



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-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.

Title:
 1000ml Media; 1/8"x1/4" C-FLEX® w/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: MCMD1000010204B
 Drawing Number: MCMD1000010204B

Revision: 1 Released By: Release Date: