

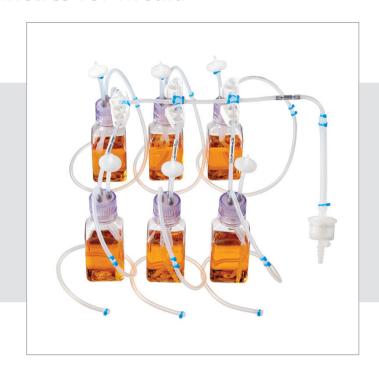
MYCAP® Manifold Bottle Assemblies for Media

Product Information

MYCAP® bottle closures are a one-piece, ready-to-use aseptic fluid transfer system.

The single-use design features integral tubing and a robust platinum-cured silicone seal.

MYCAP® bottle assemblies include Nalgene® square media bottles, Sartopore® 2 filter, are gamma irradiated and ready for immediate use.



Features	Benefits
– Ready-to-use	Save timeEliminate errors
Single-use	 No contamination risk
 One-For-All silicone closure 	 Universal validation Robust & reliable performance
- One-piece construction	 Integral closure Easy to handle
 Easy customization 	 Configure MYCAP® to your process needs
– Sartorpore [®] 2 Filter	 Sterile filter Media into media bottle

Applications

MYCAP® for Media make media distribution and storage quicker and easier than ever before. MYCAP® for Media includes six bottles.

MYCAP® for Media includes a Sartopore® 2 filter for sterile filtration directly into easy-to-handle bottles. Simply connect the filter inlet to your bulk media source to sterilize media as the bottle fills.

Cut the QUICKSEAL® aseptic disconnect after filling and your aseptic media distribution is complete, without ever having to go into a biosafety cabinet.

Choose tube welding or AseptiQuik® aseptic connector for convenient aseptic connection and dispense sterile media at point-of-use, including bioreactors.

Reduce Risk & Improve Performance With MYCAP® bottle closures, every use is its first use.

There is no risk of batch-to-batch or product-to-product cross contamination.

MYCAP® bottle assemblies pass studies of container closure by bacterial immersion.

One-For-All Closure System

At last – a truly universal bottle closure system.

All MYCAP® closure systems are constructed using the same high performance platinum-cured silicone seal regardless of the tubing material, tubing size, port configuration, bottle type or bottle size.

Ease validation and reduce regulatory work by implementing MYCAP® throughout the production process.

Easy Implementation

Alternative bottle closure systems include a number of reusable components which are costly to implement and manage.

Regulatory agencies demand a scientifically sound validation strategy for parts cleaning to avoid non-host or batch-to-batch contamination. The physical condition and life cycle of reused parts must be closely monitored to ensure cleaning efficacy.

Select pre-assembled and single-use MYCAP® bottle assemblies to bypass the burden of validation and parts management.

Reliable Performance

The platinum-cured silicone seal of MYCAP® is suitable for a range of applications.

The closure to the bottle is secure so long as the seal maintains its physical properties. Silicone maintains its elastomeric properties from -70° to 260° Celcius.

Validation demonstrates a leak-free seal when MYCAP® are installed with torqueforce from 4in-lbs to 60in-lbs.

Technical Data

Specifications

Elastomer Seal	Platinum-cured silicone
Tubing	Platinum-cured silicone or C-Flex® formulation 374
Cap Material	Polycarbonate (not fluid contact)
Bottle Style	Nalgene® PETG bottle
Tube Ports	Inlet: Sartopore® 2 Media Filter Air vent - 25 mm PES membrane MiniSart®
Fittings	Outlet: Aseptiquik® S or tube plug
Sterilization	Gamma Irradiated

Manufacturing Environment

 MYCAP° bottle closures are constructed and packaged in an ISO 7 clean room.

Monthly cleaning with a schedule of LpH®, Vesphene® and Spor Klenz® reduce and control viable organisms. Viable organisms are measured quarterly to monitor the effectiveness of the clean room management and cleaning procedures and to be compliant to EU GMPs and ISO14698.

Line clearance, weekly cleaning of equipment and work surfaces and monthly cleaning of the cleanroom reduce and control non-viable particles. Non-viable readings are taken weekly to ensure $0.5 \ \mu m/m^3$ and $5.0 \ \mu m/m^3$ particles are within the ISO Class 7 acceptance criteria, per ISO 14644-1.

