

# Tacta<sup>®</sup> Pipette Multipacks

## The Smart Solution for Your Lab



### Tacta<sup>®</sup> Mechanical Pipettes

- Reliable results, even over lengthy pipetting periods
- Comfortable to hold due to the ergonomically designed handle
- Low pipetting and tip ejection forces
- Controlled and smooth tip ejection with the new Sartorius Optiject<sup>®</sup> technology
- The unique Sartorius Optilock<sup>®</sup> system provides flexibility for volume adjustment and locking
- Calibration adjustment to provide accurate results for various liquid types
- Easy to clean 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

# Tacta® Multipacks

Multipack Order Code	Tacta® Pipettes	Optifit Tips	Accessories
Tacta® Pipette 3-pack 10 LH-729670	1-channel – 0.5-10 µl – 10-100 µl – 100-1 000 µl	Tip Racks – 96 tips – 0.1-10 µl – 0.5-200 µl – 10-1 000 µl	Linear Stand
Tacta® Pipette 3-pack 20 LH-729671	1-channel – 2-20 µl – 20-200 µl – 100-1 000 µl	Tip Racks – 96 tips – 0.5-200 µl – 10-1 000 µl	Linear Stand
Tacta® Pipette 4-pack 20 LH-729672	1-channel – 0.5-10 µl – 2-20 µl – 20-200 µl – 100-1 000 µl	Tip Racks – 96 tips – 0.1-10 µl – 0.5-200 µl – 10-1 000 µl	Linear Stand
Tacta® Pipette 4-pack 100 LH-729673	1-channel – 0.5-10 µl – 10-100 µl – 20-200 µl – 100-1 000 µl	Tip Racks – 96 tips – 0.1-10 µl – 0.5-200 µl – 10-1 000 µl	Linear Stand
Tacta® Pipette 5-pack 10 LH-729674	1-channel – 0.5-10 µl – 10-100 µl – 20-200 µl – 100-1 000 µl – 500-5 000 µl	Tip Racks – 96 tips – 0.1-10 µl – 0.5-200 µl – 10-1 000 µl Tip Rack – 50 tips – 100-5 000 µl	Linear Stand
Tacta® Pipette 5-pack 20 LH-729675	1-channel – 2-20 µl – 10-100 µl – 20-200 µl – 100-1 000 µl – 500-5 000 µl	Tip Racks – 96 tips – 0.5-200 µl – 10-1 000 µl Tip Rack – 50 tips – 100-5 000 µl	Linear Stand
Tacta® Pipette 3+1-pack LH-729676	1-channel – 0.5-10 µl – 20-200 µl – 100-1 000 µl 8-channel – 30-300 µl	Tip Racks – 96 tips – 0.1-10 µl – 0.5-200 µl – 5-350 µl – 10-1 000 µl	Linear Stand
Tacta® Pipette 3-pack max LH-729677	1-channel – 100-1 000 µl – 500-5 000 µl – 1 000-10 000 µl	Tip Racks – 10-1 000 µl – 96 tips – 100-5 000 µl – 100 tips – 100-10 000 µl – 35 tips	Linear Stand

## USA

Sartorius Corporation  
5 Orville Drive, Suite 200  
Bohemia, NY 11716  
Phone +1.631.254.4249  
Toll-free +1.800.635.2906  
www.sartorius.us

## Canada

Sartorius Canada, Inc.  
2179 Dunwin Drive, Units 4 & 5  
Mississauga, ON L5L 1X2, Canada  
Phone +1.905.569.7977  
Toll-free +1.800.668.4234  
www.sartorius.com

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

Sartorius Lab Instruments GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen  
Germany  
Phone +49.551.308.0  
www.sartorius.com

