



CIP Flush Cassettes

Flush and clean Sartorius Stedim Biotech filter holders in place



Specifications

Materials of construction

Gaskets	PVDF
Spacer	Polypropylene
Sealing Compound	Silicone grey

Available sizes

Sartorius Stedim Biotech Crossflow Cassettes are available in Standard Cassette size for pilot- | production scale and in Sartocoon Slice format for reduced volume handling.

Available filter holder

Sartorius Stedim Biotech Crossflow Cassettes are designed for Sartorius Stedim Biotech filter holders like SartocoonSlice, Sartocoon, Sartocoon 2 Plus, Sartcon 3, and different Sartoflow holders.

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 Stedim Biotech 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 Stedim Biotech filtration systems. They can be used for the following applications:

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

Product profile

CIP (Clean-In-Place) flush cassettes provide structural support when performing a cleaning process or a flush on Sartorius Stedim Biotech 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

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°, 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

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289
www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178
UK +44.1372.737159
France +33.442.845600
Italy +39.055.63.40.41
Spain +34.91.3586102
Japan +81.3.3740.5407

Specifications subject to change
without notice. Printed and copyrighted
by Sartorius Stedim Biotech GmbH
W/sart-158 · G
Publication No.: SPC2055-e08102
Order No.: 85030-531-65