

Tacta®

Mechanical Pipettes That Fits in Your Hand and Your Laboratory

Product Information

Tacta® is a premium mechanical pipette featuring superb comfort and reliability. Designed with comfort in mind, Tacta® makes pipetting effortless and safe, while producing accurate and reliable results time after time. Tacta® is made from carefully selected materials, with each component designed to meet the highest standards.

As Tacta® was designed with comfort in mind, it fits both in your hand and in your laboratory. As it requires less force than other mechanical pipettes on the market, it can be used for longer pipetting sessions without causing repetitive strain injuries, or affecting the reproducibility and reliability of the results.

Description

Tacta® is available in both single and multichannel versions in different volume ranges, 0.1 to 10,000 µL and 0.5 to 300 µL respectively. This to ensure that there is a Tacta® for your pipetting needs in your laboratory.

As a pipette, the Tacta® is comfortable to hold and use thanks to its ergonomic design, including low weight, and finger support and having the plunger, volume adjustment, volume lock, and tip ejector all close to your thumb. The Tacta® fits in your left and right hand, and can be operated with only one hand, including changing the volume.



Features

- Low pipetting cycle forces reduce the risk of strain injury
- Ergonomic design for comfortable pipetting
- Optiload feature in both single and multichannel models for easy and light tip loading with perfect tip sealing
- Optiject soft tip ejection feature detaches the tip in a smooth, controlled manner
- Optilock's dual-function volume locking prevents accidental volume changes
- Easy-to-read, four-digit volume display helps to set exact volumes
- Color-coding of volumes to ease the selection of corresponding pipette tips
- Safe-Cone Filters available for models above 10 µL with convenient built-in filter ejection mechanism

Easy Maintenance

- Fully autoclavable without disassembling
- Simple to clean and maintain with only three parts to disassemble
- Easy calibration adjustment with adjustment scale ensures accurate results, also in cases where factory calibration does not apply, for example when pipetting viscous liquids
- Materials have high chemical and UV-resistance to ensure a long life span for the pipette

Intended Use

Manual liquid-handling of volumes ranging from 0.1 µL to 10,000 µL.



Technical Data

Technical Specifications

Weight	79 g (1-ch, 1,000 µL) 138 g (8-ch, 300 µL)	Stand options	Carousel Stand for 6 pipettes Linear Stand Pipette Holder for one pipette
Length	225 mm (1-ch, 1,000 µL) 240 mm (8-ch, 300 µL)	Safe-Cone Filter ejector	Built-in ejector
Autoclavability	Fully autoclavable (121 °C, 20 min, 1 bar/100 kPa overpressure)	Limitations	The accuracy and precision values are valid under the controlled conditions described in ISO 8655. Test temperature: +15 °C to +30 °C +/-0.5 °C Relative humidity: > 50 %
UV-resistance	UV resistance: Tacta® pipettes are made of UV-resistant materials. Sartorius pipettes tolerate temporary exposure to UV radiation. Take note that prolonged or frequent exposure to UV radiation may cause yellowing and brittling of the pipette.	Instrument use outside +20 °C to +25 °C	When using instruments outside this temperature range, the correction factor for water density (Z-value µL/mg) must be used.
Pipetting principle	Air displacement	Transportation conditions	Temperature: -50 °C to +50 °C Humidity: 10–90% Shocks: <10 G
Pipetting modes	Pipetting (P) Reverse pipetting (rP) Post-delivery mixing	Long-term storage conditions	Temperature: -20 °C to +40 °C Humidity: 10–60 %
		Warranty	2 years

Volume Adjustment

The pipette calibration should be adjusted when necessary, for example when reverse pipetting, when pipetting liquids other than water, when using the pipette at high altitudes, or when the geometry of the tip you use clearly differs from standard tip geometry.

