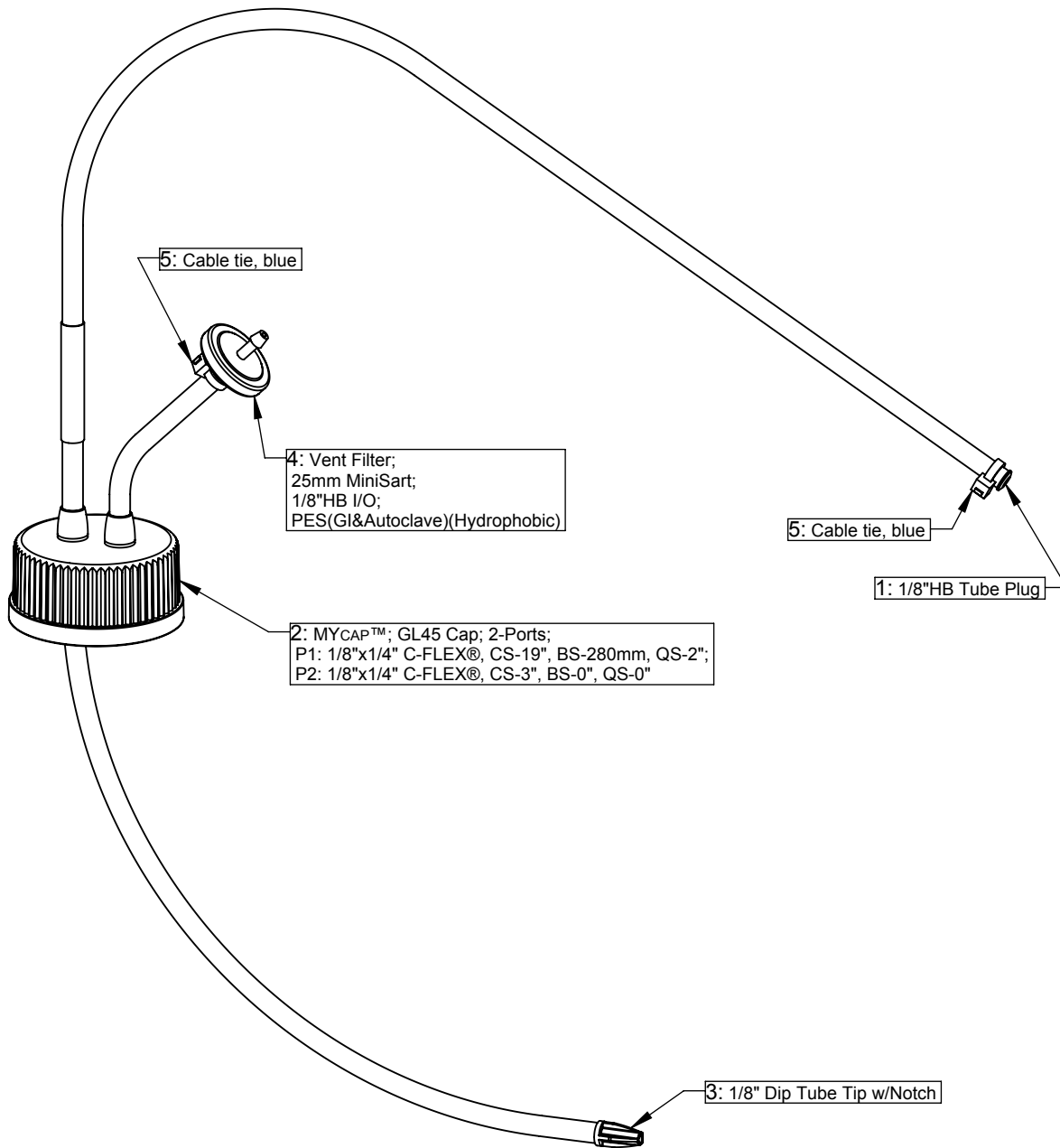


2000ml Corning Pyrex Glass, 1395-2000



5: Cable tie, blue

4: Vent Filter;
25mm MiniSart;
1/8"HB I/O;
PES(GI&Autoclave)(Hydrophobic)

2: MYCAP™; GL45 Cap; 2-Ports;
P1: 1/8"x1/4" C-FLEX®, CS-19", BS-280mm, QS-2";
P2: 1/8"x1/4" C-FLEX®, CS-3", BS-0", QS-0"

5: Cable tie, blue

1: 1/8"HB Tube Plug

3: 1/8" Dip Tube Tip w/Notch

Bill of Materials:

Item	Qty	AP Number	Material
1	1	AP-1181	-ADCF Polypropylene
2	1	AP-3062	-ADCF C-FLEX®(374) -ADCF Silicone
3	1	AP-2170	-ADCF Polypropylene
4	1	AP-2767	-Polypropylene, PES
5	2	AP-0133	-Tefzel®

Revision History:

- 0 - Initial release
- 1 - P1 CS 20.25" to 19", 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:

- No assembly required.
- Individually bagged then double bagged, 5 per bag.
- Fluid contact materials meet criteria for USP Class VI Testing. Consult Validation Guide for additional tests and details.
- 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 13952000D

Drawing Number: MYCAP 13952000D

Revision: 1

Released By:

Release Date:

Title: MYCAP™; GL45 Cap; 2-Ports;
1/8"x1/4"C-FLEX® w/Dip Tube;
QS on inlet port - 2000ml Corning Glass Bottle;
NOT IRRADIATED