

## Mline® Mechanical Pipette Multipacks Convenient Combo-Pack for Your Lab

### Mline® Mechanical Pipette

- Low pipetting forces
- Ergonomic finger support minimizes the grip force needed to hold the pipette
- Volume adjustment locking for preventing accidental volume changes
- Easy-to-read volume display
- Colour-coding of volume ranges for easy selection of corresponding pipette tips
- Easy user adjustment e.g. for different liquids
- Simple to clean and maintain with only three parts to disassemble
- Fully autoclavable

### Sartorius Optifit Tips

- High-quality standard tips
- Excellent solution for various applications
- Versatile selection of packages and purity options



# Mline® Mechanical Pipette and Tip Multipacks

Multipack Order Code	Mline® Pipettes 1-channel (except where otherwise specified)	Optifit Tips Tip Racks - 96 tips (except where otherwise specified)	Accessories
Mline® Pipette 3-pack 10 LH-725661	<ul style="list-style-type: none"> <li>▪ 0.5 - 10 µl</li> <li>▪ 10 - 100 µl</li> <li>▪ 100 - 1,000 µl</li> </ul>	<ul style="list-style-type: none"> <li>▪ 0.1 - 10 µl</li> <li>▪ 0.5 - 200 µl</li> <li>▪ 10 - 1,000 µl</li> </ul>	Linear Stand
Mline® Pipette 3-pack 20 LH-725662	<ul style="list-style-type: none"> <li>▪ 2 - 20 µl</li> <li>▪ 20 - 200 µl</li> <li>▪ 100 - 1,000 µl</li> </ul>	<ul style="list-style-type: none"> <li>▪ 0.5 - 200 µl</li> <li>▪ 10 - 1,000 µl</li> </ul>	Linear Stand
Mline® Pipette 4-pack LH-725663	<ul style="list-style-type: none"> <li>▪ 0.5 - 10 µl</li> <li>▪ 10 - 100 µl</li> <li>▪ 20 - 200 µl</li> <li>▪ 100 - 1,000 µl</li> </ul>	<ul style="list-style-type: none"> <li>▪ 0.1 - 10 µl</li> <li>▪ 0.5 - 200 µl</li> <li>▪ 10 - 1,000 µl</li> </ul>	Linear Stand
Mline® Pipette 5-pack LH-725664	<ul style="list-style-type: none"> <li>▪ 2 - 20 µl</li> <li>▪ 10 - 100 µl</li> <li>▪ 20 - 200 µl</li> <li>▪ 100 - 1,000 µl</li> <li>▪ 500 - 5,000 µl</li> </ul>	<ul style="list-style-type: none"> <li>▪ 0.5 - 200 µl</li> <li>▪ 10 - 1,000 µl</li> <li>▪ 100 - 5,000 µl (50 tips)</li> </ul>	Linear Stand
Mline® Pipette 3+1-pack LH-725665	<ul style="list-style-type: none"> <li>▪ 0.5 - 10 µl</li> <li>▪ 20 - 200 µl</li> <li>▪ 100 - 1,000 µl</li> <li>▪ 30 - 300 µl (8-channel)</li> </ul>	<ul style="list-style-type: none"> <li>▪ 0.1 - 10 µl</li> <li>▪ 0.5 - 200 µl</li> <li>▪ 5 - 350 µl</li> <li>▪ 10 - 1,000 µl</li> </ul>	Linear Stand



## Germany

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

For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)

## USA

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

## Finland & Baltics

Sartorius  
Biohit Liquid Handling Oy  
Laipatie 1  
00880 Helsinki  
Phone +358.9.755.951  
Fax +358.9.755.95.292  
[linfo.finland@sartorius.com](mailto:linfo.finland@sartorius.com)