

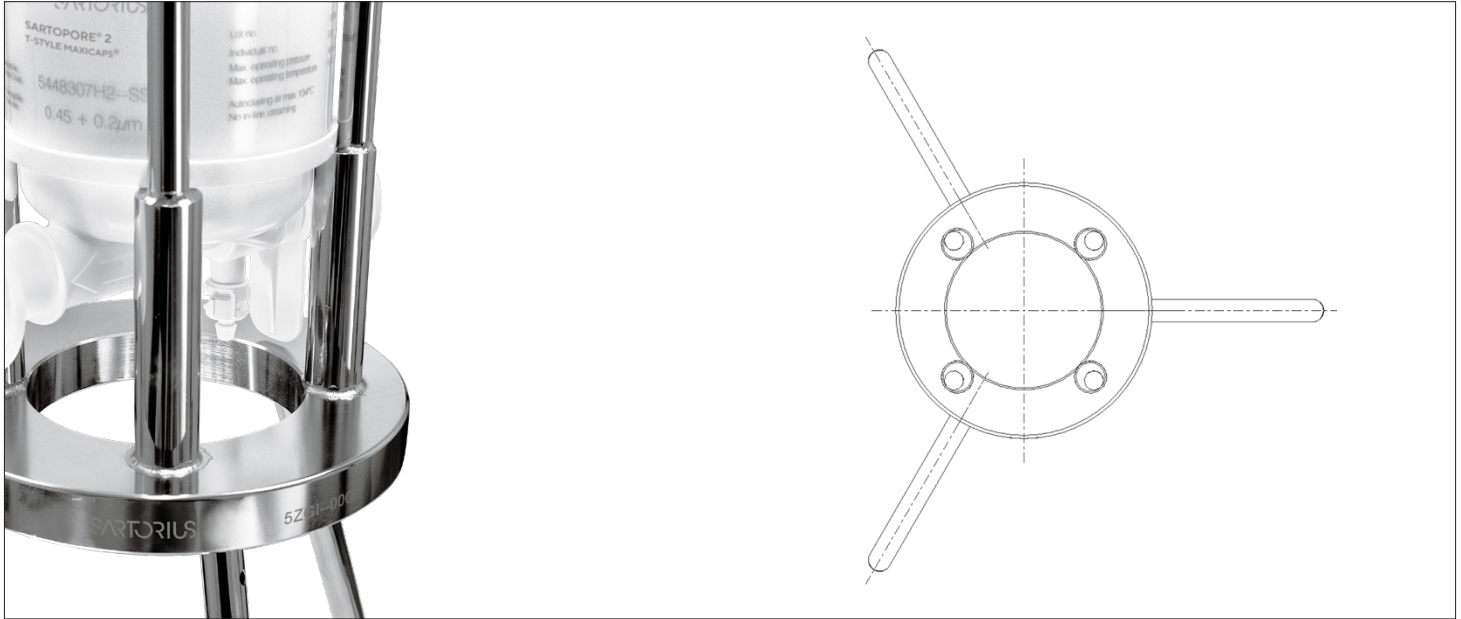
Maxicaps® T-Style Holder



Description

The Sartorius Stedim T-Style Maxicaps® holder has been specially designed to easily integrate Maxicaps® into biopharmaceutical processes. This holder is very stable, thanks to its sturdy construction, which enables a trouble-free integration of Maxicaps® 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.



Ordering Information

Ordercode	
Holder single round	5ZGI--0008
Accessories and Spareparts	
Legs (3 pcs)	7EFL--9002
TC Gasket Silicone 1.5" (2 pcs)	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