



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

11: 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"

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"

12: 1/8"IDx1/4"OD Dow Pharma50  
Silicone tube; 3"OAL

14: Cable tie,  
blue; TYP

5: 1/8"HB Cross

13: 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-0917	-PC, PE, Silicone
11	6	AP-2767	-Polypropylene, PES
12	2	AP-2291	-ADCF Silicone
13	1	AP-1488	-ADCF Polypropylene
14	26	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" ASQ

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: MCMD0500060204A  
Drawing Number: MCMD0500060204A

Revision: 1      Released By:      Release Date: