



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 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 Data

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

Pipetting Modes

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 Functions

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

* For Protocols

** Excluding 1200 µl multichannel models



Ordering Information

Picus® NxT	Picus®	Channels		Volume Range (µl)	Increment (µl)	Test Volume (µl)	Mode ^{P D}	Systematic Error ^N Limit ±		Random Error ^N Limit	
								(%)	(µl)	(%)	(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-745061	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	1.2	0.120	0.5	0.050
LH-745421	735421	12				5	P	1.5	0.075	0.8	0.040
						1	P	4.0	0.040	3.0	0.030
						0.2	P	25.0	0.050	15.0	0.030
						1	D	12.0	0.120	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	1.80	0.2	0.60
LH-745461	735461	12				150	P	0.8	1.20	0.3	0.45
						30	P	2.33	0.70	1.0	0.30
						10	P	8.0	0.80	3.0	0.30
						30	D	3.33	1.00	6.0	1.80
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
						120	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 tests per ISO 8655.

Pipettes are tested 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.

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

Sales and Service Contacts

For further contacts, visit www.sartorius.com

Europe

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen

Phone +49.551.308.0
Fax +49.551.308.3289

France & Suisse Romande

Sartorius France
2, rue Antoine Laurent de Lavoisier
ZA de la Gaudrée
91410 Dourdan

Phone +33.1.70.62.50.00
Fax +33.1.64.59.76.39

Austria

Sartorius Austria GmbH
Modecenterstrasse 22
1030 Vienna

Phone +43.1.7965760.0
Fax +43.1.7965760.24

Belgium

Sartorius Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles

Phone +32.2.756.06.90
Fax +32.2.481.84.11

Finland & Baltics

Sartorius Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki

Phone +358.9.755.951
Fax +358.9.755.95.200

Hungary

Sartorius Hungária Kft.
Kagyló u. 5.
2092 Budakeszi

Phone +3623.457.227
Fax +3623.457.147

Ireland

Sartorius Ireland Ltd.
Unit 41, The Business Centre
Stadium Business Park
Ballycoolin Road
Dublin 11

Phone +353.1.8089050
Fax +353.1.8089388

Italy

Sartorius Italy S.r.l.
Via Torino 3/5
20814 Varedo (MB)

Phone +39.0362.5557.11
Fax +39.0362.5557.99

Netherlands

Sartorius Netherlands B.V.

Phone +31.30.60.53.001
Fax +31.30.60.52.917

info.netherlands@sartorius.com

Poland

Sartorius Poland sp.z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn

Phone +48.61.6473830
Fax +48.61.6473839

Russian Federation

LLC "Sartorius RUS"
Vasilyevsky Island
5th line 70, Lit. A
199178 St. Petersburg

Phone +7.812.327.53.27
Fax +7.812.327.53.23

Spain & Portugal

Sartorius Spain, S.A.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)

Phone Spain +34.913.586.095
Phone Portugal +351.800.855.800
Fax Spain +34.913.589.623
Fax Portugal +351.800.855.799

Switzerland

Sartorius Mechatronics Switzerland AG
Ringstrasse 24a
8317 Tagelswangen (ZH)

Phone +41.44.746.50.00
Fax +41.44.746.50.50

U.K.

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9QQ

Phone +44.1372.737159
Fax +44.1372.726171

Ukraine

LLS "Sartorius RUS"
Post Box 440 "B"
01001 Kiev, Ukraine

Phone +380.44.411.4918
Fax +380.50.623.3162

Americas

USA

Sartorius Corporation
5 Orville Drive, Suite 200
Bohemia, NY 11716

Phone +1.631.254.4249
Toll-free +1.800.635.2906
Fax +1.631.254.4253

Argentina

Sartorius Argentina S.A.
Int. A. AVALOS 4251
B1605ECS Munro
Buenos Aires

Phone +54.11.4721.0505
Fax +54.11.4762.2333

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900
Fax +55.11.4362.8901

Canada

Sartorius Canada Inc
1173 North Service Road West, D4
Oakville, ON L6M 2V9

Phone +1.905.569.7977
Toll-Free +1.800.668.4234
Fax +1.905.569.7021

Mexico

Sartorius de México, S.A. de C.V.
Libramiento Norte de Tepetzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepetzotlan,
Estado de México,
C.P. 54605

Phone +52.55.5562.1102
Fax +52.55.5562.2942

leadsmex@sartorius.com

Peru

Sartorius Peru S.A.C.
Avenue Alberto del Campo 411
Floor 12 - The Office
15076 - San Isidro, Lima

Phone +51.1.441 0158
Fax +51.1.422 6100

Asia | Pacific

Australia

Sartorius Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175

Phone +61.3.8762.1800
Fax +61.3.8762.1828

China

Sartorius (Shanghai) Trading Co., Ltd.
3rd Floor, North Wing, Tower 1
No. 4560 Jinke Road
Zhangjiang Hi-Tech Park
Pudong District
Shanghai 201210, P.R. China

Phone +86.21.6878.2300
Fax +86.21.6878.2882

Hong Kong

Sartorius Hong Kong Ltd.
Unit 1012, Lu Plaza
2 Wing Yip Street
Kwun Tong
Kowloon, Hong Kong

Phone +852.2774.2678
Fax +852.2766.3526

India

Sartorius Weighing India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra,
Nelamangala Tq
562 123 Bangalore, India

Phone +91.80.4350.5250
Fax +91.80.4350.5253

Japan

Sartorius Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.3740.5408
Fax +81.3.3740.5406

Malaysia

Sartorius Malaysia Sdn. Bhd
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622
Fax +60.3.8996.0755

Singapore

Sartorius Singapore Pte. Ltd
10 Science Park Rd
The Alpha #02-13/14
Singapore Science Park II
Singapore 117684

Phone +65.6872.3966
Fax +65.6778.2494

South Korea

Sartorius Korea Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, Bundang-Gu
SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700
Fax +82.31.622.5799

Thailand

Sartorius (Thailand) Co. Ltd.
129 Rama 9 Road,
Huaykwang
Bangkok 10310

Phone +66.2643.8361-6
Fax +66.2643.8367



◀ www.sartorius.com