

Proline[®] Plus Pipette Multipacks Reliable Combo-Pack for Your Lab

Proline[®] Plus

Mechanical Pipettes

- Low pipetting forces
- Ergonomic finger hook
- Easy volume setting with click-stop mechanism
- Easy-to-read volume display
- 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



Proline® Plus Multipacks

Multipack Order Code	Proline® Plus Pipettes 1-Channel (except where otherwise specified)	Optifit Tips Tip Racks - 96 tips (except where otherwise specified)	Accessories
Proline® Plus Pipette 3-pack 10 LH-728670	<ul style="list-style-type: none"> ▪ 0.5 - 10 µl ▪ 10 - 100 µl ▪ 100 - 1,000 µl 	<ul style="list-style-type: none"> ▪ 0.1 - 10 µl ▪ 0.5 - 200 µl ▪ 10 - 1,000 µl 	Linear Stand
Proline® Plus Pipette 3-pack 20 LH-728671	<ul style="list-style-type: none"> ▪ 2 - 20 µl ▪ 20 - 200 µl ▪ 100 - 1,000 µl 	<ul style="list-style-type: none"> ▪ 0.5 - 200 µl ▪ 10 - 1,000 µl 	Linear Stand
Proline® Plus Pipette 4-pack 20 LH-728672	<ul style="list-style-type: none"> ▪ 0.5 - 10 µl ▪ 2 - 20 µl ▪ 20 - 200 µl ▪ 100 - 1,000 µl 	<ul style="list-style-type: none"> ▪ 0.1 - 10 µl ▪ 0.5 - 200 µl ▪ 10 - 1,000 µl 	Linear Stand
Proline® Plus Pipette 4-pack 100 LH-728673	<ul style="list-style-type: none"> ▪ 0.5 - 10 µl ▪ 10 - 100 µl ▪ 20 - 200 µl ▪ 100 - 1,000 µl 	<ul style="list-style-type: none"> ▪ 0.1 - 10 µl ▪ 0.5 - 200 µl ▪ 10 - 1,000 µl 	Linear Stand
Proline® Plus Pipette 5-pack 10 LH-728674	<ul style="list-style-type: none"> ▪ 0.5 - 10 µl ▪ 10 - 100 µl ▪ 20 - 200 µl ▪ 100 - 1,000 µl ▪ 500 - 5,000 µl 	<ul style="list-style-type: none"> ▪ 0.1 - 10 µl ▪ 0.5 - 200 µl ▪ 10 - 1,000 µl ▪ 100 - 5,000 µl (50 tips) 	Linear Stand
Proline® Plus Pipette 5-pack 20 LH-728675	<ul style="list-style-type: none"> ▪ 2 - 20 µl ▪ 10 - 100 µl ▪ 20 - 200 µl ▪ 100 - 1,000 µl ▪ 500 - 5,000 µl 	<ul style="list-style-type: none"> ▪ 0.5 - 200 µl ▪ 10 - 1,000 µl ▪ 100 - 5,000 µl (50 tips) 	Linear Stand
Proline® Plus Pipette 3+1-pack LH-728676	<ul style="list-style-type: none"> ▪ 0.5 - 10 µl ▪ 20 - 200 µl ▪ 100 - 1,000 µl ▪ 30 - 300 µl (8-channel) 	<ul style="list-style-type: none"> ▪ 0.1 - 10 µl ▪ 0.5 - 200 µl ▪ 5 - 350 µl ▪ 10 - 1,000 µl 	Linear Stand
Proline® Plus Pipette 3-pack max LH-728677	<ul style="list-style-type: none"> ▪ 100 - 1,000 µl ▪ 500 - 5,000 µl ▪ 1,000 - 10,000 µl 	<ul style="list-style-type: none"> ▪ 10 - 1,000 µl ▪ 100 - 5,000 µl (50 tips) ▪ 500 - 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 contacts, 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
Laipatie 1
00880 Helsinki
Phone +358.9.755.951
Fax +358.9.755.95.292
linfo.finland@sartorius.com