

Maxicaps® Holder

For “New in-line
Maxicaps® Design”



Product Information

The Sartorius Stedim Maxicaps® holder has been specially designed to integrate Maxicaps® into biopharmaceutical processes. Thanks to its sturdy construction, this holder is highly stable, thus enabling trouble-free integration of Maxicaps® into plant equipment. The open design additionally makes it easy to clean thoroughly.

The holder is designed to accommodate all different Maxicaps® heights. A few easy steps are all it takes to exchange the filter elements. Maxicaps® is attached to the holder using two knurled thumbscrews that can be tightened without the use of any tools. When Maxicaps® is mounted on the holder, all connections are readily accessible.



Technical Data

Technical Description Single Round	
Material	1.4404 (316L) 1.4301 (304) Knurled thumbscrews
Weight	3.3 Kg
Height	337 mm without Maxicaps®
Diameter	178 mm
Root circle	390 mm

Technical Description 3-round	
Material	1.4404 (316L) 1.4301 (304) Knurled thumbscrews
Weight	9.6 Kg
Height	337 mm without Maxicaps®
Diameter	338 mm
Root circle	530 mm

Ordering Information

Ordercode	
Holder single round*	5ZGI-0010
Holder 3-round*	5ZGI-0011

* technical details on request.


Accessories and Spareparts	
Knurled thumbscrew	5ZGB-V0005
Legs (3 pcs)	7EFL-9002
TC Gasket Silicone 1.5" (2 pcs)	7EDSCV0003
TC Gasket Silicone 2.0" (2 pcs)	7EDSCV0004
Clamp closure 1.5"	7ZSB-0009
Clamp closure 2.0"	7ZSB-0012
Manifold for 3-round holder (Tube version)	5ZALB-0002
Manifold for 3-round holder (Plate version)	5ZALB-0005

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

 For further contacts, visit
www.sartorius.com