# SARTURIUS

# Midi Plus Pipette Controller

## Easy and Ergonomic Pipetting of Large Volumes From Bottles or Tubes

#### Introduction

Midi Plus is a light-weight electronic pipette controller, which allows easy liquid aspiration from bottles and large tubes. Its ergonomic design helps to minimize arm and hand raises that are often required while using serological or volumetric pipettes. This eases the workload in long periods of pipetting and protects the user from Repetitive Strain Injury (RSI). 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, adapters available separately.



### Description

Midi Plus is developed for easy and safe large-volume dispensing. It provides an ergonomic alternative to the simple pipette filler with squeezable bulb, which requires extensive finger force when aspirating and dispensing liquids. The electronic operation of Midi Plus reduces the finger force needed, and the stepless speed control improves the control of the liquid flow for better accuracy. Dispensing by the force of gravity is possible by applying gentle finger pressure on 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. The improved battery enables quick charging and long battery lifetime. When the battery is low, it can be charged while working. When fully charged, the cord can be unplugged from the device, allowing safe and convenient cordless pipetting.

#### **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 (e.g., into culture media dish so as 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 detached from the table 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 pre-sterilized 0.2 µm filters prevent over-aspiration and 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

## Technical Specifications

Rechargeable Battery		
Туре	NiMH (set of 3 batteries, only for Midi Plus)	
Charging time	Approx. 3 hours	
AC-Adaptors		
Included in package	Universal adaptor with Eur, US   Jpn, UK, Chi and Aus plugs	
Weight	212 g	
Number of Pipetting Cycles	~ 2,000	
Maximum Pipetting Speed	~ 10 mL/s	
Pipette Compatibility		
Adapter set included in package	1-100 mL common glass or plastic pipettes	
Adapter set available separately	5 mL Optifit Tip	
Adapter set available separately	10 mL extra long Optifit Tip	
Gravity Dispensing	Available	
Hydrophobic Filters		
2 included in package	0.45 μm autoclavable	
Available separately	0.2 µm pre-sterilized	
Autoclavability		
Autoclavable nose cone and silicone adapters	121° C, 20 min, 1 bar	

## Ordering Information

Cat. No.	Description	Quantity
710931	Midi Plus™ Pipette Controller with universal AC-adaptor	1
LH-7129120	Replacement filter, 0.45 μm, non-sterile, autoclavable	5
LH-7129131	Replacement filter, 0.2 μm, 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
LH-712146	Nose cone (standard)	1
LH-712144	Nose cone for 5 mL Optifit Tip	1
LH-712145	Nose cone for 10 mL Optifit Tip	1
LH-712153	Silicone adapter (standard)	1
LH-712151	Silicone adapter for 5 mL Optifit Tip	1
LH-712152	Silicone adapter for 10 mL Optifit Tip	1
780300	Optifit Tip 5 mL, bulk in box	100
780308	Optifit Tip 5 mL, bulk in carton	1,000
780310	Optifit Tip 10 mL, bulk in box, (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 information, visit www.sartorius.com