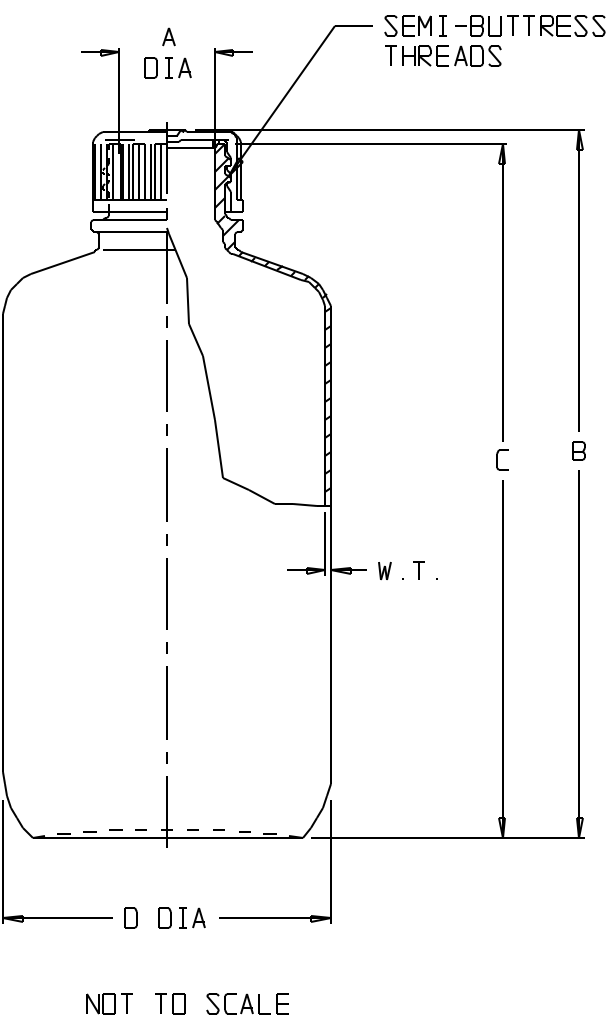


CATALOG NO.	NOMINAL CAPACITY, (oz)				MATERIALS	
					BOTTLE	CLOSURE
312097	4	8	16	32	FLPE	FLPP

SIZE CODE.....:	-0004		-0008	
SIZE (NOM VOL mL/oz):	125/4		250/8	
NOM. BRIM CAP. (mL):	140		285	
NECK FINISH (mm)....:	24-415		24-415	
THREAD (T.P.I.).....:	8		8	
NOM. WEIGHT W/CLOSURE (gms).....:	20		32	
	INCH	MM	INCH	MM
A (BTL NECK I.D.)...:	0.69 ±0.02	17.5 ±0.5	0.69 ±0.02	17.5 ±0.5
B (HT. W/CLOSURE)...:	3.97 ±0.08	100.8 ±2.0	5.22 ±0.08	132.6 ±2.0
C (HT. W/O CLOSURE):	3.88 ±0.08	98.6 ±2.0	5.13 ±0.08	130.3 ±2.0
D (DIA).....:	1.98 ±0.06	53.3 ±1.5	2.40 ±0.07	61.0 ±1.8
W.T.(NOM. WALL THK):	0.040	1.02	0.050	1.27

SIZE CODE.....:	-0016		-0032	
SIZE (NOM VOL mL/oz):	500/16		1000/32	
NOM. BRIM CAP. (mL):	525		1045	
NECK FINISH (mm)....:	28-415		38-430	
THREAD (T.P.I.).....:	6		6	
NOM. WEIGHT W/CLOSURE (gms).....:	58		100	
	INCH	MM	INCH	MM
A (BTL NECK I.D.)...:	0.84 ±0.02	21.3± 0.5	1.08 ±0.03	27.4 ±0.8
B (HT. W/CLOSURE)...:	6.71 ±0.09	170.4 ±2.3	8.50 ±0.10	215.9 ±2.5
C (HT. W/O CLOSURE):	6.59 ±0.09	167.4 ±2.3	8.35 ±0.10	212.1 ±2.5
D (DIA).....:	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

1. ALL DIMENSIONS ARE NOMINAL AND IN INCHES AND MILLIMETERS UNLESS OTHERWISE NOTED.
2. 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	PB	APP'D	PVC	PSS 312097-XXXX-3	REV. 0203
CK'D	PVC	2/25/03			
DESCRIPTION: FLUORINATED NARROW-MOUTH BOTTLES					