

# 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.





Single Round



3-round



5-round

## Technical Data

Technical Description	Single Round	3-round	5-round
Material	1.4404 (316L) 1.4301 (304) Knurled thumbscrews	1.4404 (316L) 1.4301 (304) Knurled thumbscrews	1.4404 (316L) 1.4301 (304) Knurled thumbscrews
Weight	3.3 Kg	9.6 Kg	22.8 Kg
Height	337 mm without Maxicaps®	337 mm without Maxicaps®	628 mm without Maxicaps®
Diameter	178 mm	338 mm	511 mm
Root circle	390 mm	530 mm	987 mm



## Ordering Information

### Ordercode

Holder single round*	5ZGI--0010
Holder 3-round*	5ZGI--0011
Holder 5-round*	5ZGI--0013

\* technical details on request.

### Accessories and Spareparts

Knurled thumbscrew	5ZGB-V0005
Legs (3 pcs) - for the 1- and 3-round holder	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



5ZALB-0002




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](http://www.sartorius.com)