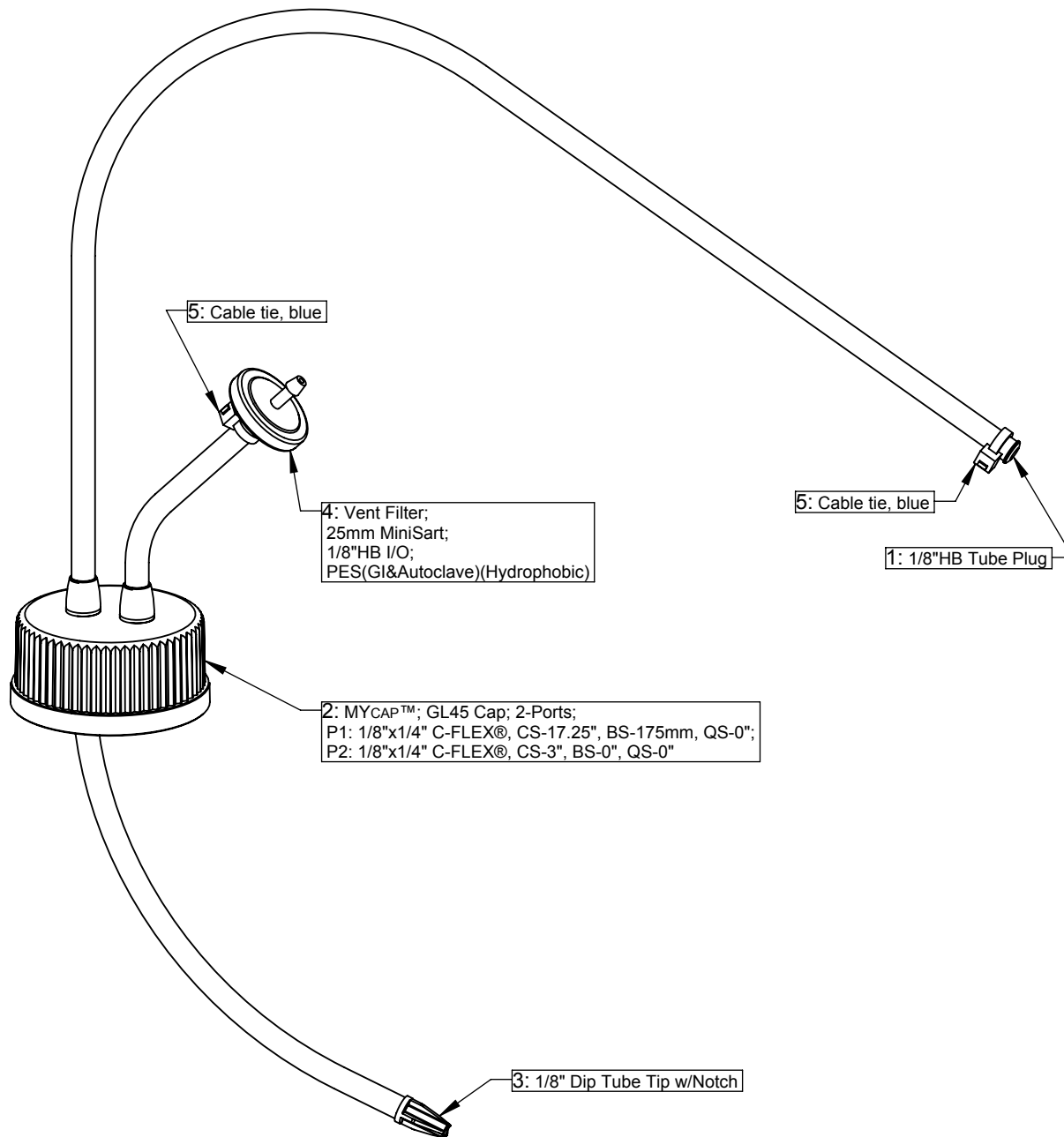


**500ml Corning Pyrex Glass, 1395-500**



**Bill of Materials:**

Item	Qty	AP Number	Material
1	1	AP-1181	-ADCF Polypropylene
2	1	AP-2808	-ADCF C-FLEX®(374) -ADCF Silicone
3	1	AP-2170	-ADCF Polypropylene
4	1	AP-2767	-Polypropylene, PES
5	2	AP-133	-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. Individually bagged then double bagged, 5 per bag.
3. Fluid contact materials meet criteria for USP Class VI Testing. Consult Validation Guide for additional tests and details.
4. Drawing not to scale.

**PROPRIETARY AND CONFIDENTIAL**

The information contained herein is the sole property of AllPure Technologies, Inc. Reproduction or distribution is prohibited without written consent from AllPure Technologies, Inc.

**Title:**

MYCAP™; GL45 Cap; 2-Ports; 1/8"x1/4"C-FLEX® w/Dip Tube - 500ml Corning Glass Bottle; NOT IRRADIATED



80 Progress Ave  
New Oxford, PA 17350  
(717) 624-3241  
(717) 624-3051 (fax)

Customer Approval: Approval Not Required

Approval Date:

Part Number: MYCAP 13950500B

Drawing Number: MYCAP 13950500B

Revision: 0

Released By:

Release Date: