

Pipette Upgrade

Upgrade Your Pipetting and Save up to 40%

NEW Picus® 2 Electronic Pipettes

Picus® 2 pipettes offer reliable performance, while also being kind to your hand. The unmatched ergonomic design not only reduces strain, but also helps to ensure reproducible pipetting results. This combined with new connectivity features and Sartorius Pipetting mobile app truly takes the productivity to the new level!

Channel	Sartorius Part no.	Vol. Range	Sugg. List	Special Price
1-Ch	LH-747021	0.2–10 µl	\$1,150.00	\$805.00
1-Ch	LH-747041	5–120 µl	\$1,150.00	\$805.00
1-Ch	LH-747061	10–300 µl	\$1,150.00	\$805.00
1-Ch	LH-747081	50–1,000 µl	\$1,150.00	\$805.00
1-Ch	LH-747101	100–5,000 µl	\$1,150.00	\$805.00
1-Ch	LH-747111	500–10,000 µl	\$1,150.00	\$805.00
8-Ch	LH-747321	0.2–10 µl	\$307.00	\$184.40
8-Ch	LH-747341	5–120 µl	\$307.00	\$184.40
8-Ch	LH-747361	10–300 µl	\$307.00	\$182.40
8-Ch	LH-747391	50–1,000 µl	\$840.00	\$504.00
12-Ch	LH-747421	0.2–10 µl	\$840.00	\$504.00
12-Ch	LH-747441	5–120 µl	\$840.00	\$504.00
12-Ch	LH-747461	10–300 µl	\$1,020.00	\$612.00
12-Ch	LH-747491	50–1,000 µl	\$1,020.00	\$612.00

Save up to 40% off



Tacta® Mechanical Pipettes

Channel	Sartorius Part no.	Vol. Range	Sugg. List	Special Price
1-Ch	LH-729010	0.1–3 µl	\$516.00	\$335.40
1-Ch	LH-729020	0.5–10 µl	\$516.00	\$335.40
1-Ch	LH-729030	2–20 µl	\$516.00	\$335.40
1-Ch	LH-729050	10–100 µl	\$516.00	\$335.40
1-Ch	LH-729060	20–200 µl	\$516.00	\$335.40
1-Ch	LH-729070	100–1000 µl	\$516.00	\$335.40
1-Ch	LH-729080	500–5000 µl	\$516.00	\$335.40
1-Ch	LH-729090	1–10 ml	\$516.00	\$335.40
8-Ch	LH-729120	0.5–10 µl	\$1,300.00	\$845.00
8-Ch	LH-729130	5–100 µl	\$1,300.00	\$845.00
8-Ch	LH-729140	30–300 µl	\$1,300.00	\$845.00
12-Ch	LH-729220	0.5–10 µl	\$1,530.00	\$994.00
12-Ch	LH-729230	5–100 µl	\$1,530.00	\$994.00
12-Ch	LH-729240	30–300 µl	\$1,530.00	\$994.00

- Comfortable to hold due to the ergonomically designed handle
- The unique Sartorius Optilock® system provides flexibility for volume adjustment and locking
- Controlled and smooth tip ejection with the new Sartorius Optiject® technology
- Calibration adjustment to provide accurate results for various liquid types
- Easy to clean, with only three parts to disassemble



Proline® Plus Mechanical Pipettes

Channel	Sartorius Part no.	Vol. Range	Sugg. List	Special Price
1-Ch	728010	0.1–3 µl	\$307.00	\$184.40
1-Ch	728020	0.5–10 µl	\$307.00	\$184.40
1-Ch	728030	2–20 µl	\$307.00	\$184.40
1-Ch	728040	5–50 µl	\$307.00	\$184.40
1-Ch	728050	10–100 µl	\$307.00	\$184.40
1-Ch	728060	20–200 µl	\$307.00	\$184.40
1-Ch	728070	100–1000 µl	\$307.00	\$184.40
1-Ch	728080	500–5000 µl	\$307.00	\$184.40
1-Ch	728090	1–10 ml	\$307.00	\$182.40
8-Ch	728120	0.5–10 µl	\$840.00	\$504.00
8-Ch	728130	10–100 µl	\$840.00	\$504.00
8-Ch	728140	30–300 µl	\$840.00	\$504.00
12-Ch	728220	0.5–10 µl	\$1,020.00	\$612.00
12-Ch	728230	10–100 µl	\$1,020.00	\$612.00
12-Ch	728240	30–300 µl	\$1,020.00	\$612.00

- Ergonomic finger support minimizes the grip force needed to hold the pipette
- Optiload mechanism in multichannel models for easy and light tip loading with perfect tip sealing
- Volume range from 3µL (5µL for fixed) to 10mL
- Color-coding of different volumes to ease the selection of matching pipette tips
- Simple to clean and maintain with only three parts to disassemble



Special offer valid only for the US from January 1, 2024 to December 31, 2024 and cannot be combined with other offers.

Contact: