

CIP Flush Cassettes

Flush and Clean
Sartorius Stedim Biotech
Filter Holders in Place



Description

Sanitization of experimental and production processes is extremely important in today's biopharmaceutical world. Federal and other institute-based standards require that drug manufacturers provide documentation of the sterility and cleanability of equipment used in the manufacturing process.

With CIP (Clean-In-Place) flush cassettes, Sartorius provides the ideal solution to the sanitization of various filter holders. The CIP system cleans in place to reduce process and validation time while lowering costs.

Applications

CIP (Clean-in-Place) flush cassettes are designed for ideal use in Sartorius filtration systems.

They can be used for the following applications:

- Flushing of filter holders
- Cleaning of filter holders
- Thermal and chemical sterilization processes

Product Information

CIP (Clean-In-Place) flush cassettes provide structural support when performing a cleaning process or a flush on Sartorius filter holders. Flush cassettes can be reused to clean multiple times without compromise of cassette integrity or performance.

Feature	Benefits
Self sealing cassette	No gaskets needed
Silicone sealing compound	No glue
Enlarged inlet and outlet holes	Lower pressure drop

Technical Data

Specifications

Materials of Construction	
Gaskets	PVDF
Spacer	Polypropylene
Sealing compound	Silicone grey

Available Sizes

Sartorius crossflow cassettes are available in standard cassette size for pilot- | production scale and in Sartocon® Slice format for reduced volume handling.

Available Filter Holder

Sartorius crossflow cassettes are designed for Sartorius filter holders like Sartocon® Slice, Sartocon®, Sartocon® 2 Plus, Sartocon® 3, and different Sartoflow® holders.

Operating Parameters	
Feed pressure, P _{in}	58 psi 4 bar maximum
Operating temperature	50 °C maximum, at 20 °C
Cleaning	NaOH, 1 M, 40 °C
Disinfection	NaOH, 1 M, 40 °C, 30 min
Storage	NaOH, 0.1 M

Sterilization

NaOH, 1 M, 40 °C, 30 min
121 °C, 30 min, steaming 121 °C,
110 min autoclaving

Regulatory Compliance

All materials have passed the current USP Biological Test.

Quality Control

Each filter cassette is individually assigned a serial number and integrity tested.

Cassettes comply with cGMP requirements for non-fiber-releasing filters and is filed under the Drug Master File Number DMF 5967 by the Food and Drug administration, Washington, DC. Validation information is available at Sartorius Stedim Biotech GmbH.

If you use holding devices from other suppliers, please contact our Applications Department. A different torque might be needed due to specific variations in design.

For further assistance, please contact your local Sartorius Stedim Biotech field engineer or our Goettingen-based Applications Department in Germany.

Technical References

Validation Guide
Publication No.: SPC5701

Directions for Use (Sartocon® Cassettes and Sartocon® Slice Cassettes) Publication No.: SPC6001

Order Information

Available types and order numbers


Cassette type	Order no.
Sartocon® Cassettes	302 000 00 00E--SG
Sartocon® Slice Cassettes	305 000 00 00E--SG

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