



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

- Colour-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 ml.



Technical Data

Technical Specifications

Weight	79 g (1-ch, 1000 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, 1000 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)	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.

For accurate adjustment results, please refer to the table below.

Volume Adjustment

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
100-1000	1	1	-45	-10	-5	-1	0	1	5	10	45
500-5000	1	5	-225	-50	-25	-5	0	5	25	50	225
1000-10000	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	
					(%)	(µl)	(%)	(µl)
LH-729010	1	0.1 – 3	0.002	3	1.4	0.042	0.8	0.024
				1.5	2.6	0.039	1.6	0.024
				0.3	10.0	0.030	6.0	0.018
LH-729020	1	0.5 – 10	0.01	10	1.0	0.100	0.6	0.060
				5	1.5	0.075	1.0	0.050
				1	3.0	0.030	2.0	0.020
LH-729030	1	2 – 20	0.02	20	1.0	0.200	0.5	0.100
				10	1.4	0.140	0.9	0.090
				2	4.0	0.080	3.0	0.060
LH-729050	1	10 – 100	0.10	100	0.8	0.80	0.2	0.20
				50	1.0	0.50	0.4	0.20
				10	3.0	0.30	1.0	0.10
LH-729060	1	20 – 200	0.20	200	0.6	1.20	0.2	0.40
				100	0.8	0.80	0.3	0.30
				20	2.3	0.46	0.9	0.18
LH-729070	1	100 – 1,000	1.00	1,000	0.7	7.0	0.2	2.0
				500	0.8	4.0	0.2	1.0
				100	2.5	2.5	0.6	0.6
LH-729080	1	500 – 5,000	10.0	5,000	0.6	30	0.2	10
				2,500	0.7	17.5	0.3	7.5
				500	2.4	12	0.6	3
LH-729090	1	1,000 – 10,000	20.0	1,000	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	0.150	1.0	0.100
LH-729220	12			5	2.5	0.125	2.0	0.100
				1	5.5	0.055	4.0	0.040
LH-729130	8	5 – 100	0.10	100	0.9	0.90	0.4	0.40
LH-729230	12			50	1.2	0.60	0.7	0.35
				10	4.0	0.40	2.0	0.20
LH-729140	8	30 – 300	0.20	300	0.6	1.80	0.25	0.75
LH-729240	12			150	1.0	1.50	0.5	0.75
				30	2.5	0.75	1.0	0.30

N Note: The listed systematic and random error values can be achieved only under strictly controlled conditions during type tests per ISO 8655. Due to the continuous product development by Sartorius, the systematic and random error values are subject to change without prior notice.

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