

Picus® & Picus® Nxt Electronic Pipettes

The Most Sophisticated
and Ergonomic Pipettes Ever!

Product Information

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).



Description

The Picus® family pipettes are kind to your hand with unbeatable ergonomic design that ensures reliable and repeatable experiment results. Repeatable pipetting results are guaranteed with the electronic piston control and brake, raising all users to expert 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

- 2 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.

Fully electronic liquid handling in the volume range of 0.2 µL to 10 mL.

Technical Data

Technical Specifications	
Rechargeable battery	Li-Polymer with protection circuit
Charging time	Approx. 1 hour
Charger	Universal charger with EU, US JPN, UK, CI IN, 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	>1,000
Volume range	1-ch: 0.2 - 10,000 µL 8- & 12 - ch: 0.2 - 1200 µL
Pipetting modes	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	Picus®	Channels	Volume Range (µL)	Increment (µL)	Test Volume (µL)	Mode ^{P/D}	Systematic Error ^N Limit ± (%) (µL)		Random Error ^N Limit (%) (µL)		
LH-745021	735021	1	■	0.2 - 10	0.01	10	P	1.0	0.100	0.4	0.040
						5	P	1.2	0.060	0.7	0.035
						1	P	3.0	0.030	2.0	0.020
						0.2	P	17.5	0.035	10	0.020
						1	D	6.0	0.060	7.0	0.070
LH-745041	735041	1	■	5 - 120	0.10	120	P	0.5	0.60	0.15	0.18
						60	P	0.7	0.42	0.2	0.12
						12	P	2.0	0.24	1.0	0.12
						5	P	5.5	0.275	2.5	0.125
						12	D	4.0	0.48	4.0	0.48
LH-725061	735061	1	■	10 - 300	0.20	300	P	0.5	1.50	0.15	0.45
						150	P	0.6	0.90	0.2	0.30
						30	P	1.5	0.45	0.8	0.24
						10	P	5.0	0.50	2.4	0.24
						30	D	3.0	0.90	3.0	0.90
LH-745081	735081	1	■	50 - 1,000	1.00	1,000	P	0.45	4.5	0.15	1.5
						500	P	0.6	3.0	0.2	1.0
						100	P	2.0	2.0	0.5	0.5
						50	P	4.0	2.0	1.0	0.5
						100	D	2.5	2.5	2.0	2.0
LH-745101	735101	1	■	100 - 5,000	5.00	5,000	P	0.5	25	0.15	7.5
						2,500	P	0.7	17.5	0.2	5
						500	P	1.6	8	0.4	2
						100	P	8.0	8	2.0	2
						500	D	2.4	12	2.4	12
LH-745111	735111	1	■	500-10,000	10.00	10,000	P	0.6	60	0.2	20
						5,000	P	0.9	45	0.3	15
						1,000	P	3.0	30	0.6	6
						500	P	7.0	35	1.2	6
						1,000	D	4.0	40	2.4	24
LH-745321	735321	8	■	0.2 - 10	0.01	10	P	0.6	0.72	0.5	0.050
LH-745421	735421	12				5	P	0.8	0.48	0.8	0.040
						1	P	2.5	0.30	3.0	0.030
						0.2	P	6.0	0.30	15.0	0.030
						1	D	4.5	0.54	15.0	0.150
LH-745341	735341	8	■	5 - 120	0.10	120	P	0.6	0.72	0.3	0.36
LH-745441	735441	12				60	P	0.8	0.48	0.4	0.24
						12	P	2.5	0.30	1.67	0.20
						5	P	6.0	0.30	4.0	0.20
						12	D	4.5	0.54	8.0	0.96
LH-745361	735361	8	■	10 - 300	0.20	300	P	0.6	0.150	1.0	0.100
LH-745461	735461	12				150	P	0.8	0.125	2.0	0.100
						30	P	2.33	0.055	4.0	0.040
						10	P	8.0			
						30	D	3.33			
LH-745391	735391	8	■	50-1,200	1.00	1,200	P	0.6	7.2	0.2	2.4
LH-745491	735491	12				600	P	1.0	6.0	0.3	1.8
						120	P	2.5	3.0	1.0	1.2
						50	P	8.0	4.0	2.4	1.2
						20	D	3.33	4.0	3.33	4.0

^N Note: The listed systematic and random error values can be achieved only under strictly controlled conditions during type test per ISO 8655. The best compatibility is achieved when combining Sartorius pipettes and Sartorius tips. The systematic error and random error results, in tests, have been achieved using Sartorius Optifit tips at factory default speed settings. Due to the continuous product development by Sartorius, the systematic and random error values are subject to change without prior notice.

^P P = Pipetting Mode

^D D = Multi-dispensing mode. The listed systematic and random error values are of 10 measurements at 10 % of the nominal volume.

Germany


Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

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
Laippatie 1
00880 Helsinki
Phone +358 9 755 951

 For further information, visit
www.sartorius.com