

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 <ul style="list-style-type: none"> ▪ 0.5–10 µL ▪ 10–100 µL ▪ 100–1,000 µL 	Tip Racks – 96 tips <ul style="list-style-type: none"> ▪ 0.1–10 µL ▪ 0.5–200 µL ▪ 10–1,000 µL 	Linear Stand
Tacta® Pipette 3-pack 20 LH-729671	1-channel <ul style="list-style-type: none"> ▪ 2–20 µL ▪ 20–200 µL ▪ 100–1,000 µL 	Tip Racks – 96 tips <ul style="list-style-type: none"> ▪ 0.5–200 µL ▪ 10–1,000 µL 	Linear Stand
Tacta® Pipette 4-pack 20 LH-729672	1-channel <ul style="list-style-type: none"> ▪ 0.5–10 µL ▪ 2–20 µL ▪ 20–200 µL ▪ 100–1,000 µL 	Tip Racks – 96 tips <ul style="list-style-type: none"> ▪ 0.1–10 µL ▪ 0.5–200 µL ▪ 10–1,000 µL 	Linear Stand
Tacta® Pipette 4-pack 100 LH-729673	1-channel <ul style="list-style-type: none"> ▪ 0.5–10 µL ▪ 10–100 µL ▪ 20–200 µL ▪ 100–1,000 µL 	Tip Racks – 96 tips <ul style="list-style-type: none"> ▪ 0.1–10 µL ▪ 0.5–200 µL ▪ 10–1,000 µL 	Linear Stand
Tacta® Pipette 5-pack 10 LH-729674	1-channel <ul style="list-style-type: none"> ▪ 0.5–10 µL ▪ 10–100 µL ▪ 20–200 µL ▪ 100–1,000 µL ▪ 500–5,000 µL 	Tip Racks – 96 tips <ul style="list-style-type: none"> ▪ 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 <ul style="list-style-type: none"> ▪ 2–20 µL ▪ 10–100 µL ▪ 20–200 µL ▪ 100–1,000 µL ▪ 500–5,000 µL 	Tip Racks – 96 tips <ul style="list-style-type: none"> ▪ 0.5–200 µL ▪ 10–1,000 µL Tip Rack – 50 tips <ul style="list-style-type: none"> ▪ 100–5,000 µL 	Linear Stand
Tacta® Pipette 3+1-pack LH-729676	1-channel <ul style="list-style-type: none"> ▪ 0.5–10 µL ▪ 20–200 µL ▪ 100–1,000 µL 8-channel <ul style="list-style-type: none"> ▪ 30–300 µL 	Tip Racks – 96 tips <ul style="list-style-type: none"> ▪ 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 <ul style="list-style-type: none"> ▪ 100–1,000 µL ▪ 500–5,000 µL ▪ 1,000–10,000 µL 	Tip Racks <ul style="list-style-type: none"> ▪ 10–1,000 µL – 96 tips ▪ 100–5,000 µL – 50 tips ▪ 100–10,000 µL – 35 tips 	Linear Stand



Germany

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

For further information, visit
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
Laippatie 1
00880 Helsinki
Phone +358.9.755.951
Fax +358.9.755.95.292
linfo.finland@sartorius.com