

# Midi Plus™ Pipette Controller

Easy and ergonomic pipetting of large volumes from bottles or tubes



## Technical Specifications

|                                              |                                                           |
|----------------------------------------------|-----------------------------------------------------------|
| <b>Rechargeable battery</b>                  |                                                           |
| Type                                         | NiMH (set of 3 batteries, only for Midi Plus)             |
| Charging time                                | Approx. 3 hours                                           |
| <b>AC-adaptors</b>                           |                                                           |
| Included in package                          | Universal adaptor with Eur, US/Jpn, UK, Chi and Aus plugs |
| Weight                                       | 212 g                                                     |
| Number of pipetting cycles                   | ~2000                                                     |
| Maximum pipetting speed                      | ~10 ml / s                                                |
| <b>Pipette compatibility</b>                 |                                                           |
| 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                              |
| <b>Gravity Dispensing</b>                    | Available                                                 |
| <b>Hydrophobic Filters</b>                   |                                                           |
| 2 included in package                        | 0.45 µm autoclavable                                      |
| available separately                         | 0.2 µm pre-sterilized                                     |
| <b>Autoclavability</b>                       |                                                           |
| Autoclavable nose cone and silicone adapters | 121°C, 20 min, 1 bar                                      |

## 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.

## 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 MidiPlus™ 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 life-time. 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, 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 detached from the table surface to avoid contamination (fold-out stand not available in North America).
- 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.

## Ordering Information

| Order Code | Description                                                                                   | Quantity |
|------------|-----------------------------------------------------------------------------------------------|----------|
| 710931     | Midi Plus™ Pipette Controller with Universal AC-adaptor                                       | 1        |
| 710929     | Midi Plus™ Pipette Controller (order code for North America)                                  | 1        |
| LH-7129120 | Replacement filter, 0.45 µm, non-sterile, autoclavable                                        | 5        |
| LH-7129130 | 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                                                              | 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  
 Fax +358.9.755.95.220

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](http://www.sartorius.com)



Specifications subject to change without notice.  
 Printed and copyrighted by  
 Sartorius Biohit Liquid Handling Oy  
 Publication No.:SUL2001-e131102  
 Order No.: 85037-542-21  
 Ver. 11|2013