# SARTURIUS

# Prospenser and Prospenser Plus Bottle-top Dispensers



#### Product Information

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.

# Description | Intended Use

Prospenser and Prospenser Plus are premium bottle-top dispensers available in a volume range from 0.2 to 60 mL, used for manual dispensing of liquids.

### **Features**

- Extremely high chemical and thermal resistance due to a ceramic piston, glass cylinder, and highly resistant fluoroplastic parts
- The incremental volume setting ensures accurate and repeatable results
- Smooth and light to operate
- Easy to attach aspiration tube with a secure threaded connection
- User-replaceable valves
- Wide range of accessories
- Fully autoclavable

# Additional Prospenser Plus Features

- 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

## Technical Data

Weight	200 g (5 mL, Prospenser)   388 g (5 mL, Prospenser Plus)				
Length	143 mm (5 mL, Prospenser)   175 mm (5 mL, Prospenser Plus)				
Autoclavability	Fully autoclavable (121 °C, 20 min, 1 bar/100 kPa)				
Threaded bottle fitting	Prospenser: A32   Prospenser Plus: A45				
Bottle adapters included	Prospenser: A28, S40, A45   Prospenser Plus: A32, A38, S40				
Limitations	The accuracy and precision values are valid under the controlled conditions described in ISO 8655. Test temperature: $15 ^{\circ}\text{C}$ to $30 ^{\circ}\text{C}$ +/-0.5 $^{\circ}\text{C}$ . 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: Highly pure Al2O3-ceramic (99.7%)   Cylinder: Borosilicate glass 3.3   Valve block: Halar (ECTFE) Valve ball: Borosilicate glass 3.3   Valve seat: Borosilicate glass 3.3   Valve spring: Hastelloy (99.7%)				
Warranty	2 years				

# Ordering Information

Propenser	Propenser Plus	Volume Range (mL)	Increment (mL)	Test Volume (mL)	Systematic Error* (%)	Systematic Error* (mL)	Random Error* (%)	Random Error* (mL)
LH-723060	LH-723070	0.2-1	0.05	1	0.6	0.006	0.2	0.002
LH-723061	LH-723071	0.4-2	0.05	2	0.6	0.012	0.2	0.004
LH-723062	LH-723072	1-5	0.10	5	0.6	0.03	0.2	0.01
LH-723063	LH-723073	2-10	0.25	10	0.6	0.06	0.2	0.02
LH-723064	LH-723074	5-30	0.50	30	0.6	0.18	0.2	0.06
LH-723065	LH-723075	10-60	1.00	60	0.6	0.36	0.2	0.12

<sup>\*</sup> 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.





#### Germany

Sartorius Lab Instruments GmbH & Co. KG Otto-Brenner-Strasse 20 37079 Göttingen 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

