

Tacta[®]

Mechanical Pipettes

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.



Description

Tacta[®] is a premium mechanical pipette available in volume range from 0.1 to 10,000 μL in single channel models, and from 0.5 μL to 300 μL in multichannel models.

Features

- Low pipetting cycle forces reduce the risk of WRULD (Work Related Upper Limb Disorder)
- 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
- 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 (included in the pipette package)
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 calibration of Tacta® is based on ISO 8655-6: gravimetric test method for volumetric instruments. It has been factory-checked using the forward-pipetting technique and certified according to ISO 3696 at 22°C using grade-3 distilled water.

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
5-100	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 (%)	Systematic Error ^N Limit ± (µL)	Random Error ^N Limit (µL)
LH-729010	1	0.1–3	0.002	3	1.4	0.8	0.042	0.024
				1.5	2.6	1.6	0.039	0.024
				0.3	10.0	6.0	0.030	0.018
LH-729020	1	0.5–10	0.01	10	1.0	0.6	0.100	0.060
				5	1.5	1.0	0.075	0.050
				1	3.0	2.0	0.030	0.020
LH-729030	1	2–20	0.02	20	1.0	0.5	0.200	0.100
				10	1.4	0.9	0.140	0.090
				2	4.0	3.0	0.080	0.060
LH-729050	1	10–100	0.10	100	0.8	0.2	0.80	0.20
				50	1.0	0.4	0.50	0.20
				10	3.0	1.0	0.30	0.10
LH-729060	1	20–200	0.20	200	0.6	0.2	1.20	0.40
				100	0.8	0.3	0.80	0.30
				20	2.3	0.9	0.46	0.18
LH-729070	1	100–1,000	1.00	1,000	0.7	0.2	7.0	2.0
				500	0.8	0.2	4.0	1.0
				100	2.5	0.6	2.5	0.6
LH-729080	1	500–5,000	10.0	5,000	0.6	0.2	30	10
				2,500	0.7	0.3	17.5	7.5
				500	2.4	0.6	12	3
LH-729090	1	1,000–10,000	20.0	10,000	0.6	0.2	60	20
				5,000	1.2	0.3	60	15
				1,000	3.0	0.6	30	6
LH-729120	8	0.5–10	0.01	10	1.5	1.0	0.150	0.100
LH-729220	12			5	2.5	2.0	0.125	0.100
				1	5.5	4.0	0.055	0.040
LH-729130	8	5–100	0.10	100	0.9	0.4	0.90	0.40
LH-729230	12			50	1.2	0.7	0.60	0.35
				10	4.0	2.0	0.40	0.20
LH-729140	8	30–300	0.20	300	0.6	0.25	1.80	0.75
LH-729240	12			150	1.0	0.5	1.50	0.75
				30	2.5	1.0	0.75	0.30

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.

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 contacts, visit
www.sartorius.com