



**Bill of Materials:**

Item	Qty	AP Number	Material
1	1	AP-1181	-ADCF Polypropylene
2	1	AP-3212	-ADCF C-FLEX®(374) -ADCF Silicone
3	1	AP-1048	-HDPE
4	1	AP-2170	-ADCF Polypropylene
5	1	AP-0639	-Polystyrene
6	1	AP-2767	-Polypropylene, PES
7	2	AP-0133	-Tefzel®

**Revision History:**

0 - Initial release  
 1 - P1: CS 13.25" to 13.5"; BS 124mm to 117mm;  
 QS 2.25" to 2.5"

**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.



AllPure Technologies, LLC  
 80 Progress Ave  
 New Oxford, PA 17350

biotech  
 www.sartorius.com  
 www.allpureinc.com

Customer Approval: Approval Not Required

Approval Date:

Part Number: MYCAP CPS G0015D

Drawing Number: MYCAP CPS G0015D

Revision: 1

Released By:

Release Date:

Title:

MYCAP™: 15ml PS Centrifuge; 2-Ports;  
 1/8"x1/4" C-FLEX® w/Dip Tube Tip; QS on inlet