Quality Assurance

MYCAP® bottle assemblies are manufactured in accordance with the key principles of cGMP.

Each MYCAP® bottle closure is issued with a unique lot number on its' Certificate of Compliance. The lot number provides full traceability to all components used and operations performed.

The Certificate of Conformance for MYCAP® bottle assemblies reports compliance to:

- USP <87>: Biological reactivity tests, in Vitro
- USP <88>: Biological reactivity tests, in Vivo
- USP <788>: Particulate matter
- USP <85>: Endotoxin
- Sterility Assurance (SAL) 10⁻⁶ per ISO11137
- USP <381>: Extractables
- 21CFR177.2600: Extractables

Shelf Life

MYCAP® bottle closures, components and component resins have shelf-life that exceeds 2 years.

TSE BSE Risk

The components and materials used in construction of MYCAP® bottles assemblies are treated at high temperatures and undergo chemical reactions which far exceed the stringent requirements laid down in the documents CPMP|BWP|1230|98 and WHO | CPS | VPH | 95.145 and | or EMA | 410 | 01, rev 3. These processing conditions are considered sufficient to inactivate TSE & BSE transmitters.

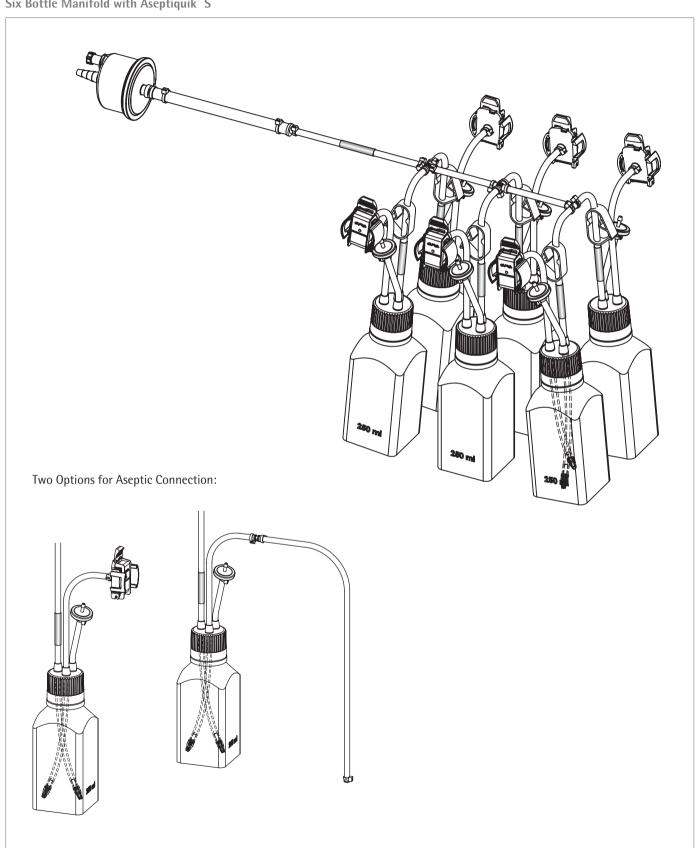
C-Flex® is a registered trademark of Saint-Gobain Performance Plastics Corporation Luer® is a registered trademark of Becton Dickinson LpH®, Vesphene® Spor Klenz® are registered trademarks of STERIS Corporation National Registeries and Property of Theorems Fisher Scientific

Nalgene® is a registered trademark of ThermoFisher Scientific Sartopore®, Minisart®, QUICKSEAL® and MYCAP® is a registered trademark of Sartorius-Stedim Biotech

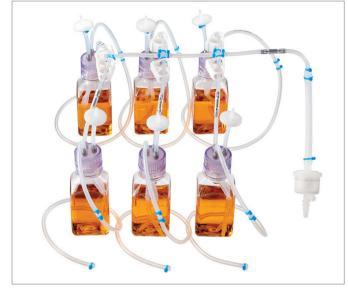
AseptiQuik®, is a registered trademark of Colder Products Company

Ordering Information

Six Bottle Manifold with Aseptiquik® S







Before Media Fill After Media Fill

Six Bottle Manifold with Aseptiquik® S on C-Flex® for Aseptic Connections

Universal Design Features

- Optimized Tube Lengths including dip tube for fluid transfer
- Sartopore[®] 2 0.2 μm Filter at Manifold Inlet for Sterile Media Fill
- Includes Nalgene® PETG Media Bottle
- Gamma Irradiated
- Individually Double Bagged

