

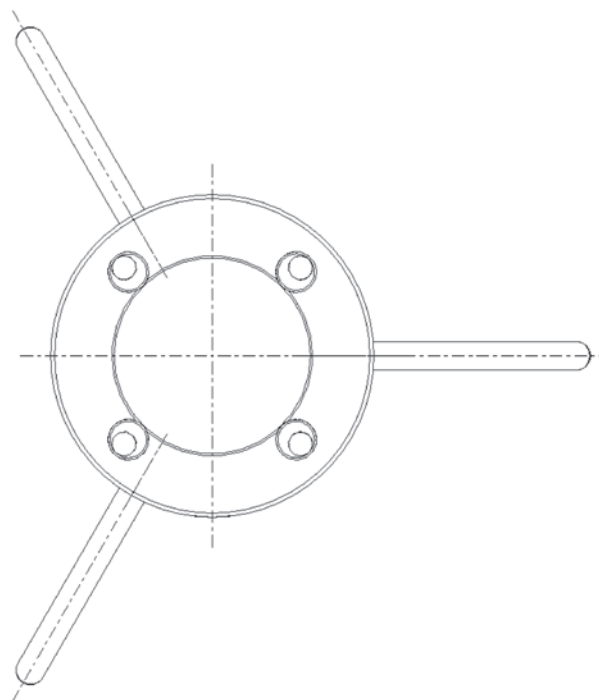
# Maxicaps<sup>®</sup> T-Style Holder

## Description

The Sartorius Stedim T-Style Maxicaps<sup>®</sup> holder has been specially designed to easily integrate Maxicaps<sup>®</sup> into biopharmaceutical processes. This holder is very stable, thanks to its sturdy construction, which enables a trouble-free integration of Maxicaps<sup>®</sup> into existing plant equipment. Its open design makes it easy to clean thoroughly.



The holder is designed to accommodate all 3 T-style Maxicap sizes. Its Stainless steel surfaces are polished to acceptable standards. A few easy steps is all it takes to install & remove the T-style capsules. No special tooling is needed during installation. All connections are readily accessible with the filter mounted in.



## Ordercode

Holder single round	5ZGI--0008
---------------------	------------

## Accessories and Spareparts

Legs (3pcs)	7EFL-9002
TC Gasket Silicone 1.5" (2pcs)	7EDSCV0003
Clamp closure 1.5"	7ZSB-0009

## Technical Description

Material	316 L
Weight	5,1 Kg
Height	620 mm without Maxicaps®
Diameter	165 mm
Root circle	370 mm

### 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)

Specifications subject to change without notice.  
© 2021 Sartorius Stedim Biotech GmbH, August-Spindler-Strasse 11, 37079 Goettingen, Germany

Publication No.: S--2063-e | Order No.: 85037-539-23 | Status: 01 | 29 | 2021