



Directions for use

Sartoclear® F&B Sheets

1. Introduction

The purpose of this document is to outline the appropriate method for identifying, installing and using a Sartorius Stedim Biotech Sartoclear® F&B Sheets.

Sartoclear® F&B sheets are intended to be used for the clarification fluids in the Foods & Beverages Industries, e.g. wines and beers. The use of this product for other applications or in an other manner as described in this manual may result in personal injury and malfunction of the filter. Please follow the user manual for your own safety and to obtain optimal filtration results.

2. Identification

This document pertains to Sartorius Stedim Biotech Sartoclear® F&B filter sheets with the following article numbers:

40 x 40 cm sheets
290S7-040040FN
290S5-040040FN
290S3-040040FN
290F7-040040FN
290F4-040040FN
290C8-040040FN
290C3-040040FN

60 x 60 cm sheets
290S7-060060FK
290S5-060060FK
290S3-060060FK
290F7-060060FK
290F4-060060FK
290C8-060060FK
290C3-060060FK.

3. Installation

During unpacking and installation it is recommended to wear gloves in order to prevent contamination of the sheets. Please verify that the correct depth filter grade is taken for the application.

Note:

Do not use cracked or broken depth filter sheets.

Insert the sheets with the smooth side facing the filtrate frame in the downstream direction. The rough side of the filter sheet should face the plate in the upstream direction. Please make sure that all edge sealing's cover the edges of the filter sheets. When all the filter sheets are properly sealed the filter unit may be clamped to approximately 50% of its maximal clamping strength. In this manner there is room left for swelling of the sheets due to wetting which secures a proper sealing. The wetting can now be started with the outlet slightly closed so that a backpressure of approximately 20 kPa is created. After 5 minutes of wetting the filter sheets are swollen and the filter unit can be closed completely.

4. Pre-rinsing of the Sartoclear® F&B sheets

Prior to use, Sartoclear® F&B depth filter sheets must be pre-rinsed or flushed with at least 100 l/m² of WFI or buffer in the direction of filtration. The flushing flow rate should be 1,25 times the flow of filtration. In case of a required steam sterilization it is recommended to flush again prior starting the filtration.

5. Sterilization | Sanitization

The filter press and filters may be sterilized by steam sterilization or by using a hot water flush.

5.1 Steam sterilization

The filtration system can be sterilized with steam of up to 134°C (273°F) at a saturated steam pressure of 200 kPa (29 Psi). Before applying steam to the filter pack it is important to loosen the tightened filter pack and slightly open all the vent and drain valves. In this way the steam penetration is optimized and steam shocks will be prevented.

Note:

Please, take care of using particle free steam. The end user is responsible for the effectiveness of the applied steam sterilization procedure in the used equipment.

Note:

Re-tighten the filter press after the system completely cooled down.

5.2. Hot water sanitization

For hot water sanitization it is recommended to use demineralised water which is free of particles. The max allowed temperature is 85°C. The recommended flush rate is 1,5 times the applied filtration speed. The hot water flux should last at least 30 minutes.

Note:

Re-tighten the filter press after the system completely cooled down

6. Filtration

The maximum allowed differential pressure is 1,5 Bar (21,75 psi)

Note:

Constant flow is the preferred filtration mode, where the flow rate through the sheets is held constant during the entire filtration process.

7. Disposal

All components of the Sartoclear® F&B sheets are non-hazardous and can be disposed through normal waste removal methods, in accordance with the local legislation. The disposal method may depend on the nature of the residual material originating from the filtered material.

8. Return of used Filter Elements

All used filter elements should be properly sterilized prior to shipping to Sartorius Stedim Biotech. This allows our staff to handle them with minimal risk during the inspection of the filter elements. The German law requires that a return shipment form (available through your local Sartorius Stedim Biotech representative) must be completed prior to shipping of used filter materials.

9. Liability

Sartorius Stedim Biotech can not assume liability if Sartoclear® F&B Sheets subjected to improper use. In the interest of product development we reserve the right to make changes. Thank you for working with Sartorius Stedim Biotech, we appreciate your business.

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0
www.sartorius.com

© 2021
Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany
MB | Publication No.: SR-6001-e110802
Status: 08 | 2011

List of Sartorius material numbers applying to EPA-FIFRA

290AC4060060-K	290F4H060060-K	290S7-040040BN
290C3-040040BN	290F4H100200-GSC	290S7-040040FN
290C3-040040FN	290F5-040040BN	290S7-060060-K
290C3-060060FK	290F5-060060BK	290S7-060060BK
290C5-040040FN	290F7-040040BNDH	290S7-060060FK
290C5-060060BK	290F7-040040FN	J1-M503LP
290C5-060060FK	290F7-060060BK	U1290F4H040040-N
290C5H060060-K	290F7-060060FK	U1290F7H040040-N
290C5H060060DK	290F7-100200-G	U1290S3-040040-N
290C7-060060-K	290F7H040040-NDH	
290C8-040040-NDH	290K-060120-G	
290C8-040040FN	290K-060120-GSC	
290C8-060060-K	290K-080160-G	
290C8-060060FK	290K-100200-G	
290C8H040040-N	290K-100200-GSC	
290C8H040040DN	290S3-040040-NDH	
290C8H060060-K	290S3-040040FN	
290C8H060060DK	290S3-060060BK	
290C8H100200-GSC	290S3-060060DK	
290C9-060060BK	290S3-060060FK	
290F4-040040-NDH	290S5-040040FN	
290F4-040040FN	290S5-060060BK	
290F4-060060FK	290S5-060060FK	
290F4H040040-NDH	290S7-040040-NDH	