

## mLINE® Pipette Multipacks



# Our Most Ergonomic Mechanical Pipettes in an Affordable Package of Your Choice!

- 5 mLINE® Multipack options available for various needs with a great value
- A set of 3 to 5 mLINE® mechanical pipettes, a Linear Stand and racks of matching tips included
  - Renew your existing pipettes with a set of new, highly ergonomic pipettes with excellent accuracy and precision
  - Set up new workstations for every researcher with new, high quality pipettes

Order Code  
**LH-725661**

Multipack Contents  
**mLINE® Pipette 3-pack 10**

- mLINE® pipettes 0.5-10 µl, 10-100 µl, 100-1000 µl
- Sartorius Optifit Tip tray box 0.2-10 µl, 0.5-200 µl, 10-1000 µl
- Linear Stand



Order Code  
**LH-725662**

Multipack Contents  
**mLINE® Pipette 3-pack 20**

- mLINE® pipettes 2-20 µl, 20-200 µl, 100-1000 µl
- Sartorius Optifit Tip tray box 0.5-200 µl, 10-1000 µl
- Linear Stand



Order Code  
**LH-725663**

Multipack Contents  
**mLINE® Pipette 4-pack**

- mLINE® pipettes 0.5-10 µl, 10-100 µl, 20-200 µl, 100-1000 µl
- Sartorius Optifit Tip tray box 0.2-10 µl, 0.5-200 µl, 10-1000 µl
- Linear Stand



## mLINE® - Effortless Accuracy

Sartorius' most advanced mechanical pipette in the Biohit*family* - the mLINE® - offers excellent ergonomics, performance and safety. It is the best choice for long-lasting and repetitive manual pipetting in order to prevent the risk of work related hand, arm or shoulder disorders.

### Ergonomics

- Exceptionally light pipetting and tip ejection forces protect you from repetitive strain injuries (RSI)
- Ergonomic finger support allows the pipette to rest effortlessly in your hand

### Performance

- Thermal insulation of internal components for steady sampling temperature provides you with highly accurate and repeatable results also in long pipetting series
- Easy to use calibration allows quick re-calibration for liquids with various consistencies assuring better accuracy across a variety of liquids

### Safety

- Minimized risk of contamination with Safe-Cone Filters and full autoclavability (steam sterilization)
- Volume-lock prevents accidental volume changes while pipetting for error free results



Order Code    Multipack Contents  
**LH-725665**    **mLINE® Pipette 3+1-pack**

- mLINE® pipettes 0.5-10 µl, 20-200 µl, 100-1000 µl and 8-channel 30-300 µl
- Sartorius Optifit Tip tray box 0.2-10 µl, 0.5-200 µl, 5-350 µl, 10-1000 µl
- Linear Stand



Order Code    Multipack Contents  
**LH-725664**    **mLINE® Pipette 5-pack**

- mLINE® pipettes 2-20 µl, 10-100 µl, 20-200 µl, 100-1000 µl, 500-5000 µl
- Sartorius Optifit Tip tray box 0.5-200 µl, 10-1000 µl, 5000 µl
- Linear Stand

## Technical Specifications

Item	Channels	Volume	Increment	Test volume	Systematic error	Random error
mLINE® Mechanical Pipette	1	0.5-10 µl	0.01 µl	10 µl	1.00 %	0.60 %
				5 µl	1.50 %	1.00 %
				1 µl	2.50 %	1.50 %
mLINE® Mechanical Pipette	1	2-20 µl	0.02 µl	20 µl	0.90 %	0.40 %
				10 µl	1.20 %	1.00 %
				2 µl	3.00 %	2.00 %
mLINE® Mechanical Pipette	1	10-100 µl	0.10 µl	100 µl	0.80 %	0.15 %
				50 µl	1.00 %	0.40 %
				10 µl	2.00 %	1.00 %
mLINE® Mechanical Pipette	1	20-200 µl	0.20 µl	200 µl	0.60 %	0.15 %
				100 µl	0.80 %	0.30 %
				20 µl	2.00 %	0.80 %
mLINE® Mechanical Pipette	1	100-1000 µl	1.00 µl	1000 µl	0.60 %	0.20 %
				500 µl	0.60 %	0.20 %
				100 µl	1.00 %	0.40 %
mLINE® Mechanical Pipette	1	500-5000 µl	10.0 µl	5000 µl	0.50 %	0.20 %
				2500 µl	0.60 %	0.30 %
				500 µl	2.00 %	0.60 %
mLINE® Mechanical Pipette	8	30-300 µl	0.20 µl	300 µl	0.60 %	0.25 %
				150 µl	1.00 %	0.50 %
				30 µl	2.00 %	1.00 %

The data for systematic error (inaccuracy) and random error (imprecision), according to EN ISO 8655, is valid only when Sartorius pipette tips are used

### Contact Details

**Sartorius Corporation**  
5 Orville Drive, Suite 200  
Bohemia, NY 11716  
Phone | 800.635.2906  
Fax | 631.254.4253  
www.sartorius.us

**Sartorius Canada**  
2179 Dunwin Drive #4  
Mississauga, Canada ON L5L 1X2  
Phone +1.800.668.4234  
Fax +1-905.569.7021  
www.sartorius.com

**Sartorius Biohit Liquid Handling Oy**  
Laippatie 1  
00880 Helsinki  
Phone +358.9.755.951  
Fax +358.9.755.95.292  
lhinfo.finland@sartorius.com

**Sartorius Lab Instruments GmbH & Co. KG**  
Weender Landstrasse 94-108  
37075 Goettingen  
Phone +49.551.308.0  
Fax +49.551.308.3289



◀ www.sartorius.com