



Tacta[®] Mechanical
Pipettes
Perfectly Balanced

Simplifying Progress

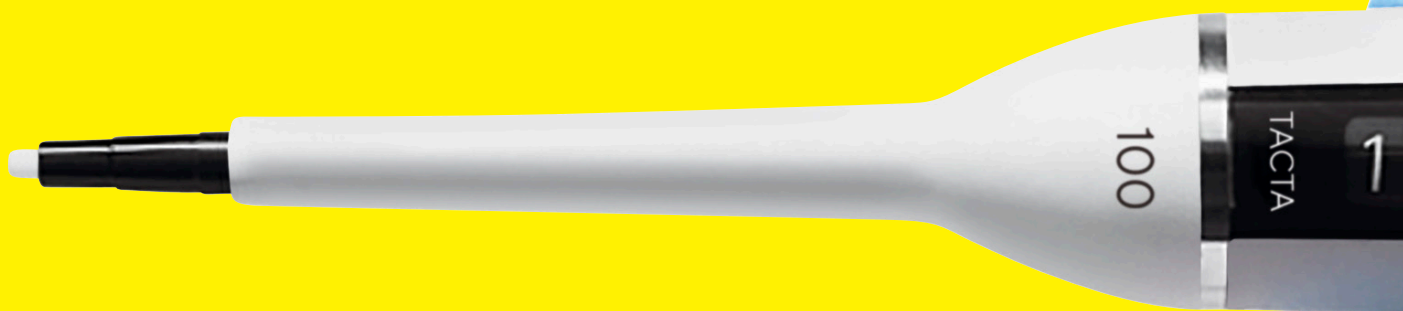
SARTORIUS

Tacta[®] Mechanical Pipettes

Have you ever considered the overall user experience in pipetting? We have.

Sartorius proudly presents Tacta[®], the premium mechanical pipette with superb comfort and reliability. Tacta[®] makes pipetting effortless and safe, while producing accurate and reliable results time after time.

At Sartorius, we listened to our customers and combined their views with our extensive experience, and state-of-the-art R&D, to create the exceptionally balanced design of Tacta[®]. Tacta[®] is designed to feel great in your hand, and it is easy and comfortable to use. All materials used for Tacta[®] have been carefully selected, and each component is designed to meet the highest standards.





The Tacta[®] Experience - Perfectly Balanced

- Comfortable to hold due to the ergonomically designed handle
- Low pipetting and tip ejection forces that reduce the risk of strain injury
- The unique Sartorius Optilock system provides flexibility for volume adjustment and locking
- Controlled and smooth tip ejection with the new Sartorius Optiject technology
- Calibration adjustment to provide accurate results for various liquid types
- Easy to clean, with only three parts to disassemble
- Reliable results, even over lengthy pipetting periods





SARTORIUS

200.0

TACTA

200

Feel the Comfort

Superior Ergonomics

Tacta® rests lightly in your hand thanks to its ergonomically designed handle and finger hook. Tacta® is effortless to use and exceptionally comfortable to hold. Tacta® eases your workload and protects you from strain, even when you pipette for extended periods of your working day. The unique handle and finger hook design lets the pipette rest lightly in your hand, with no need to grip the handle tightly.



🌐 For more information, please visit
www.sartorius.com/en/products/pipetting-and-dispensing/mechanical-pipettes/tacta

Volume Adjustment

Use the Sartorius Optilock feature to prevent accidental volume changes during pipetting. Either hold down the volume lock button with one hand and adjust the volume with the other, then release to lock, or with just one hand, slide the volume lock up, adjust the volume and slide down to lock again. The choice is in your hands.



Optilock

Large, Clear Display

The large, easy-to-read display helps you to see all four digits of the set volume from a distance without straining your eyes. The volume is easy to read even when the pipette is angled – eliminating the need to turn your head into an uncomfortable position.



The volume is easy to read even when the pipette is angled.

Calibration Adjustment

Tacta® is very easy to adjust, for various types of liquids, using a simple adjustment key. The adjustment scale shows the degree of adjustment. By noting this value for a specific liquid, you can return to that setting any time. Easy calibration adjustment ensures accurate results with a variety of liquids.



Calibration Adjustment

Low Pipetting, Tip Loading and Ejection Forces

The forces needed for a pipetting cycle, from tip attachment to pipetting and finally tip ejection, are exceptionally low with Tacta®, protecting you from strain injury.

The Optiload feature, with spring-loaded tip cones in both the single and multichannel models, secures tip loading with perfect sealing

and minimal force. Consequently, Optiload also ensures light tip ejection to protect your hand.

The light tip ejection is further enhanced by Sartorius Optiject, the soft, levered tip ejection feature that lets the tip detach in a smooth, controlled way with little force.

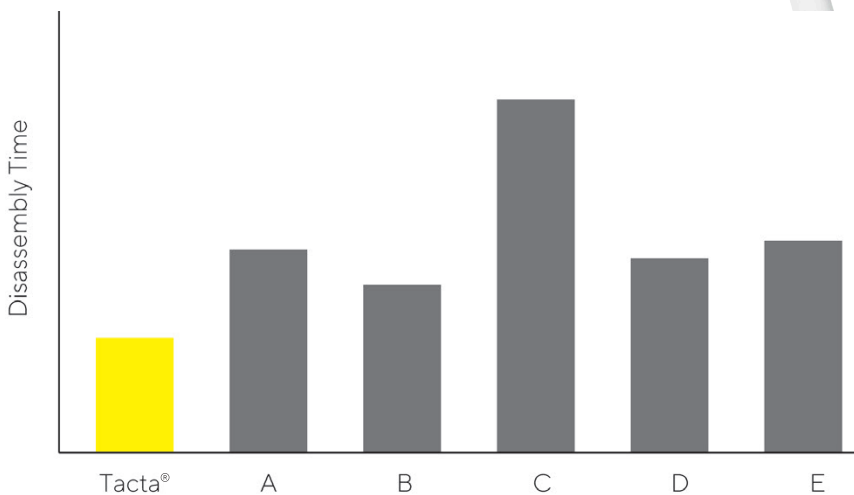


Optiload

Safe from Contamination

Easy Cleaning

Tacta® is the quickest and easiest pipette, on the market, to clean, with only three parts to be disassembled for cleaning and no tools are needed for disassembly. Tacta® can also be steam-sterilized or autoclaved as-is, without disassembly. It has also high UV and chemical resistance.



The clever design of the Tacta® requires only three parts to be disassembled for cleaning, making the Tacta®'s disassembly time significantly shorter than its competitors'. This makes pipette cleaning an easy and convenient task for anyone and simplifies routine maintenance. Regular cleaning will ensure reliable results and prolong the life of the pipette.

The test was carried out using 1000 µl pipettes.



Safe-Cone Filters

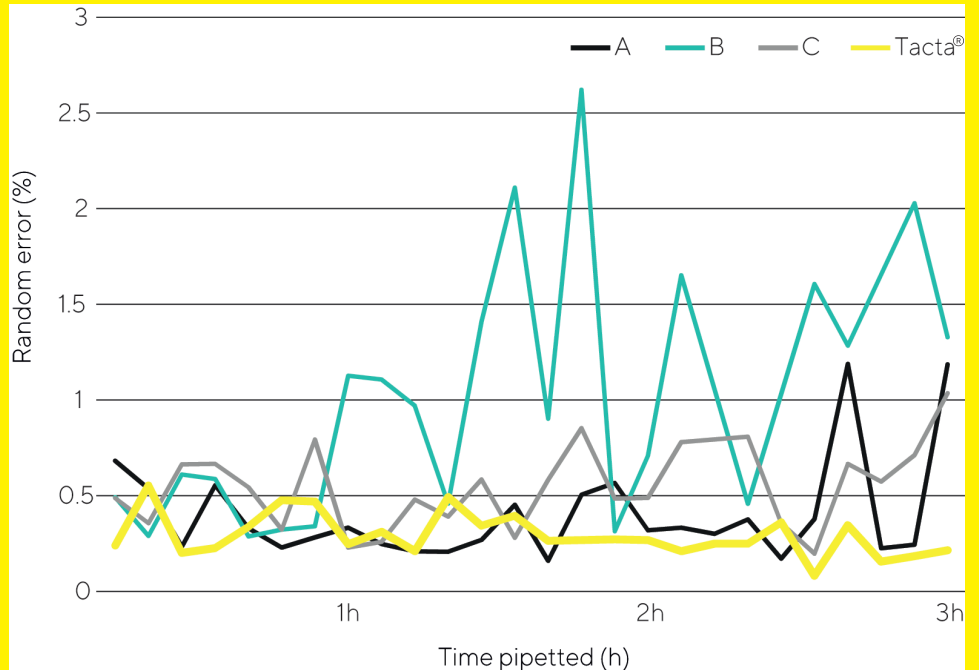
The exchangeable Safe-Cone Filters, used in pipette tip cones, act as barriers to prevent sample aerosols and fluids from contaminating the internal components of the pipette. The unique Safe-Cone Filter ejector enables the removal of used filters without human contact with contaminated filters.

Safe-Cone Filters are available for all Tacta® models greater than 10 µl and offer a cost-effective method to reduce contamination. Filters should be changed regularly, and always after over-aspiration.

Achieve Reliable Results

Proven Precision

The following graph illustrates how the precision of pipetting results is impacted over time when using various pipettes. Tacta[®], which is designed down to the last detail to create a comfortable pipetting experience, produced consistently precise results during a three-hour test period.



Precision of results tested over a three-hour period using four pipettes from various manufacturers. The test was performed by an independent laboratory in accordance with the ISO 8655 standard using 1000 µl pipettes with 1000 µl tips from the corresponding manufacturers.

Sartorius Pipette Tips and Services

Complete your Tacta[®] experience with original Sartorius pipette tips. Our Optifit Tips and Safetyspace[®] Filter Tips are designed to fit your Tacta[®] pipette perfectly, enabling maximum accuracy and precision.

Sartorius also provides pipette maintenance, repair and calibration services globally. Contact your nearest Sartorius service centre or office for services.



Ordering Information

Tacta®

Order Code	Channels	Color Code	Volume Range (µl)	Safe-Cone Filters	Standard Plus	Optifit Tips* (µl)	Safetyspace® Filter Tips* (µl)
LH-729010	1	■	0.3-3	-	-	10	10
LH-729020	1	■	1-10	-	-	10	10
LH-729030	1	■	2-20	721014	-	200	20
LH-729050	1	■	10-100	721008	721018	200	120
LH-729060	1	■	20-200	721007	721017	200	200
LH-729070	1	■	100-1,000	721006	721016	1,000	1,000
LH-729080	1	■	500-5,000	721005	721015	5,000	5,000
LH-729090	1	■	1,000-10,000	721005	721015	10,000	-
LH-729120	8	■	1-10	-	-	10	10
LH-729130	8	■	10-100	721008	721018	200	120
LH-729140	8	■	30-300	721007	721017	350	300
LH-729220	12	■	1-10	-	-	10	10
LH-729230	12	■	10-100	721008	721018	200	120
LH-729240	12	■	30-300	721007	721017	350	300

* Tips are available in various packaging formats. Please contact your Sartorius representative or visit our websites for further information.

Storing the Pipettes

When Tacta® is not in use, it should be stored in an upright position in order to avoid contamination from work surfaces. The compact Carousel Stand, designed for six pipettes, is ideal for saving bench space.

The Linear Stand is designed for all Sartorius mechanical and electronic pipettes. The simplest solution of all is to attach pipette holders to the front edge of a shelf. All Tacta®s are supplied with a pipette holder.

Pipette Stands

Order Code	Item
LH-725630	Carousel Stand for 6 mechanical pipettes
725620	Linear Stand
LH-727640	Pipette Holder



🌐 For more information, please visit www.sartorius.com/en/products/pipetting-and-dispensing/mechanical-pipettes/tacta

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 Liquid Handling Oy
Tulppatie 1
00880 Helsinki
Email: lhinfo.finland1@sartorius.com

 For further contacts, visit
www.sartorius.com