

Biotrate

Digital Burette

Product Information

Sartorius Biotrate is a premium digital burette equipped with sophisticated functions for easy titration. Biotrate offers high chemical resistance and it is a smart solution for titrating various liquids accurately and safely.



Description

Biotrate is a premium digital burette for convenient titration.

Features

- Extremely high chemical resistance
- Smooth and light to operate with soft touch operating wheels
- Media recirculation system that allows unused media to be recovered from the device and returned to the bottle
- 360° rotating dispensing head
- Universal connector at the back of the unit for drying tubes and filters
- Easy to attach aspiration tube with a secure threaded connection
- User replaceable valves
- Supplied with UV-protected replacement window for light sensitive media protection
- Long life CR2032 battery
- Easy custom adjustment for various liquids
- Wide range of accessories
- Autoclavable liquid pathway

Intended Use

Manual titration of liquids.

Technical Data

Technical Specifications	
Weight	844 g (20 ml)
Length	260 mm (20 ml)
Autoclavability	Autoclavable liquid pathway (121°C, 20 min, 1 bar/100 kPa)
Threaded bottle fitting	A45
Bottle adapters included	A32, A38, S40
Battery	CR2032
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%
Transportation conditions	Temperature: 0°C to +40°C

Long-term storage conditions	Temperature: 0°C to +40°C
Chemical resistance	Information available on request
Materials	Piston: Compound Halar (ECTFE) Cylinder: Borosilicate glass 3.3 Valve block: Halar (ECTFE) Valve mantle: Halar (ECTFE) Valve ball: Borosilicate glass 3.3 Valve seat: Borosilicate glass 3.3 Valve spring: Hastelloy Tubes: FEP Sealing rings: PTFE
Warranty	2 years

Ordering Information

Order Code	Nominal Volume [ml]	Readability [ml]	Test Volume [ml]	Systematic Error* [%]	Systematic Error* [ml]	Random Error* [%]	Random Error* [ml]
Biotrate							
LH-723080	10	0,01	10	0,3	0,03	0,1	0,01
LH-723081	20	0,01	20	0,2	0,04	0,1	0,02
LH-723082	50	0,01	50	0,2	0,10	0,1	0,05

* The listed systematic and random error values can be achieved under strictly controlled conditions during type tests as per ISO 8655. Due to our continuous product development, the systematic and random error values may change without prior notice.



Sartorius Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki, Finland
Phone +358 9 755 951
Fax +358 9 755 95 200

Sartorius Lab Instruments
GmbH & Co. KG
Weender Landstrasse 94-108
37075 Goettingen, Germany
Phone +49 551 3080
Fax +49 551 308 3289

www.sartorius.com