

eLINE® Electronic Pipettes

Proven Convenience, Reliability and Speed



Description

The electronic eLINE® pipette family offers a great price-performance ratio for any user. Its ergonomic design eases the users workload and reduces the risk of strain injuries. The fully electronic operation secures accurate and precise results independent of the user. In addition, the versatile selection of liquid handling modes supports and speeds up the user's work. One eLINE® can be used for several liquid handling tasks from regular pipetting to diluting with mixing and automatic repetitive dispensing.

Features

For convenience:

- Fully electronic operation minimizes the force needed to operate the pipette
- Electronic tip ejection ensures light removal of tips also with multichannel models
- The ergonomic design of the handle, and the easy-to-reach operating buttons, allow for convenient single-handed operation for both right and left-handed users
- The convenient finger hook provides proper support to minimize grip force

For reliable results:

- Electronic piston movement, electronic brake and piston control guarantee accurate and reliable pipetting results, independent of the user

- The Optoload feature permits tip loading with perfect tip sealing onto every tip cone for secured accuracy
- Safe-Cone Filters prevent the risk of contamination cost-effectively
- Unique Super Pipetting feature with the 0,1–5 µl model shoots the last droplets out – no need to touch the vessel wall

For speed:

- The versatile pipetting modes reduce the needed work stages, enabling significantly faster liquid handling than a mechanical pipette
- Multichannel models enable pipetting on the microwell plates considerably faster than the mechanical pipettes
- Speed setting available for both aspiration and dispensing

Applications

- PCR and other DNA/RNA techniques
- ELISA
- Protein analysis
- Cell culture

Intended use

For fully electronic liquid dispensing, mixing and diluting in volumes under 5 ml.

Technical Specifications

Rechargeable battery	NiMH
Charging time	Approx. 12 hours (4 hours quick charging)
Charger	Universal charger with EU, US JPN, UK, CHN, AUS and KOR plugs
Weight	170 g (1-ch, 300 µl) 223 g (8-ch, 300 µl)
Length	241 mm (1-ch, 300 µl) 240 mm (8-ch, 300 µl)
Number of pipetting cycles	>1000
Volume range	1-ch: 0.1–5000 µl 8- and 12-ch: 0.2–1200 µl
Pipetting modes	7+3
DC-motor concept	Piston control Electronic brake
Memory places	6
Tip ejection	Electronic
Spring loaded tip cones	Optoload feature in multichannel models
Filters	Safe-Cone Filters in all models >10 µl
Autoclavable lower parts*	121°C, 20 min, 1 bar
Charging Stands, available separately	Charging Stand for 1 pipette, Charging Carousel for 4 pipettes
Warranty	2 years, possibility for 1 year extended warranty

*Excluding 1200 µl multichannel models

Pipetting Modes and Mode Additions

- Pipetting (P)
- Reverse pipetting (rP)
- Pipetting with mixing (P mix)
- Pipetting with cycle counter (P count)
- Manual pipetting (P man), with 1-ch models
- Multiple dispensing (d)
- Sequential dispensing (Sd), excl. 0.1–0.5 µl model
- Automatic multi dispensing (Ad)
- Diluting (dd)
- Diluting with mixing (dd mix)
- Multi-aspirating (SA)
- Super Pipetting (SP), with 0.1–5 µl model

Ordering Information

Order Code	Channels		Volume Range (μl)	Increment (μl)	Test Volume (μl)	System. Error (%)	Random Error (%)	Safe-Cone Standard	Filters Plus	Optifit Tips** (μl)	SafetySpace Filter Tips** (μl)
730011	1		0.1-5	0.05	5 2.5 0.5	1.20 2.00 12.00	0.70 1.20 8.00	-	-	10	10
730021	1		0.2-10	0.05	10 5 1	1.00 1.40 4.00	0.40 0.70 2.30	-	-	10	10
730041	1		5-120	0.50	120 60 12	0.60 0.90 3.00	0.20 0.20 1.00	721008	721018	200 350	120
730061	1		10-300	1.00	300 150 30	0.60 0.60 1.50	0.15 0.20 0.80	721007	721017	350	300
730081	1		50-1000	5.00	1000 500 100	0.40 0.60 1.60	0.15 0.20 0.50	721006	721016	1000	1000
730101*	1		100-5000	10.0	5000 2500 500	0.50 0.80 1.60	0.17 0.20 0.40	721006	721016	5000	-
730321	8		0.2-10	0.05	10 5 1	0.90 1.50 4.00	0.60 0.90 3.50	-	-	10	10
730341	8		5-120	0.50	120 60 12	0.80 0.80 3.00	0.30 0.35 1.50	721008	721018	200 350	120
730361	8		10-300	1.00	300 150 30	0.70 0.90 3.00	0.20 0.30 1.20	721007	721017	350	300
730391	8		50-1200	5.00	1200 600 120	0.90 1.20 2.50	0.30 0.30 0.80	721006	721016	1200	1200
730421	12		0.2-10	0.05	10 5 1	0.90 1.50 4.00	0.60 0.90 3.50	-	-	10	10
730441	12		5-120	0.50	120 60 12	0.80 0.80 3.00	0.30 0.35 1.50	721008	721018	200 350	300
730461	12		10-300	1.00	300 150 30	0.70 0.90 3.00	0.20 0.30 1.20	721007	721017	350	300
730491	12		50-1200	5.00	1200 600 120	0.90 1.20 2.50	0.30 0.30 0.80	721006	721016	1200	1200

All pipettes include universal charger (EU, US | JPN, UK, CHN, AUS, KOR plugs)

*NOTE: Min. volume in P-mode is 500 µl. 100 µl is possible in d-mode.

**NOTE: Low Retention Tips are available in volumes up to 1200 µl.

Sartorius Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki, Finland
Phone +358.9.755.951
Fax +358.9.755.95.200

Sartorius Lab Instruments GmbH & Co. KG
Weender Landstrasse 94-108
37075 Goettingen, Germany
Phone +49.551.3080
Fax +49.551.308.3289

www.sartorius.com

Specifications subject to change without notice.
Printed and copyrighted by
Sartorius Biohit Liquid Handling Oy
Publication No.: SUL2002-e140102
Order No.: 85037-543-80
Ver. 2|2015

