



Midi Plus

Pipetting Controller

Product Information

Midi Plus is a light-weight electronic pipette controller, which allows easy liquid aspiration from bottles and large tubes. Its ergonomic design helps to minimise arm and hand raises that are often required while using serological or volumetric pipettes. This eases the workload in long periods of pipetting, and helps to protect the user from Work Related Upper Limb Disorder (WRULD).



Description

Midi Plus is developed for easy and safe large-volume dispensing. It provides an ergonomic alternative to the simple pipette filler with a squeezable bulb, which requires significant finger force when aspirating and dispensing liquids. The electronic operation of Midi Plus reduces the needed force, and the stepless speed control improves the control of the liquid flow for better accuracy. Dispensing by using the force of gravity is possible, by gently pressing the operating button. Pipetting accuracy depends on the user's ability to set the meniscus of the liquid at the graduation marks on a serological pipette.

Midi Plus features convenient cordless pipeting. The battery can be charged quickly and has a long life-time. When the battery is low, it can be charged while working.

Midi Plus can be used with commonly used glass or plastic pipettes in the 1 to 100 ml volume range, and also with 5 ml and 10 ml Sartorius Optifit Tips.

Features

- Ergonomic handle design with electronic operating buttons for reduced finger force
- Stepless speed control for accurate dispensing
- Possibility of dispensing with the force of gravity for slow, drop-by-drop dispensing, eg. into culture media dish not to break the

fine surface of the medium

- Fold-out stand saves bench space when Midi Plus is not in use and keeps the pipette free from the table's surface to avoid contamination.
- Wide compatibility with glass or plastic pipettes, also compatible with Sartorius 5 ml Optifit Tips and 10 ml extra-long Optifit Tips (adapter sets for 5 ml and 10 ml tips available separately)
- Hydrophobic 0,45 µm and 0,2 µm (sterile) filters help to prevent over-aspiration, and thus contamination of the device
- Direct charging cable enables operation while charging
- Low battery indicator

Applications

- Cell culture
- Liquid preparation in large volumes

Intended Use

For large volume pipetting from bottles or large tubes in volume range of 1 to 100 ml.

Technical Data

Technical Specifications	
Weight	212 g
Autoclavability	Autoclavable nose cone and silicone adapters (121°C, 20 min, 1 bar/100 kPa)
Battery (included in package)	Rechargeable NiMH battery (set of three batteries)
Charger (included in package)	Universal charger with EU, US/ JPN, UK, CHN, KOR and AUS plugs
Charging time	Approx. 3 hours
Battery charge	~2000* pipetting cycles (when fully charged)
Maximum pipetting speed	~10 ml/s
Dispensing method	Motor driven dispensing Gravity dispensing
Compatibility with volumetric pipettes	1-100 ml common glass or plastic pipettes

Compatibility with tips (adapter set available separately)	5 ml Sartorius Optifit Tips 10 ml Sartorius extra long Optifit Tips
Hydrofobic filters	Pore size 0,45 µl, autoclavable (2 included in package) Pore size 0,2 µl, sterile (available separately)
Recommended operating temperature	Normal room temperature
Transportation conditions	Temperature: -50°C to +50°C Humidity: 10-90% Shocks: <10 G
Long-term storage conditions	Temperature: -5°C to +35°C Humidity: 10-60%
Warranty	2 years

*Depends on dispensing method, used liquid and volumes.

Ordering Information

Order Code	Description	Quantity
710931	Midi Plus Pipette Controller with Universal Charger	1
LH-7129120	Replacement filter, 0,45 µm, non-sterile, autoclavable	5
LH-7129130	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 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, Bulk in Box	100
780308	Optifit Tip 5 ml, Bulk in Box	1000
780310	Optifit Tip 10 ml, Bulk in Box, (extra long tip for Midi Plus)	100

Sartorius Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki, Finland
Phone +358.9.755.951

Sartorius Lab Instruments GmbH & Co. KG
Weender Landstrasse 94-108
37075 Goettingen, Germany
Phone +49.551.3080
Fax +49.551.308.3289



www.sartorius.com