



1: Sartopore® 2 capsule filter; 0.2µm; 300cm²; 1/4"-1/2" HB I/O (GI)

3: 1/4"x1/8"HB Reducer

9: Ratchet clamp; 1/4"OD tubing

10: 1/8"HB Straight Connector

13: Vent Filter; 25mm MiniSart; 1/8"HB I/O; PES(GI/Autoclave)(Hydrophobic)

6: MYCAP™; 38-430 Cap; 3-Ports;
 P1: 1/8"x1/4" Dow Pharma 50 Sil, CS-11", BS-175mm, QS-2";
 P2: 1/8"x1/4" Dow Pharma 50 Sil, CS-11", BS-175mm, QS-0";
 P3: 1/8"x1/4" Dow Pharma 50 Sil, CS-3", BS-0", QS-0"

11: 1/8"IDx1/4"OD C-FLEX® 374 tube; 18"OAL

12: Colder AseptiQuik® S; 1/8"HB; Not Autoclaveable

2: 1/4"IDx7/16"OD Dow Pharma 50 Silicone tube; 6"OAL

4: 1/8"IDx1/4"OD Dow Pharma 50 Silicone tube; 6"OAL w/QS collar at 3"

14: 1/8"IDx1/4"OD Dow Pharma50 Silicone tube; 3"OAL

5: 1/8"HB Cross

15: 1/8"HB Tee

7: 1/8" Dip Tube Tip w/Notch

8: 500ml PETG Bottle

Bill of Materials:

Item	Qty	AP Number	Material
1	1	AP-3311	-PP, PES, Polyester -Silicone
2	1	AP-2119	-ADCF Silicone
3	1	AP-0029	-ADCF Polypropylene
4	1	AP-0827	-ADCF Silicone
5	2	AP-0638	-ADCF Polypropylene
6	6	AP-3312	-ADCF Silicone
7	12	AP-2170	-ADCF Polypropylene
8	6	AP-0298	-PETG
9	6	AP-0118	-Polypropylene
10	6	AP-0089	-ADCF Polypropylene
11	6	AP-0378	-ADCF C-FLEX®(374)
12	6	AP-0917	-PC, PE, Silicone
13	6	AP-2767	-Polypropylene, PES
14	2	AP-2291	-ADCF Silicone
15	1	AP-1488	-ADCF Polypropylene
16	38	AP-0133	-Tefzel®

Revision History:

- 0 - Initial release
- 1 - Adjusted Sartopore® filter (AP-3311) for improved image scale

Dimensional Tolerances:

- Tubing Lengths:
- 0-12" ±0.25"
- 0-36" ±0.5"
- 36"+ ±1.0"

Notes:

1. No assembly required.
2. Certificate of Conformance & Certificate of Irradiation provided for each lot.
3. Individually double bagged.
4. Fluid contact materials meet criteria for USP Class VI Testing. Consult Validation Guide for additional tests and details.
5. Gamma Irradiated.
6. Drawing not to scale.

PROPRIETARY AND CONFIDENTIAL

The information contained herein is the sole property of AllPure Technologies, LLC. Reproduction or distribution is prohibited without written consent from AllPure Technologies, LLC.

Title:
6x500ml Media; 1/8"x1/4" C-FLEX® w/ASQ

sartorius stedim
 biotech
 AllPure Technologies, LLC
 80 Progress Ave
 New Oxford, PA 17350
 www.sartorius.com
 www.allpureinc.com

Customer Approval: Approval Not Required
 Approval Date:
 Part Number: MCMD0500060204B
 Drawing Number: MCMD0500060204B

Revision: 1 Released By: Release Date: