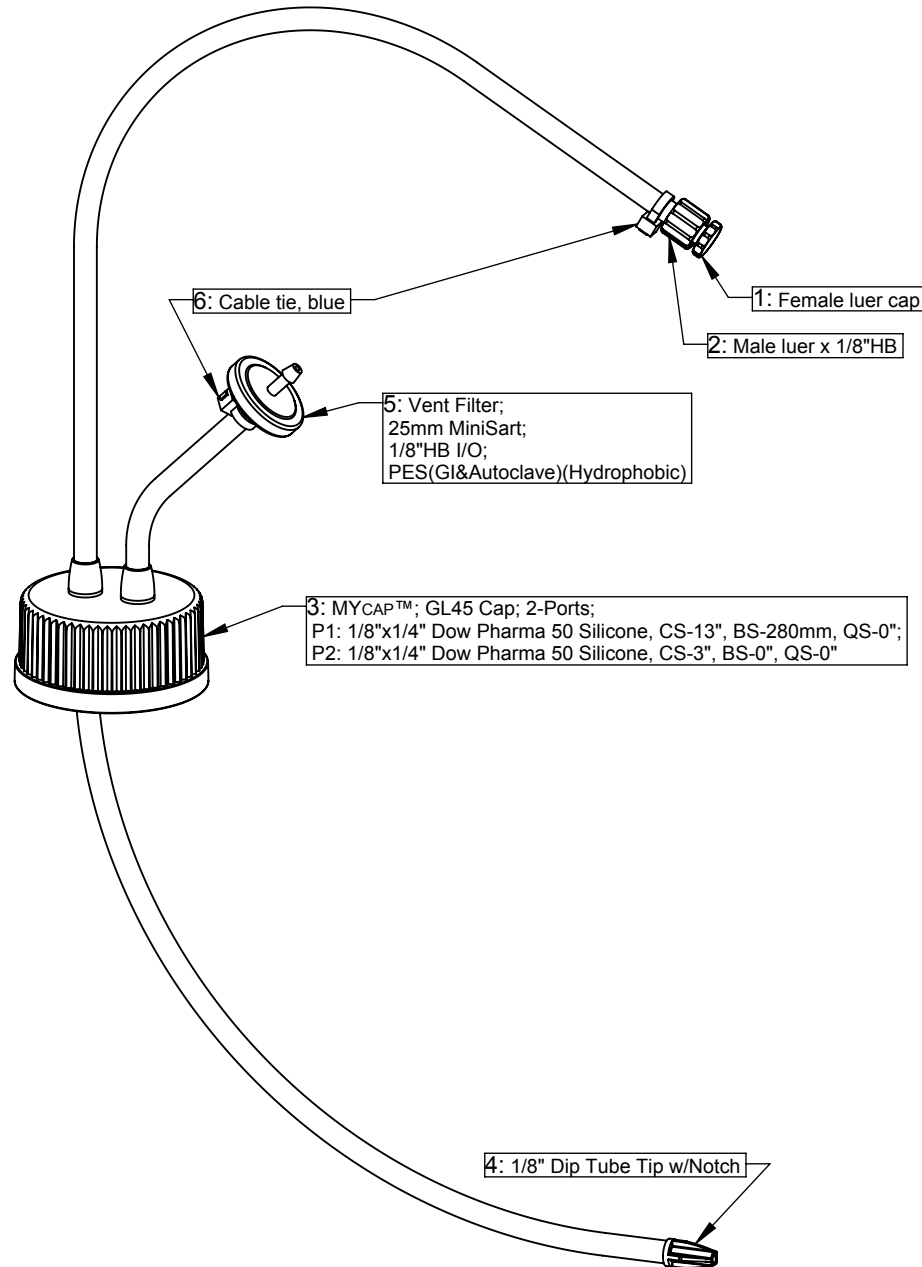


2000ml Corning Pyrex Glass, 1395-2000



Bill of Materials:

Item	Qty	AP Number	Material
1	1	AP-0099	-ADCF Polypropylene
2	1	AP-0031	-ADCF Polypropylene
3	1	AP-2776	-ADCF Silicone
4	1	AP-2170	-ADCF Polypropylene
5	1	AP-2767	-Polypropylene, PES
6	2	AP-0133	-Tefzel®

Revision History:

- 0 - Initial release
- 1 - P1 CS 14" to 13", BS 252mm to 280mm
- Update template to include Sartorius name/logo

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 Sartorius Stedim North America. Reproduction or distribution is prohibited without written consent from Sartorius Stedim North America.



Sartorius Stedim North America, Inc.
 80 Progress Ave
 New Oxford, PA 17350

biotech
 www.sartorius.com

Customer Approval: Approval Not Required

Approval Date:

Part Number: MYCAP 13952000A

Drawing Number: MYCAP 13952000A

Revision: 1

Released By:

Release Date:

Title:

MYCAP™; GL45 Cap; 2-Ports; 1/8"x1/4" Silicone w/Dip Tube - 2000ml Corning Glass Bottle;
 NOT IRRADIATED