Model [µL]	Channels	Increment [µL]	-45	-10	-5	-1	0	1	5	10	45
0.1–3	1	0.002	-0.09	-0.02	-0.01	-0.002	0	0.002	0.01	0.02	0.09
0.5–10	1	0.01	-0.45	-0.1	-0.05	-0.01	0	0.01	0.05	0.1	0.45
2–20	1	0.02	-0.9	-0.2	-0.1	-0.02	0	0.02	0.1	0.2	0.9
10–100	1	0.1	-4.5	-1	-0.5	-0.1	0	0.1	0.5	1	4.5
20–200	1	0.2	-9	-2	-1	-0.2	0	0.2	1	2	9
100–1,000	1	1	-45	-10	-5	-1	0	1	5	10	45
500–5,000	1	5	-225	-50	-25	-5	0	5	25	50	225
1,000–10,000	1	10	-450	-100	-50	-10	0	10	50	100	450
0.5–10	8 & 12	0.01	-0.45	-0.1	-0.05	-0.01	0	0.01	0.05	0.1	0.45
10–100	8 & 12	0.1	-4.5	-1	-0.5	-0.1	0	0.1	0.5	1	4.5
30–300	8 & 12	0.2	-9	-2	-1	-0.2	0	0.2	1	2	9

Ordering Information

Tacta [®] Order Code	Channels		Volume Range (μL)	Increment (μL)	Test Volume (μL)	Systematic Error ^N Limit ± (%)	Random Error ^N Limit (μL)		
LH-729010	1		0.1–3*	0.002	3	1.6	0.048	0.8	0.024
					1.5	3.0	0.045	1.6	0.024
					0.3	12.0	0.036	6.0	0.018
LH-729020	1		0.5–10*	0.01	10	1.0	0.10	0.6	0.06
					5	1.5	0.075	1.0	0.05
					1	3.0	0.03	2.0	0.02
LH-729030	1		2–20	0.02	20	1.0	0.2	0.4	0.08
					10	1.5	0.15	0.7	0.07
					2	6.0	0.12	3.0	0.06
LH-729050	1		10–100	0.10	100	0.8	0.8	0.2	0.2
					50	1.0	0.5	0.3	0.15
					10	3.0	0.3	1.0	0.1
LH-729060	1		20–200	0.20	200	0.6	1.2	0.2	0.4
					100	1.0	1	0.3	0.3
					20	2.5	0.5	0.9	0.18
LH-729070	1		100–1,000	1.00	1,000	0.7	7	0.2	2
					500	0.9	4.5	0.2	1
					100	3.0	3	0.6	0.6
LH-729080	1		500–5,000	5.00	5,000	0.6	30	0.2	10
					2,500	0.7	17.5	0.25	6.25
					500	2.4	12	0.6	3
LH-729090	1		1,000–10,000	10.00	10,000	0.6	60	0.2	20
					5,000	1.2	60	0.3	15
					1,000	3.0	30	0.6	6
LH-729120	8		0.5–10*	0.01	10	1.5	0.15	1.0	0.1
LH-729220	12				5	2.5	0.125	2.0	0.1
					1	5.5	0.055	4.0	0.04
LH-729130	8		10–100	0.10	100	0.9	0.9	0.4	0.4
LH-729230	12				50	1.2	0.6	0.6	0.3
					10	4.0	0.4	2.0	0.2
LH-729140	8		30–300	0.20	300	0.6	1.8	0.25	0.75
LH-729240	12				150	1.0	1.5	0.4	0.6
					30	3.0	0.9	1.0	0.3

Note: The values listed for the systematic and random errors are valid for listed pipettes used in combination with suitable Sartorius Optifit non-sterile tips and can only be achieved under strictly controlled conditions during type examinations in accordance with ISO 8655. The listed pipettes with other Sartorius tips fulfill specifications listed in ISO 8655.

*The ISO 8655 does not recognize using the pipette below 10% of its nominal (maximum) volume and therefore it is not recommended although the pipette's volume can be adjusted below 10% of nominal volume.

Germany

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

USA

Sartorius Corporation
3874 Research Park Drive
Ann Arbor, MI 48108
Phone +1 734 769 1600

 For further contacts, visit
www.sartorius.com