



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

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

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

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

5: 1/8"HB Cross

9: Ratchet clamp; 1/4"OD tubing

10: 1/8"HB Straight Connector

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

12: 1/8"HB Tube Plug

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

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

15: 1/8"HB Tee

16: Cable tie, blue; TYP

6: MYCAP™; 38-430 Cap; 3-Ports; P1: 1/8"x1/4" Dow Pharma 50 Sil, CS-9.5", BS-215mm, QS-1.5"; P2: 1/8"x1/4" Dow Pharma 50 Sil, CS-9.5", BS-215mm, QS-0"; P3: 1/8"x1/4" Dow Pharma 50 Sil, CS-3", BS-0", QS-0"

**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-3110	-ADCF Silicone
7	12	AP-2170	-ADCF Polypropylene
8	6	AP-1330	-PETG
9	6	AP-0118	-Polypropylene
10	6	AP-0089	-ADCF Polypropylene
11	6	AP-0378	-ADCF C-FLEX®(374)
12	6	AP-1181	-ADCF Polypropylene
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:**

6x1000ml Media; 1/8"x1/4" C-FLEX®



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: MCMD1000060204W

Drawing Number: MCMD1000060204W

Revision: 1

Released By:

Release Date: