



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-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"

6: 1000ml PETG Bottle

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

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

10: 1/8"HB Tube Plug

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-3110	-ADCF Silicone
5	2	AP-2170	-ADCF Polypropylene
6	1	AP-1330	-PETG
7	1	AP-2767	-Polypropylene, PES
8	1	AP-0089	-ADCF Polypropylene
9	1	AP-0378	-ADCF C-FLEX®(374)
10	1	AP-1181	-ADCF Polypropylene
11	7	AP-0133	-Tefzel®

Revision History:

- 0 - Initial release
- 1 - Adjusted Sartopore® filter (AP-2365) for improved image scale

Dimensional Tolerances:

- 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	Title: 1000ml Media; 1/8"x1/4" C-FLEX®	
Approval Date:		
Part Number: MCMD1000010204W		
Drawing Number: MCMD1000010204W		
Revision: 1	Released By:	Release Date: