



**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-0917	-PC, PE, Silicone
9	5	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.

**Title:**

1000ml 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: MCMD1000010204A

Drawing Number: MCMD1000010204A

Revision: 1

Released By:

Release Date: