

Picus® & Picus® NxT Electronic Pipettes

The Most Sophisticated and Ergonomic Pipettes Ever!



Description

Sartorius Picus® & Picus® NxT are the most sophisticated and ergonomic electronic pipettes on the market. These exceptionally compact and lightweight pipettes have been specially designed to ease the user's workload and to protect the user from repetitive strain injury (RSI). The newest technology and fully electronic operation assure high accuracy and precision of results, regardless of user skill level. Picus® NxT provides distinct advantages for highly regulated laboratories.

Features

Picus® & Picus® NxT

- Highest level of ergonomics provided by the uniquely low weight, light electronic tip ejection and comfortable handle design
- Extensive range of pipetting modes reduces the needed pipetting steps and speeds up work
- Electronic brake and piston control system provide outstanding accuracy and repeatability of pipetting results, independent of the user
- Intuitive user interface in five language options: English, French, German, Russian and Chinese, enables ease of use
- Adjustment wheel offers extremely fast volume setting and menu navigation
- Optiload enables perfect tip sealing for accurate delivery from each channel
- Safe-Cone Filters prevent the risk of contamination cost-effectively

- Microwell plate tracker guides the user to pipette into the correct wells
- Calibration adjustment in 1, 2 or 3 points

Picus® NxT

- Certificate of accredited 3-point calibration (per ISO 17025 and ISO 8655) delivered with the product at no extra charge
- User programmable pipetting protocols enable the storage of three frequently needed pipetting workflows; easily activated when needed.
- 2-level password protection for stored programs to prevent unauthorized changes (optional)
- Pipette locking, e.g. in case of contamination, increases lab safety by disabling the pipette from use.
- Service and calibration reminders help the users to remember important service dates.
- Repeated blow-out helps to dispense the last droplets of e.g. viscous liquids

Applications

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

Intended use

Fully electronic liquid handling in the volume range of 0.2 µl to 10 ml.

Technical Specifications

Rechargeable battery	Li-Polymer with protection circuit
Charging time	Approx. 1 hour
Charger	Universal charger with EU, US JPN, UK, CHN, AUS and KOR plugs
Weight	100 g (1-ch, 300 µl) 160 g (8-ch, 300 µl)
Length	210 mm (1-ch, 300 µl) 216 mm (8-ch, 300 µl)
Number of pipetting cycles	>1000
Volume range	1-ch: 0.2–10 000 µl 8- & 12-ch: 0.2–1200 µl
Pipetting modes	Picus®: 8 + 6 Picus® NxT: 9 + 7
DC-motor concept	Electronic piston control Electronic brake
Memory places	Picus®: 10 Picus® NxT: 3* + 10
Tip ejection	Electronic
Spring loaded tip cones	Optiload 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

*For Protocols

**Excluding 1200 µl multichannel models

Pipetting Modes	Advanced Functions
Pipetting	Tracker, Mixing, Counter, Repeated Blow-out*
Reverse Pipetting	Tracker, Counter, Excess Volume Adjustment
Manual Pipetting	Repeated Blow-out*
Multi-Dispensing	Tracker, Excess Volume Adjustment, Auto-Dispensing
Diluting	Mixing, Repeated Blow-out*
Sequential Dispensing	Excess Volume Adjustment
Multi-Aspiration	Repeated Blow-out*
Titrate	Fast Dispensing
Protocol*	All additional modes

*Advanced function, Repeated Blow-out, and pipetting mode, Protocol, are only available in Picus® NxT models.

Ordering Information

Picus® NxT Order Code	Picus® 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)
LH-745021	735021	1 	0,2-10	0,01	10 5 1	0,90 1,00 2,50	0,40 0,70 1,50	-	-	10	10
LH-745041	735041	1 	5-120	0,10	120 60 12	0,40 0,60 2,00	0,15 0,20 1,00	721008	721018	200 350	120
LH-745061	735061	1 	10-300	0,20	300 150 30	0,40 0,60 1,50	0,15 0,20 0,80	721007	721017	350	300
LH-745081	735081	1 	50-1000	1,00	1000 500 100	0,40 0,60 1,50	0,15 0,20 0,50	721006	721016	1000	1000
LH-745101	735101	1 	100-5000	5,00	5000 2500 500	0,50 0,80 1,00	0,15 0,20 0,40	721005	721015	5000	-
LH-745111	735111	1 	500-10000	10,00	10000 5000 1000	0,60 1,20 3,00	0,20 0,30 0,60	721005	721015	10000	-
LH-745321	735321	8 	0,2-10	0,01	10 5 1	0,90 1,50 4,00	0,50 0,80 3,00	-	-	10	10
LH-745341	735341	8 	5-120	0,10	120 60 12	0,50 0,70 2,00	0,20 0,30 1,50	721008	721018	200 350	120
LH-745361	735361	8 	10-300	0,20	300 150 30	0,50 0,70 2,00	0,20 0,30 1,00	721007	721017	350	300
LH-745391	735391	8 	50-1200	1,00	1200 600 120	0,50 1,00 2,50	0,20 0,30 1,00	721006	721016	1200	1200
LH-745421	735421	12 	0,2-10	0,01	10 5 1	0,90 1,50 4,00	0,50 0,80 3,00	-	-	10	10
LH-745441	735441	12 	5-120	0,10	120 60 12	0,50 0,70 2,00	0,20 0,30 1,50	721008	721018	200 350	120
LH-745461	735461	12 	10-300	0,20	300 150 30	0,50 0,70 2,00	0,20 0,30 1,00	721007	721017	350	300
LH-745491	735491	12 	50-1200	1,00	1200 600 120	0,50 1,00 2,50	0,20 0,30 1,00	721006	721016	1200	1200

*The listed systematic and random error values are valid for Pipetting mode only, and achieved under strictly controlled conditions during type tests per ISO 8655. Due to the continuous product development by Sartorius, the systematic and random error values may change without prior notice.

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

All pipettes are supplied with a universal charger (EU, UK, US | JPN, KOR, AUS and CHN plugs).

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

