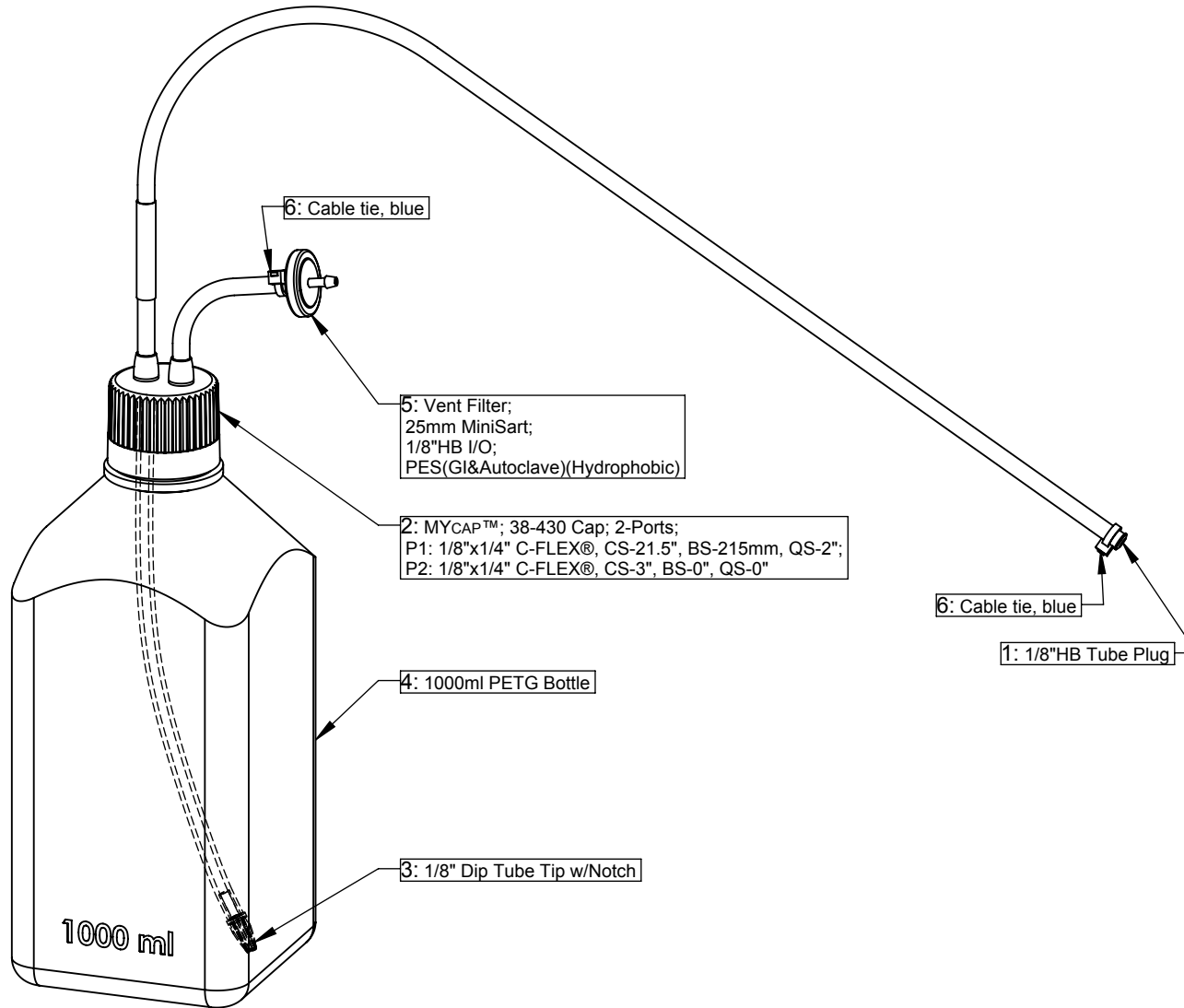


1000ml Nalgene Square Media Bottle, PETG



6: Cable tie, blue

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

2: MYCAP™; 38-430 Cap; 2-Ports;
P1: 1/8"x1/4" C-FLEX®, CS-21.5", BS-215mm, QS-2";
P2: 1/8"x1/4" C-FLEX®, CS-3", BS-0", QS-0"

4: 1000ml PETG Bottle

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

6: Cable tie, blue

1: 1/8"HB Tube Plug

Bill of Materials:

Item	Qty	AP Number	Material
1	1	AP-1181	-ADCF Polypropylene
2	1	AP-3091	-ADCF C-FLEX®(374) -ADCF Silicone
3	1	AP-2170	-ADCF Polypropylene
4	1	AP-1330	-PETG
5	1	AP-2767	-Polypropylene, PES
6	2	AP-0133	-Tefzel®

Revision History:

- 0 - Initial release
- 1 - P1 QS 1.5" to 2"
- Update drawing 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. 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 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 NGPG1000D

Drawing Number: MYCAP NGPG1000D

Revision: 1

Released By:

Release Date:

Title:

MYCAP™; 1000ml PETG; 2-Ports;
1/8"x1/4" C-FLEX® w/Dip Tube;
QS on inlet port