

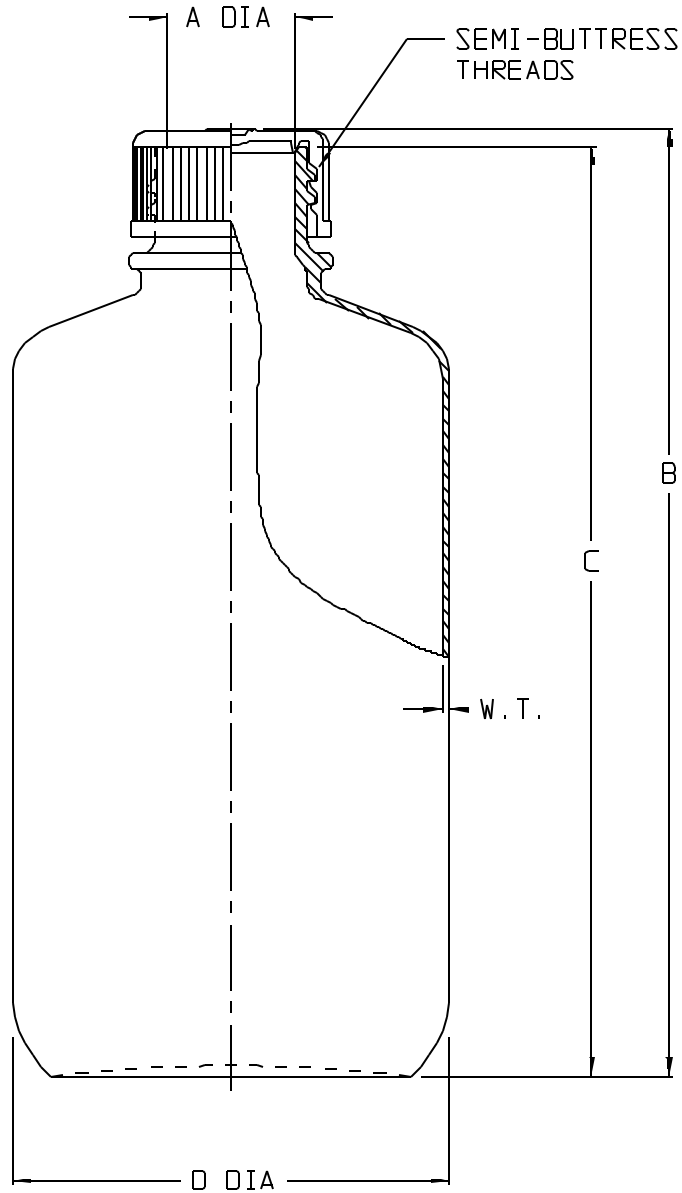
CATALOG NO.	NOMINAL CAPACITY, (oz)						MATERIALS	
							BOTTLE	CLOSURE
312084	1	2	4	8	16	32	TRANSLUCENT AMBER HDPE	OPAQUE AMBER PP
312087	1	2	4	8	16	32	PPCO	PP
312088	1	2	4	8	16	32	LDPE	PP
312089	1	2	4	8	16	32	HDPE	PP

SIZE CODE.....	-0001		-0002		-0004	
SIZE (NOM VOL mL/oz)	30/1		60/2		125/4	
NOM. BRIM CAP. (mL)	34		64		143	
NECK FINISH (mm)...	20-415		20-415		24-415	
THREAD (T.P.I.).....	8		8		8	
NOM. WEIGHT W/CLOSURE (gms).....	8		12		16	
	INCH	MM	INCH	MM	INCH	MM
A (NECK I.D.).....	0.54 ±0.02	13.7 ±0.5	0.54 ±0.02	13.7 ±0.5	0.69 ±0.02	17.5 ±0.5
B (HT. W/CLOSURE)...	2.40 ±0.07	61.0 ±1.8	3.36 ±0.08	85.3 ±2.0	3.97 ±0.08	100.8 ±2.0
C (HT. W/O CLOSURE):	2.30 ±0.07	58.4 ±1.8	3.25 ±0.08	82.6 ±2.0	3.87 ±0.08	98.3 ±2.0
D (DIA).....	1.35 ±0.05	34.3 ±1.3	1.52 ±0.06	38.61 ±1.5	1.99 ±0.06	50.5 ±1.5
W.T.(NOM. WALL THK):	0.035	0.89	0.035	0.89	0.030	0.76

SIZE CODE.....	-0008		-0016		-0032	
SIZE (NOM VOL mL/oz)	250/8		500/16		1000/32	
NOM. BRIM CAP. (mL)	289		555		1090	
NECK FINISH (mm)...	24-415		28-415		38-430	
THREAD (T.P.I.).....	8		6		6	
NOM. WEIGHT W/CLOSURE (gms).....	26		40		78	
	INCH	MM	INCH	MM	INCH	MM
A (NECK I.D.).....	0.69 ±0.02	17.5 ±0.5	0.84 ±0.02	21.3 ±0.5	1.080 ±0.025	27.4 ±0.6
B (HT. W/CLOSURE)...	5.23 ±0.08	132.8 ±2.0	6.73 ±0.09	170.9 ±2.3	8.51 ±0.10	216.2 ±2.5
C (HT. W/O CLOSURE):	5.14 ±0.08	130.6 ±2.0	6.61 ±0.09	167.9 ±2.3	8.37 ±0.10	212.6 ±2.5
D (DIA).....	2.40 ±0.07	61.0 ±1.8	2.87 ±0.08	72.9 ±2.0	3.60 ±0.08	91.4 ±2.0
W.T.(NOM. WALL THK):	0.030	0.76	0.040	1.02	0.035	0.89

2. This is not a production drawing, but is for informational purposes only,
and is subject to change without notice.

1. All dimensions are nominal and in inches and millimeters unless otherwise noted.



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 CK'D	MCC DM	APP'D GRP	PSS 31-208X-XXXX-3	REV. 1107
DESCRIPTION: NARROW MOUTH, ROUND PACKAGING BOTTLES				