



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

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

8: 1/8"HB Straight Connector

1: Sartopore® 2 capsule filter; 0.2µm; 150cm²; 1/4"-1/2" HB I/O; (GI)

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

11: Cable tie, blue; TYP

4: 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"

6: 500ml PETG Bottle

5: 1/8" Dip Tube Tip w/Notch

9: 1/8"IDx1/4"OD C-FLEX® 374 tube; 18"OAL

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

**Bill of Materials:**

Item	Qty	AP Number	Material
1	1	AP-2365	-PP, PES, Polyester -Silicone
2	1	AP-2119	-ADCF Silicone
3	1	AP-0029	-ADCF Polypropylene
4	1	AP-3312	-ADCF Silicone
5	2	AP-2170	-ADCF Polypropylene
6	1	AP-0298	-PETG
7	1	AP-2767	-Polypropylene, PES
8	1	AP-0089	-ADCF Polypropylene
9	1	AP-0378	-ADCF C-FLEX®(374)
10	1	AP-0917	-PC, PE, Silicone
11	7	AP-0133	-Tefzel®

**Revision History:**

- 0 - Initial release
- 1 - Adjusted Sartopore® filter (AP-2365) 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.

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

Customer Approval: Approval Not Required	Revision: 1	Released By:	Release Date:
Approval Date:			
Part Number: MCMD0500010204B			
Drawing Number: MCMD0500010204B			

Title:  
 500ml Media; 1/8"x1/4" C-FLEX® w/ASQ