

Jumbo Star The New Filter Generation



Modular Filter Cartridge Construction

Structual integrity

In addition to ease of handling, the new Jumbo construction provides better mechanical stability of the filter system and can withstand back-flushing. The adapter joints between the modules are constructed to meet the most stringent requirements of Food and Beverage applications.

Jumbo Star filter and FDA requirements

All materials used for the construction of Jumbo Star filter meet FDA requirements for use in food and beverage applications. You can choose a wide selection of filters for the different application.

Applications

Filtration of wine and sparkling wine

Jumbo Star filter cartridges are ideally suited for the filtration of wine and sparkling wine.

Typical applications for Jumbo Star in wineries:

- Prefiltration step (last cellar filtration) with Sartoclean SC 30
- Prefiltration with Sartoclean SC 12 upstream of membrane filtration
- Clarification with Sartopure PP2

Benefits:

- Easier to handle and less expensive than pad filters
- Reduces the cost of your operations
- Ensures exact retention ratings
- Requires less rinse water
- Compact and closed filter system
- Easy to sterilize

Filtration of water

Sartorius Stedim Biotech offers a variety of Jumbo Star filter cartridges made to deliver the specific results desired for the filtration of water. Two types of Jumbo filter cartridges are available for particle removal specifically in the range from 0.65 µm to 20 µm and for clarification and reduction of suspended particles.

- Sartopure PP2
- Sartopure GF Plus

Both types do the "heavy work" of removing particles, ensuring an optimal service life of the downstream membrane filter system.

Filtration of beer

Sartorius Stedim Biotech provides a Jumbo Star filter cartridge optimally designed for beer filtration.

- Sartopure PP2

Sartopure PP2 is the ideal choice for particle removal downstream of diatomaceous earth | PVPP filters. This cartridge is available in retention rating of 8 and 20 μ m for reliable particle removal. Even under the most severe conditions, they trap any particles that might have broken through upstream filtration stages, thus ensuring a particle-free filtrate.





The materials of construction of the Jumbo Star filter offer excellent chemical compatibility and thus allow CIP cycles.

Filtration of spirits

Sartorius Stedim Biotech supplies two Jumbo Star filter types for the filtration of spirits:

- Sartopure PP2
 Specifically for particle removing filtration
- Sartopure GF Plus
 For clarification and chill haze removal

Jumbo Star filter cartridges are widely used to clarify cloudy spirits. The closed filter system avoids the oxidation of your product.

Air Venting for reservoir and storage tank

The Jumbo Star Sartopure GA is designed for all venting applications, that do not necessarily require an integrity testable membrane filter cartridge.

Jumbo Star Sartopure GA offers an optimised ratio of air flow rate and retention as well as a great hydrophobic.

Featuring a retention of 0.2 µm for gas Jumbo Star Sartopure GA efficiently protects stored products in the food & beverage industry. Special reservoir and storage tank with a great tankage.







Sartorius Stedim Biotech – Your Knowledgeable Partner for Filtration

Quality from start to finish!All filter cartridges are required to pass stringent in-process quality control tests.

Completely traceable documentation of the entire manufacturing process of Jumbo cartridges is guaranteed by our stringent quality assurance policy that is certified for compliance with ISO 9001. This results in consistently high product quality rangingfrom raw materials to the finished product. Establishing a good partnership with our customers is our No. 1 priority. That's why Sartorius Stedim Biotech offers not only excellent products, but also an international customer support service network that will make your daily work easy.

Our experienced engineers are ready to assist you in solving your filtration problems.





Ordering information

Cartridge type	Adapter type	Retention rating (μm)	1-high (10")	2-high (20")	3-high (30")	4-high (40")
Sartoclean SC	40	1.20	5634003J1	5634003J2	5634003J3	5634003J4
	40	3.00	5634002J1	5634002J2	5634002J3	5634002J4
Sartopure PP2	40	0.65	5594005J1	5594005J2	5594005J3	5594005J4
	40	1.20	5594003J1	5594003J2	5594003J3	5594003J4
	40	3.00	5594002J1	5594002J2	5594002J3	5594002J4
	40	8.00	5594001J1	5594001J2	5594001J3	5594001J4
	40	20.00	5594020J1	5594020J2	5594020J3	5594020J4
Sartopure GF Plus	40	0.65	5554005J1	5554005J2	5554005J3	5554005J4
	40	1.20	5554003J1	5554003J2	5554003J3	5554003J4
Sartopure GA	40	0.20	5594007J1GA		5594007J2GA	



Jumbo Star Filter Cartridge ...the New Filter Generation

Features and benefits

- Highest flow rates
- Highest dirt holding capacityimproved service life
- Modular filter design for easy adaptation and handling
- Improved filter construction for highest thermal and mechanical stability
- Able to handle a broad range of flow requirements

Improved overall design and construction

The large filter area and proprietary pleat design of Jumbo Star provides highest flow rates and total throughputs. Capable of flow rates ranging from 10.000 to 40.000 liter per hour, Jumbo Star filters are especially designed to match requirements for any large volume food and beverage filtration process. Jumbo Star filters feature up to 46 m² of filter area and thus can replace up to twenty four standard filter cartridges. The new modular filters are available in 10", 20", 30" and 40" height allow for easy adaptation into the system.

The rugged construction of Jumbo Star features unique combinations of polypropylene and glass

fiber filter layers. This prevents breakthrough of particles, colloids and bio burden – providing maximum protection for downstream standard filters.

The unique, modular construction of the Jumbo Star filter requires much less set-up work and time than standard filter cartridges technology.

The closed filter system of Jumbo Star allows ease of use and hygienic processing. Product losses and oxidation are also minimized as a result of this design.





Sales and Service Contacts

For further contacts, visit www.sartorius-stedim.com

Europe

Germany

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

Phone +49.551.308.0 Fax +49.551.308.3289

www.sartorius-stedim.com

Sartorius Stedim Systems GmbH Schwarzenberger Weg 73–79 34212 Melsungen

Phone +49.5661.71.3400 Fax +49.5661.71.3702

www.sartorius-stedim.com

France

Sartorius Stedim Biotech S.A. Zl Les Paluds Avenue de Jouques – BP 1051 13781 Aubagne Cedex

Phone +33.442.845600 Fax +33.442.845619

Sartorius Stedim France SAS ZI Les Paluds Avenue de Jouques – CS 71058 13781 Aubagne Cedex

Phone +33.442.845600 Fax +33.442.846545

Austria

Sartorius Stedim Austria GmbH Franzosengraben 12 A-1030 Vienna

Phone +43.1.7965763.18 Fax +43.1.796576344

Belgium

Sartorius Stedim Belgium N.V. Leuvensesteenweg, 248/B 1800 Vilvoorde

Phone +32.2.756.06.80 Fax +32.2.756.06.81

Denmark

Sartorius Stedim Nordic A/S Hoerskaetten 6D, 1. DK-2630 Taastrup

Phone +45.7023.4400 Fax +45.4630.4030

Hungary

Sartorius Stedim Hungária Kft Kagyló u. 5 2092 Budakeszi

Phone +36.23.457.227 Fax +36.23.457.147

Italy

Sartorius Stedim Italy S.p.A. Via dell'Antella, 76/A 50012 Antella-Bagno a Ripoli (FI)

Phone +39.055.63.40.41 Fax +39.055.63.40.526

Netherlands

Sartorius Stedim Netherlands B.V. Edisonbaan 24 3439 MN Nieuwegein

Phone +31.30.6025080 Fax +31.30.6025099

Poland

Sartorius Stedim Poland Sp. z o.o. ul. Wrzesinska 70 62-025 Kostrzyn

Phone +48.61.647.38.40 Fax +48.61.879.25.04

Spain

Sartorius Stedim Spain SA C/Isabel Colbrand 10, Oficina 70 Poligono Industrial de Fuencarral 28050 Madrid

Phone +34.90.2110935 Fax +34.91.3589623

Switzerland

Sartorius Stedim Switzerland AG Ringstr. 24 a 8317 Tagelswangen

Phone +41.52.354.36.36 Fax +41.44.52.354.36.46

U.K.

Sartorius Stedim UK Limited Longmead Business Park Blenheim Road, Epsom Surrey KT19 9 QQ

Phone +44.1372.737159 Fax +44.1372.726171

America

USA

Sartorius Stedim North America Inc. 5 Orville Drive Bohemia, NY 11716

Toll-Free +1.800.368.7178 Fax +1.631.254.4253

Sartorius Stedim SUS Inc. 1910 Mark Court Concord, CA 94520

Phone +1.925.689.6650 Toll Free +1.800.914.6644 Fax +1.925.689.6988

Sartorius Stedim Systems Inc. 201 South Ingram Mill Road Springfield, MO 65802

Phone +1.417.873.9636 Fax +1.417.873.9275

Argentina

Sartorius Argentina S.A. Int. A. Avalos 4251 B1605ECS Munro Buenos Aires

Phone +54.11.4721.0505 Fax +54.11.4762.2333

Brazil

Sartorius do Brasil Ltda Av. Dom Pedro I, 241 Bairro Vila Pires Santo André São Paulo Cep 09110-001

Phone +55.11.4451.6226 Fax +55.11.4451.4369

Mexico

Sartorius de México S.A. de C.V. Circuito Circunvalación Poniente No. 149 Ciudad Satélite 53100 Naucalpan, Estado de México

Phone +52.5555.62.1102 Fax +52.5555.62.2942

Asia | Pacific

Australia

Sartorius Stedim Australia Pty. Ltd. Unit 5, 7-11 Rodeo Drive Dandenong South Vic 3175

Phone +61.3.8762.1800 Fax +61.3.8762.1828

China

Sartorius Stedim Beijing Representative Office No. 33, Yu'an Road, Airport Industrial Zone B, Shunyi District Beijing 101300

Phone +86.10.80426516 Fax +86.10.80426580

Sartorius Stedim Shanghai Represantative Office Room 618, Tower 1, German Centre, Shanghai, PRC., 201203

Phone +86.21.28986393 Fax +86.21.28986392.11

Sartorius Stedim Guangzhou Office Room 704, Broadway Plaza, No. 233–234 Dong Feng West Road Guangzhou 510180

Phone +86.20.8351.7921 Fax +86.20.8351.7931

India

Sartorius Stedim India Pvt. Ltd. #69/2-69/3, Jakkasandra Kunigal Road, Nelamangala Tq Bangalore – 562 123

Phone +91.80.4350.5361 Fax +91.80.4350.5253

Japan

Sartorius Stedim Japan K.K. KY Building, 8–11 Kita Shinagawa 1-chome Shinagawa-ku Tokyo 140-0001

Phone +81.3.3740.5407 Fax +81.3.3740.5406

Malaysia

Sartorius Stedim Malaysia Sdn. Bhd. Lot L3-E-3B, Enterprise 4 Technology Park Malaysia Bukit Jalil 57000 Kuala Lumpur

Phone +60.3.8996.0622 Fax +60.3.8996.0755

Singapore

Sartorius Stedim Singapore Pte. Ltd. 1 Science Park Road, The Capricorn, #05-08A, Singapore Science Park 2 Singapore 117528

Phone +65.6872.3966 Fax +65.6778.2494