

Prospenser and Prospenser Plus – Superior Bottle-top Dispensers



Prospenser & Prospenser Plus – Dispensing Made Easy

Prospenser and Prospenser Plus are high quality bottle-top dispensers with excellent chemical resistance. They are safe, convenient, and customizable, making them perfect for a wide range of dispensing needs.

Superior Chemical Resistance

- Both the ceramic piston and all parts in the liquid pathway are highly chemical resistant, guaranteeing versatility for a wide range of applications.
- High quality materials guarantee a long lifetime.

High Accuracy & Precision

- The incremental volume setup ensures accurate and repeatable results.
- High quality parts and assembly ensure smooth piston movement.

Easy to Use

- Easy and comfortable dispensing for various applications.
- Smooth and light to operate.



Prospenser & Prospenser Plus – Features

Chemical Resistance

Prospenser and Prospenser Plus offer excellent chemical resistance thanks to their high quality materials. The ceramic piston is chemically stable and compatible with an extremely wide range of liquids.

Light to Use

The operating handle is designed to be comfortable to hold while aspirating and light to press while dispensing. The piston moves smoothly, guaranteeing excellent results in everyday use.

Easy to Attach Aspiration Tube

The aspiration tube is easy to attach with a secure threaded connection. This prevents air entering the system and ensures consistent results.

Easily Replaceable Valves

Valves can be easily replaced by the user so there is no need to send the unit to be serviced.

Volume Setting

Prospenser and Prospenser Plus have an incremental volume setting enabling you to set the exact volume every time.

Versatile Accessories

There is a wide range of accessories available including bottle adapters, dispensing heads, and aspiration tubes.

Fully Autoclavable

Prospenser and Prospenser Plus are easy to disassemble for cleaning and are fully autoclavable.



Prospenser Plus – Advanced Features



Media Recirculation System

Prospenser Plus is equipped with a media recirculation system that allows unused media to be recovered from the device and returned to the bottle. This ensures easy priming and minimizes media loss.



Smart 360° Rotating Dispensing Head

The dispensing head can be turned to the required position without needing to turn the bottle. This ensures the bottle label is always visible to the user.



Connector for Drying Tube and Filters

Prospenser Plus has a luer connection at the back of the unit for drying tubes and filters. Drying tubes are recommended for humidity sensitive liquids.





Prospenser and Prospenser Plus – Technical Details

Features	Prospenser	Prospenser Plus
Volume ranges	0,2–1 ml 0,4–2 ml 1–5 ml 2–10 ml 5–30 ml 10–60 ml	0,2–1 ml 0,4–2 ml 1–5 ml 2–10 ml 5–30 ml 10–60 ml
360° rotating dispensing head	No	Yes
Media recirculation system	No	Yes
Connector for drying tube and filters	No	Yes
Threaded bottle fitting	A32	A45
Bottle adapters included	A28, S40, A45	A32, A38, S40
Extremely high chemical and thermal resistance	Ceramic piston, glass cylinder, and highly resistant fluoroplastic parts	Ceramic piston, glass cylinder, and highly resistant fluoroplastic parts
Repeatable volume setting	Yes	Yes
Operating force	Low	Low
Aspiration tube fitting	Threaded	Threaded
Delivered with quality certificate	Yes	Yes
Easily replaceable valves	Yes	Yes
Easy to dismantle for cleaning	Yes	Yes
Fully autoclavable	121°C, 1 bar	121°C, 1 bar

Ordering Information

Bottle-top Dispensers

Prospenser

Order Code	Item	Volume Range
LH-723060	Prospenser	0,2–1 ml
LH-723061	Prospenser	0,4–2 ml
LH-723062	Prospenser	1–5 ml
LH-723063	Prospenser	2–10 ml
LH-723064	Prospenser	5–30 ml
LH-723065	Prospenser	10–60 ml

Prospenser Plus

Order Code	Item	Volume Range
LH-723070	Prospenser Plus	0,2–1 ml
LH-723071	Prospenser Plus	0,4–2 ml
LH-723072	Prospenser Plus	1–5 ml
LH-723073	Prospenser Plus	2–10 ml
LH-723074	Prospenser Plus	5–30 ml
LH-723075	Prospenser Plus	10–60 ml

Note: Bottle is not included



Accessories

Order Code	Item	Material	Prospenser Plus	Prospenser	Prospenser
			All models	1, 2, 5 & 10 ml	30 & 60 ml
LH-721647	Dispensing head, standard	FEP	•		
LH-721655	Dispensing head, standard	FEP		•	
LH-721656	Dispensing head, standard	FEP			•
LH-721648	Dispensing head, 1,5 mm spiral	FEP	•		
LH-721649	Dispensing head, 3,0 mm spiral	FEP	•		
LH-721657	Dispensing head, spiral	FEP		•	
LH-721658	Dispensing head, spiral	FEP			•
LH-721653	Dispensing head, luer lock	FEP/PP	•		
LH-721654	Dispensing head, luer lock	FEP/PFA	•		
LH-721659	Dispensing head, luer lock	FEP/PP			•
LH-721660	Dispensing head, luer lock	FEP/PFA			•
LH-721678	Aspiration tube, 310 mm	FEP	•		•
LH-721680	Aspiration tube, 310 mm	FEP		•	
LH-721679	Aspiration tube, 400 mm	FEP	•		•
LH-721681	Aspiration tube, 400 mm	FEP		•	
LH-721682	Recirculation tube, 70 mm	FEP	•		
LH-721736	Bottle adapter, A25	PP		•	•
LH-721737	Bottle adapter, A28	PP		•	•
LH-721689	Bottle adapter, A28	ETFE	•		
LH-721684	Bottle adapter, A28	PP	•		
LH-721683	Bottle adapter, A32	PP	•		
LH-721688	Bottle adapter, A32	ETFE	•		
LH-721686	Bottle adapter, A38	PP	•		
LH-721733	Bottle adapter, A38	ETFE	•		
LH-721738	Bottle adapter, A38	PP		•	•
LH-721685	Bottle adapter, A38/430	PP	•		
LH-721732	Bottle adapter, A38/430	ETFE	•		
LH-721741	Bottle adapter, A45	ETFE		•	•
LH-721739	Bottle adapter, A45	PP		•	•
LH-721687	Bottle adapter, S40	PP	•		
LH-721734	Bottle adapter, S40	ETFE	•		
LH-721740	Bottle adapter, S40	PP		•	•
LH-721742	Bottle adapter, S40	ETFE		•	•
LH-721735	Bottle adapter, NS29/32	Silicone	•		
LH-721743	Bottle support, for bottles 75 - 120 mm	Silicone	•	•	•

Sartorius Minisart® Syringe Filters

Sartorius Minisart® filters fit Prospenser and Prospenser Plus dispensers perfectly. Minisart® filters offer superior flow rate, high total throughput volumes, and exceptional quality.



Prospenser and Prospenser Plus can be equipped with Minisart® filters for media filtration. Dispensing heads with luer lock connection are available separately.



For filtering replacement air, Prospenser Plus can be equipped with a Minisart® filter.

Sartorius Pipettes, Pipette Tips, and Services

Our leading range of superior electronic and mechanical pipettes are the perfect solution for your liquid handling needs. Sartorius pipettes and tips fit seamlessly together maximizing pipetting performance and ensuring accuracy and precision. For more information, visit www.sartorius.com.

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



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