Article Number	Description	Bottle Size	Port 1	Port 2	Port 3	Qty/Pack
MCMD1000060204B	6×1000 mL Square Media Bottle; 1/8×1/4 C-Flex® Tubing w/Aseptiquik® S	1000 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; 1/4 – 1/2 barb inlet; QUICKSEAL®; dip tube to bottom	Aseptiquik®; dip tube to bottom	Minisart® hydropho- bic PES vent filter	1
MCMD0500060204B	6×500 mL Square Media Bottle; 1/8×1/4 C-Flex® Tubing w/Aseptiquik® S	500 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; 1/4 – 1/2 barb inlet; QUICKSEAL®; dip tube to bottom	Aseptiquik®; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0250060204B	6×250 mL Square Media Bottle; ½8×½ C-Flex [®] Tubing w/Aseptiquik [®] S	250 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; ¹ / ₄ – ¹ / ₂ barb inlet; QUICKSEAL®; dip tube to bottom	Aseptiquik®; dip tube to bottom	Minisart® hydropho- bic PES vent filter	1

Six Bottle Manifold with Aseptiquik® S for Aseptic Connections

Universal Design Features

- Optimized Tube Lengths including dip tube for fluid transfer
- Sartopore® 2 0.2 μm Filter at Manifold Inlet for Sterile Media Fill
- Includes Nalgene® PETG Media Bottle
- Gamma Irradiated
- Individually Double Bagged

Article Number	Description	Bottle Size	Port 1	Port 2	Port 3	Qty/Pack
MCMD1000060204A	6×1000 mL Square Media Bottle; ½8×¼ Silicone Tubing; Aseptiquik® S	1000 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; 1/4 – 1/2 barb inlet; QUICKSEAL®; dip tube to bottom	Aseptiquik®; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0500060204A	6×500 mL Square Media Bottle; 1/8×1/4 Silicone Tubing; Aseptiquik® S	500 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; 1/4 – 1/2 barb inlet; QUICKSEAL®; dip tube to bottom	Aseptiquik [®] ; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0250060204A	6×250 mL Square Media Bottle; ½8×½ Silicone Tubing; Aseptiquik® S	250 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; 1/4 – 1/2 barb inlet; QUICKSEAL®; dip tube to bottom	Aseptiquik®; dip tube to bottom	Minisart® hydropho- bic PES vent filter	1

Six Bottle Manifold with C-Flex® for Aseptic Connections via Tube Welding

Universal Design Features

- Optimized Tube Lengths including dip tube for fluid transfer
- Sartopore® 2 0.2 μm Filter at Manifold Inlet for Sterile Media Fill
- Includes Nalgene® PETG Media Bottle
- Gamma Irradiated
- Individually Double Bagged

Article Number	Description	Bottle Size	Port 1	Port 2	Port 3	Qty/Pack
MCMD1000060204W	6 × 1000 mL Square Media Bottle; 1/8 × 1/4 C-Flex® Tubing	1000 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; 1/4 – 1/2 barb inlet; QUICKSEAL®; dip tube to bottom	Tube Plug; dip tube to bottom	Minisart® hydropho- bic PES vent filter	1
MCMD0500060204W	6×500 mL Square Media Bottle; 1/8×1/4 C-Flex® Tubing	500 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; 1/4 – 1/2 barb inlet; QUICKSEAL®; dip tube to bottom	Tube Plug; dip tube to bottom	Minisart® hydrophobic PES vent filter	1
MCMD0250060204W	6×250 mL Square Media Bottle; ½8×½ C-Flex® Tubing	250 mL	(Manifold Inlet) Size 5 Sartopore® 2 capsule filter; 0.2 µm; 1/4 – 1/2 barb inlet; QUICKSEAL®; dip tube to bottom	Tube Plug; dip tube to bottom	Minisart® hydropho- bic PES vent filter	1

Europe

Germany Sartorius Lab Instruments GmbH & Co. KG Otto-Brenner-Strasse 20 37079 Goettingen

Phone +49.551.308.0

France & Suisse Romande

Sartorius France 2. rue Antoine Laurent de Lavoisier ZA de la Gaudrée 91410 Dourdan

