

Tacta® Pipette Multipacks The Smart Solution for Your Lab

Tacta® 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® technology
- The unique Sartorius Optilock® 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 1 – 10 µL 10 – 100 µL 100 – 1,000 µL	Tip Racks – 96 tips 10 µL 200 µL 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 200 µL 1,000 µL	Linear Stand
Tacta® Pipette 4-pack 20 LH-729672	1-channel 1 – 10 µL 2 – 20 µL 20 – 200 µL 100 – 1,000 µL	Tip Racks – 96 tips 10 µL 200 µL 1,000 µL	Linear Stand
Tacta® Pipette 4-pack 100 LH-729673	1-channel 1 – 10 µL 10 – 100 µL 20 – 200 µL 100 – 1,000 µL	Tip Racks – 96 tips 10 µL 200 µL 1,000 µL	Linear Stand
Tacta® Pipette 5-pack 10 LH-729674	1-channel 1 – 10 µL 10 – 100 µL 20 – 200 µL 100 – 1,000 µL 500 – 5,000 µL	Tip Racks – 96 tips 10 µL 200 µL 1,000 µL Tip Rack – 50 tips 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 200 µL 1,000 µL Tip Rack – 50 tips 5,000 µL	Linear Stand
Tacta® Pipette 3+1-pack LH-729676	1-channel 1 – 10 µL 20 – 200 µL 100 – 1,000 µL 8-channel 30 – 300 µL	Tip Racks – 96 tips 10 µL 200 µL 350 µL 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 1,000 µL – 96 tips 5,000 µL – 50 tips 10,000 µL – 35 tips	Linear Stand



Germany

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

USA

Sartorius Corporation
3874 Research Park Drive
Ann Arbor, MI 48108
Phone +1 734 769 1600

For further information, visit
www.sartorius.com