



Bill of Materials:

| Item | Qty | AP Number | Material |
|------|-----|-----------|--------------------------------------|
| 1 | 1 | AP-1181 | -ADCF Polypropylene |
| 2 | 1 | AP-3218 | -ADCF C-FLEX®(374) -ADCF Silicone |
| 3 | 1 | AP-2170 | -ADCF Polypropylene |
| 4 | 1 | AP-0202 | -Polypropylene |
| 5 | 1 | AP-2767 | -Polypropylene, PES |
| 6 | 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 G0050B

Drawing Number: MYCAP CPP G0050B

Revision: 0

Released By:

Release Date:

Title:

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