Phone +33.1.70.62.50.00

Austria

Sartorius Austria GmbH Modecenterstrasse 22 1030 Vienna

Phone +43 1 7965760 0

Belgium

Sartorius Belgium N.V. Rue Colonel Bourg 105 1030 Bruxelles

Phone +32.2.756.06.90

Finland & Baltics

Sartorius Biohit Liquid Handling Oy Lainnatie 1 00880 Helsinki

Phone +358.9.755.951

Hungary

Sartorius Hungária Kft. Kagyló u. 5. 2092 Budakeszi

Phone +3623.457.227

Ireland

Sartorius Ireland Ltd. Unit 41, The Business Centre Stadium Business Park Ballycoolin Road Dublin 11

Phone +353.1.8089050

Italy

Sartorius Italy S.r.l. Via Torino 3/5 20814 Varedo (MB)

Phone +39.0362.5557.11

Netherlands

Sartorius Netherlands B.V.

Phone +31.30.60.53.001 Fax +31.30.60.52.917

info.netherlands@sartorius.com

Poland

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

Phone +48.61.6473830

Russian Federation

LLC "Sartorius RUS' Vasilyevsky Island 5th line 70, Lit. A 199178 St. Petersburg

Phone +7.812.327.53.27

Spain & Portugal

Sartorius Spain, S.A. Avda. de la Industria, 32 Edificio PAYMA 28108 Alcohendas (Madrid)

Phone Spain +34.913.586.095 Phone Portugal +351.800.855.800

Switzerland

Sartorius Mechatronics Switzerland AG Ringstrasse 24a 8317 Tagelswangen (ZH)

Phone +41.44.746.50.00

Sartorius UK Ltd. Longmead Business Centre Blenheim Road, Epsom Surrey KT19 9QQ

Phone +44.1372.737159

Ukraine

LLS "Sartorius RUS" Post Box 440 "B" 01001 Kiev, Ukraine

Phone +380.44.411.4918

Americas

IISA

Sartorius Corporation 5 Orville Drive, Suite 200 Bohemia, NY 11716

Phone +1.631.254.4249 Toll-free +1.800.635.2906

Argentina

Sartorius Argentina S.A. Int. A. Ávalos 4251 B1605ECS Munro **Buenos Aires**

Phone +54.11.4721.0505

Brazil

Sartorius do Brasil Ltda Avenida Senador Vergueiro 2962 São Bernardo do Campo CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900

Canada

Sartorius Canada Inc 1173 North Service Road West, D4 Oakville, ON L6M 2V9

Phone +1.905.569.7977 Toll-Free +1.800.668.4234

Mexico

Sartorius de México, S.A. de C.V. Libramiento Norte de Tepotzotlan s/n, Colonia Barrio Tlacateco, Municipio de Tepotzotlan, Estado de México, C.P. 54605

Phone +52.55.5562.1102 leadsmex@sartorius.com

Peru

Sartorius Peru S.A.C. Avenue Alberto del Campo 411 Floor 12 - The Office 15076 - San Isidro, Lima

Phone +51.1.441 0158

Asia | Pacific

Australia

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

Phone +61.3.8762.1800

China

Sartorius (Shanghai) Trading Co., Ltd. 3rd Floor, North Wing, Tower 1 No. 4560 Jinke Road Zhangjiang Hi-Tech Park **Pudong District** Shanghai 201210, P.R. China

Phone +86.21.6878.2300

Hong Kong

Sartorius Hong Kong Ltd. Unit 1012, Lu Plaza 2 Wing Yip Street Kwun Tong Kowloon, Hong Kong

Phone +852.2774.2678

Sartorius Weighing India Pvt. Ltd. #69/2-69/3, NH 48, Jakkasandra, Nelamangala Tg 562 123 Bangalore, India

Phone +91.80.4350.5250

Japan

Sartorius Japan K.K. 4th Fl., Daiwa Shinagawa North Bldg. 8-11, Kita-Shinagawa 1-chome Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.3740.5408

Malaysia

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

Phone +60.3.8996.0622

Singapore

Sartorius Singapore Pte. Ltd 10 Science Park Rd The Alpha #02-13/14 Singapore Science Park II Singapore 117684

Phone +65.6872.3966

South Korea

Sartorius Korea Ltd. 8th Floor, Solid Space B/D, PanGyoYeok-Ro 220, BunDang-Gu SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700

Thailand

Sartorius (Thailand) Co. Ltd. 129 Rama 9 Road, Huaykwang Bangkok 10310

Phone +66.2643.8361-6