including nose-cone and the silicone adapter	1
LH-711017	Adapter set for 5 ml pipette tip, autoclavable, including nose-cone and the silicone adapter	1
LH-711018	Adapter set for 10 ml pipette tip, autoclavable, including nose-cone and the silicone adapter	1
780300	Optifit Tip 5 ml	100
780308	Optifit Tip 5ml	1000
780310	Optifit Tip 10 ml (extra long tip for Midi Plus)	100



Germany

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

USA

Sartorius Corporation 565 Johnson Avenue Bohemia, NY 11716 Phone +1 631 254 4249 Toll-free +1 800 635 2906

For further contacts, visit www.sartorius.com