

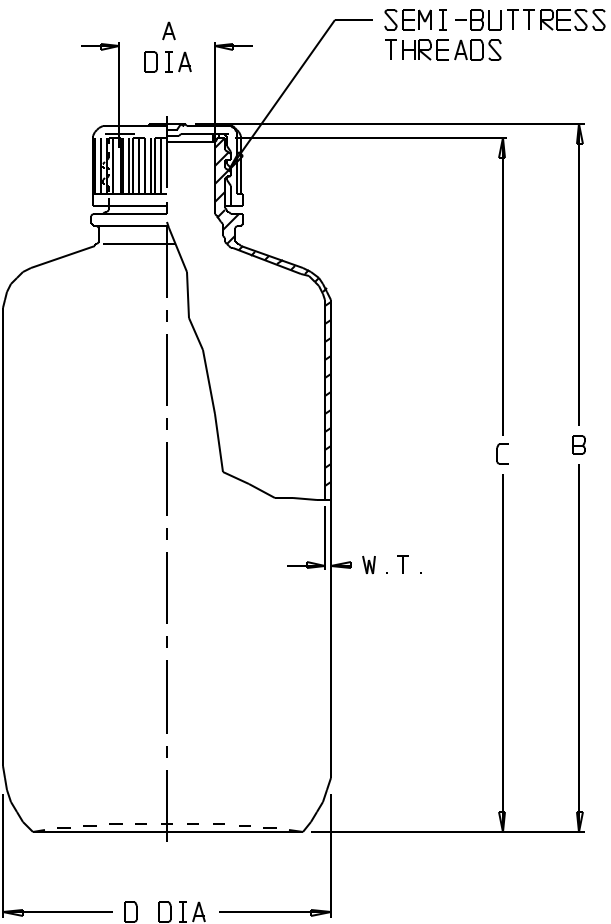
CATALOG NO.	NOMINAL CAPACITY, (oz)					MATERIALS	
						BOTTLE	CLOSURE
002097	8	16	32	64	1 GAL	FLPE	FLPP

SIZE CODE.....:	-0008		-0016		-0032	
SIZE (NOM VOL mL/oz):	250/8		500/16		1000/32	
NOM. BRIM CAP. (mL):	285		525		1045	
NECK FINISH (mm)....:	24-415		28-415		38-430	
THREAD (T.P.I.).....:	8		6		6	
NOM. WEIGHT W/CLOSURE (gms).....:	32		58		100	
	INCH	MM	INCH	MM	INCH	MM
A (BTL NECK I.D.)...:	0.69 ±0.02	17.5 ±0.5	0.84 ±0.02	21.3± 0.5	1.08 ±0.03	27.4 ±0.8
B (HT. W/CLOSURE)...:	5.22 ±0.08	132.6 ±2.0	6.71 ±0.09	170.4 ±2.3	8.50 ±0.10	215.9 ±2.5
C (HT. W/O CLOSURE):	5.13 ±0.08	130.3 ±2.0	6.59 ±0.09	167.4 ±2.3	8.35 ±0.10	212.1 ±2.5
D (DIA).....:	2.40 ±0.07	61.0 ±1.8	2.86 ±0.08	72.6 ±2.0	3.60 ±0.08	91.4 ±2.0
W.T.(NOM. WALL THK):	0.050	1.27	0.050	1.27	0.050	1.27

SIZE CODE.....:	-0005		-0010	
SIZE (NOM VOL mL/oz)	2000/64		4000/1gal	
NOM. BRIM CAP. (mL):	2100		4160	
NECK FINISH (mm)....:	38-430		38-430	
THREAD (T.P.I.).....:	6		6	
NOM. WEIGHT W/CLOSURE (gms).....:	160		370	
	INCH	MM	INCH	MM
A (BTL NECK I.D.)...:	1.08 ±0.03	27.4 ±0.8	1.02 ±0.03	25.9 ±0.8
B (HT. W/CLOSURE)...:	9.85 ±0.12	250.2 ±3.0	13.14 ±0.16	333.8 ±4.1
C (HT. W/O CLOSURE):	9.68 ±0.12	245.9 ±3.0	12.96 ±0.16	329.2 ±4.1
D (DIA).....:	4.74 ±0.08	120.4 ±2.0	6.02 ±0.10	152.9 ±2.5
W.T.(NOM. WALL THK):	0.050	1.27	0.050	1.27

NOTES:

- ALL DIMENSIONS ARE NOMINAL AND IN INCHES AND MILLIMETERS UNLESS OTHERWISE NOTED.
- THIS IS NOT A PRODUCTION DRAWING, BUT IS FOR INFORMATIONAL PURPOSES ONLY, AND IS SUBJECT TO CHANGE WITHOUT NOTICE.



SPECS ON BACK

NOT TO SCALE

CONFIDENTIAL

THIS PRODUCT SPECIFICATION SHEET CONTAINS CONFIDENTIAL AND/OR PROPRIETARY INFORMATION AND IS TO BE USED ONLY FOR THE EXPRESSED PURPOSES AGREED UPON BY AN AUTHORIZED OFFICER OF THERMO FISHER SCIENTIFIC.

ThermoFisher
SCIENTIFIC

Thermo Fisher Scientific
75 Panorama Creek Drive
Rochester, NY 14625-2385, USA

MADE	MG	APP'D	PVC	PSS 00-2097-XXXX-1	REV. 0499
CK'D	PVC		4/28/99		

DESCRIPTION:

FLUORINATED NARROW-MOUTH BOTTLES