

## 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     | \$1,440.00 | \$1,008.00    |
| 1-Ch    | LH-747041          | 5–120 µl      | \$1,440.00 | \$1,008.00    |
| 1-Ch    | LH-747061          | 10–300 µl     | \$1,440.00 | \$1,008.00    |
| 1-Ch    | LH-747081          | 50–1,000 µl   | \$1,440.00 | \$1,008.00    |
| 1-Ch    | LH-747101          | 100–5,000 µl  | \$1,440.00 | \$1,008.00    |
| 1-Ch    | LH-747111          | 500–10,000 µl | \$1,440.00 | \$1,008.00    |
| 8-Ch    | LH-747321          | 0.2–10 µl     | \$2,450.00 | \$1,715.00    |
| 8-Ch    | LH-747341          | 5–120 µl      | \$2,450.00 | \$1,715.00    |
| 8-Ch    | LH-747361          | 10–300 µl     | \$2,450.00 | \$1,715.00    |
| 8-Ch    | LH-747391          | 50–1,000 µl   | \$2,450.00 | \$1,715.00    |
| 12-Ch   | LH-747421          | 0.2–10 µl     | \$3,060.00 | \$2,142.00    |
| 12-Ch   | LH-747441          | 5–120 µl      | \$3,060.00 | \$2,142.00    |
| 12-Ch   | LH-747461          | 10–300 µl     | \$3,060.00 | \$2,142.00    |
| 12-Ch   | LH-747491          | 50–1,000 µl   | \$3,060.00 | \$2,142.00    |

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    | \$644.00   | \$418.60      |
| 1-Ch    | LH-729020          | 0.5–10 µl   | \$644.00   | \$418.60      |
| 1-Ch    | LH-729030          | 2–20 µl     | \$644.00   | \$418.60      |
| 1-Ch    | LH-729050          | 10–100 µl   | \$644.00   | \$418.60      |
| 1-Ch    | LH-729060          | 20–200 µl   | \$644.00   | \$418.60      |
| 1-Ch    | LH-729070          | 100–1000 µl | \$644.00   | \$418.60      |
| 1-Ch    | LH-729080          | 500–5000 µl | \$644.00   | \$418.60      |
| 1-Ch    | LH-729090          | 1–10 ml     | \$644.00   | \$418.60      |
| 8-Ch    | LH-729120          | 0.5–10 µl   | \$1,620.00 | \$1,053.00    |
| 8-Ch    | LH-729130          | 5–100 µl    | \$1,620.00 | \$1,053.00    |
| 8-Ch    | LH-729140          | 30–300 µl   | \$1,620.00 | \$1,053.00    |
| 12-Ch   | LH-729220          | 0.5–10 µl   | \$1,910.00 | \$1,241.50    |
| 12-Ch   | LH-729230          | 5–100 µl    | \$1,910.00 | \$1,241.50    |
| 12-Ch   | LH-729240          | 30–300 µl   | \$1,910.00 | \$1,241.50    |

- 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    | \$384.00   | \$230.40      |
| 1-Ch    | 728020             | 0.5–10 µl   | \$384.00   | \$230.40      |
| 1-Ch    | 728030             | 2–20 µl     | \$384.00   | \$230.40      |
| 1-Ch    | 728040             | 5–50 µl     | \$384.00   | \$230.40      |
| 1-Ch    | 728050             | 10–100 µl   | \$384.00   | \$230.40      |
| 1-Ch    | 728060             | 20–200 µl   | \$384.00   | \$230.40      |
| 1-Ch    | 728070             | 100–1000 µl | \$384.00   | \$230.40      |
| 1-Ch    | 728080             | 500–5000 µl | \$384.00   | \$230.40      |
| 1-Ch    | 728090             | 1–10 ml     | \$384.00   | \$230.40      |
| 8-Ch    | 728120             | 0.5–10 µl   | \$1,050.00 | \$630.00      |
| 8-Ch    | 728130             | 10–100 µl   | \$1,050.00 | \$630.00      |
| 8-Ch    | 728140             | 30–300 µl   | \$1,050.00 | \$630.00      |
| 12-Ch   | 728220             | 0.5–10 µl   | \$1,280.00 | \$768.00      |
| 12-Ch   | 728230             | 10–100 µl   | \$1,280.00 | \$768.00      |
| 12-Ch   | 728240             | 30–300 µl   | \$1,280.00 | \$768.00      |

- 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 3 µL (5 µL for fixed) to 10 mL
- 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 Canada from January 1, 2024 to December 31, 2024 and cannot be combined with other offers.

Contact:

Your Logo