



**Bill of Materials:**

Item	Qty	AP Number	Material
1	1	AP-0099	-ADCF Polypropylene
2	1	AP-0031	-ADCF Polypropylene
3	1	AP-3220	-ADCF Silicone
4	1	AP-2170	-ADCF Polypropylene
5	1	AP-0202	-Polypropylene
6	1	AP-2767	-Polypropylene, PES
7	2	AP-0133	-Tefzel®

**Revision History:**

0 - Initial release

**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 CPP G0050A

Drawing Number: MYCAP CPP G0050A

Revision: 0

Released By:

Release Date:

**Title:**

MYCAP™: 50ml PP Centrifuge; 2-Ports;  
 1/8"x1/4" Silicone w/Dip Tube Tip