



Picus[®] and
Picus[®] NxT
Electronic Pipettes

Simplifying Progress

SARTORIUS



Picus[®] and Picus[®] NxT

Pipetting simplified

The Picus[®] family pipettes, Picus[®] and Picus[®] NxT, with Sartorius pipette tips are the new standard for lab professionals.

The Picus family pipettes are kind to your hand with unbeatable ergonomic design that ensures reliable and repeatable experiment results. Picus is the right tool for everybody from intern to lab expert, being as easy to start with as a mechanical pipette and offering options for even the most advanced user. Picus saves you time in the lab with pipetting modes and customizable programs for every need. The pipette is easy to get comfortable with and quick to use. Just pick it up, switch it on and pipette!

Picus[®] NxT takes on where Picus[®] left with added features like user customizable protocols, password protection, and calibration reminders. Especially when conformity to the strictest regulations is needed by selecting the advanced Picus[®] NxT you can take your compliance to a new level with added certification and safety features.

Both Picus[®] and Picus[®] NxT, are available in single-channel models, covering a volume range of 0.2-10 000 µl and in multi-channel models from 0.2 µl to 1200 µl.

Choose the tools that any professional deserves.



Be kind to your hand

- Effortless pipetting through the uniquely lightweight (only 100 grams) and streamlined design
- Muscle strain minimized with conveniently located soft-touch operating button and electronic tip ejection
- Comfortable handle design and finger hook let each Picus[®] pipette rest lightly in your hand



Speed up your work

- Extremely quick volume setting and menu navigation using the unique adjustment wheel
- The extensive range of pipetting modes, from diluting and titrating to serial dispensing, makes pipetting quick and easy
- Intuitive user interface with a range of language options: English, French, German, Russian, and Chinese.
- With Picus[®] NxT user-defined protocols further speed up routine pipetting



Achieve reliable results

- Repeatable pipetting results are guaranteed with the electronic piston control and brake, raising all users to expert level
- Unique plate tracker increases reliability in microwell plate dispensing
- The option to use Safe-Cone Filters provide cost-effective contamination prevention
- With Picus[®] NxT the repeated blow-out function helps dispense every last droplet, ensuring complete sample recovery



Picus[®] NxT: Conform to Lab regulations

- Certificate of accredited 3-point calibration (ISO 17025 and ISO 8655) conforms to the strictest regulations
- Get alerts for maintenance and calibration reminders when service is required
- Advanced password protection for settings and pipetting protocol memory
- Possibility to lock the pipette, when its use is to be prevented, e.g. in case of contamination

Ergonomic Award Winning Design



reddot design award
winner 2012



Large easy-to-read display

All information of your active pipetting step available at a glance.



Soft Keys for Programming and a Hotkey

Fast and easy access to up to 10 of your most used pipetting settings by saving them with hotkey.



Electronic Tip Ejection

Protects you and your hand from Work-Related Upper Limb Disorder by removing the tips effortlessly with an electronic system.



Adjustment Wheel

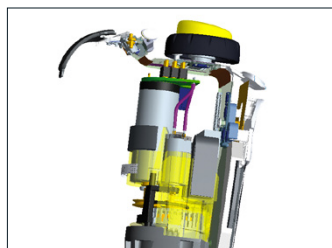
Offers extremely fast volume setting and easy menu navigation. Enables ergonomic one-handed operation. Accurately control manual pipetting and titrating speed with just a light touch of the thumb.

Reliable Pipetting



Microwell Plate Tracker

Never lose track of the next well on a 96 or 384 well plates with the unique well plate tracker functionality that improves work efficiency and reliability of your results. Tracker is an optional feature used in conjunction with the Pipetting, Reverse Pipetting and Multi Dispensing modes.



Advanced precision with sophisticated piston control

Achieve low dispensing variation with Sartorius pipette tips especially in serial dispensing with the optical sensor and the electronic brake precisely controlling the piston movement.



Optiload for Perfect Tip Sealing

Load tips with an equal constant force to every channel. Perfect tip sealing ensured to every individual tip cone when using Sartorius pipette tips. Optiload available in all multichannel versions.



Safe-Cone Filters and Autoclavable Lower Parts

Protect your samples and pipette from getting contaminated with exchangeable Safe-Cone Filters.

Picus® NxT: Regulatory compliance

Speed Up Your Work

Many workflows are designed to use multiple different pipetting modes for optimal results. The Picus® NxT pipette allows you to simplify your workflow by allowing you to create three custom pipetting protocols with up to 10 steps each. This reduces errors, speeds up the work and lets you to focus on pipetting by eliminating the need to change modes and volume settings back and forth in the midst of the workflow.

3-Point Calibration Certificate

It is essential to have pipettes calibrated in compliance with ISO 17025 for all laboratories that depend on the quality and reliability of results. Every Picus® NxT is ready for work directly from the box as they are delivered with a certificate of accredited 3-point calibration (ISO 17025 and ISO 8655).



Two-level Password Protection

Separate administrator and user profiles in Picus® NxT prevent unauthorized modifications to pipette settings and can be used to restrict access to specific individuals or groups. This feature ensures that your work is done exactly as the method, regulations or laboratory guideline dictates. Passwords are an optional feature and can be set as inactive when extra protection is not necessary.

Pipette Locking

Have you had sticky notes stating “Do not use, waiting for service!” attached to your old pipettes? With the pipette locking feature you can prevent the pipette from being used when it is awaiting service or cleaning. If password protection is enabled, the administrator password is required to unlock the pipette.

Service and Calibration Reminders

Your pipettes as precision measuring instruments must be serviced and calibrated regularly, according to GLP guidelines and ISO standards. The Picus® NxT can be programmed with service and calibration reminders according to time and/or cycle intervals. Maintaining and calibrating your pipettes regularly is critical to ensure the quality and accuracy of your results.

Choosing between Picus® and Picus® NxT

Picus is the choice when you want to upgrade your pipetting and leave the variables of manual pipetting behind and pipette with the touch of a button. When you wish to optimize your workflow and regulatory compliance, select Picus® NxT and take advantage of the user programmable protocols and safety features. In the table below the pipette functions are introduced with Picus® NxT specific ones highlighted in yellow.

Pipetting Modes	Advanced Functions (ADV)*						
	Plate Tracker	Mixing	Counter	Excess Volume Adjustment	Auto-Dispensing	Fast Dispensing	Repeated Blow-out
Pipetting	■	■	■				■
Reverse Pipetting	■		■	■			
Manual Pipetting							■
Multi-Dispensing	■			■	■		
Diluting		■					■
Sequential Dispensing				■			
Multi-Aspiration							■
Titration						■	
Protocols**		■		■	■	■	■

*Advanced functions are used in conjunction with the pipetting modes. **The availability of advanced functions varies according to the pipetting mode selected. Highlighted available in Picus® NxT models.

Remember also these Picus family features

Quick charge battery lasts for the whole day.

Fully recharge your pipette in an hour.

Plug-and-pipette

Starting to use a Picus is easy thanks to the intuitive user interface. The pipettes are ready to use directly from the package.

Liquid Adjustment Feature

Enables maximal repeatability even with difficult liquids. Combine with Low Retention tips for optimal performance.

Picus® Charging Options

The electronic pipettes can be charged with a single pipette charging stand, a carousel stand for 4 pipettes, or with a USB cable that you can also use while pipetting. The compact design of the carousel stand is ideal for saving bench space in the laboratory. It features a rotating head that provides easy access to all of your pipettes.

















Sartorius Optifit and SafetySpace™ Filter tips

To achieve the performance specifications of Picus® and Picus® NxT choose Sartorius pipette tips for your lab. Also remember to calibrate your pipette using the same pipette tips you are working with.

Pipette Stands

Order Code	Item
730981	Charging Stand for one pipette
730991	Charging Carousel for 4 pipettes
725620	Linear Stand (non-charging)

Ordering Information: Picus® and Picus® NxT

Picus® NxT Order Codes	Picus® Order Codes	Channels	Volume Range (µl)	Safe-Cone Standard	Filters Plus	Optifit Tips* (µl)	SafetySpace Filter Tips* (µl)
LH-745021	735021	1 	0.2 - 10	-	-	10	10
LH-745041	735041	1 	5 - 120	721008	721018	200	120
LH-745061	735061	1 	10 - 300	721007	721017	350	300
LH-745081	735081	1 	50 - 1,000	721006	721016	1,000	1,000
LH-745101	735101	1 	100 - 5,000	721005	721015	5,000	5,000
LH-745111	735111	1 	500 - 10,000	721005	721015	10,000	-
LH-745321	735321	8 	0.2 - 10	-	-	10	10
LH-745341	735341	8 	5 - 120	721008	721018	200	120
LH-745361	735361	8 	10 - 300	721007	721017	350	300
LH-745391	735391	8 	50 - 1,200	721006	721016	1,200	1,200
LH-745421	735421	12 	0.2 - 10	-	-	10	10
LH-745441	735441	12 	5 - 120	721008	721018	200	120
LH-745461	735461	12 	10 - 300	721007	721017	350	300
LH-745491	735491	12 	50 - 1,200	721006	721016	1,200	1,200

* Low Retention Tips are available in volumes up to 1,200 µl.

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 sartorius.com

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-Free +1 800 368 7178

Finland & Baltics

Sartorius
Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki
Phone +358 9 755 951