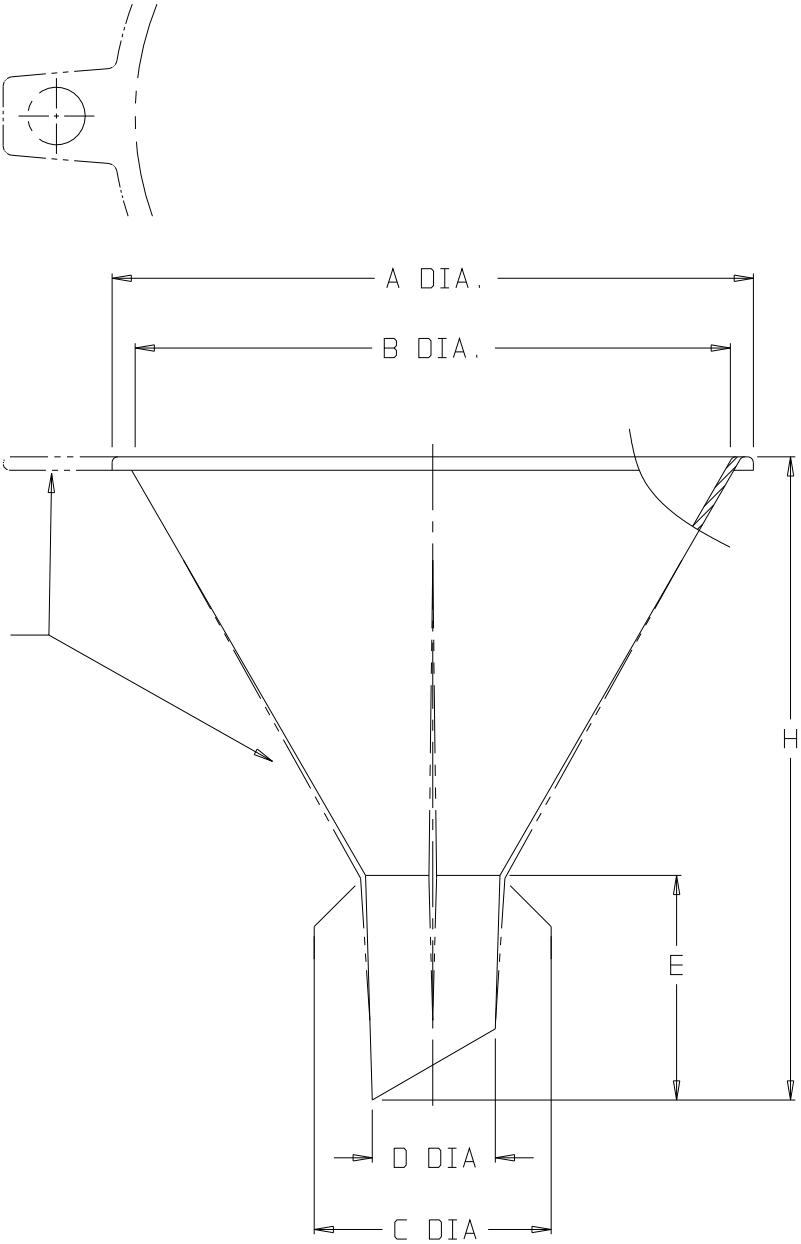


CATALOG NUMBER	4260-0040		4260-0060		4260-0100	
MATERIAL	LDPE		LDPE		LDPE	
NOM WT (gms)	52		154		385	
	INCH	MM	INCH	MM	INCH	MM
A DIA	4.34 ±0.04	110.2 ±1.0	6.6 ±0.1	167 ±3	10.5 ±0.3	267 ±8
B (TOP I.D.)	3.96 ±0.04	100.6 ±1.0	5.9 ±0.1	150 ±3	9.9 ±0.3	251 ±8
C DIA	0.84 ±0.02	21.3 ±0.5	0.69 ±0.02	17.5 ±0.5	1.11 ±0.04	28.2 ±1.0
D DIA	0.46 ±0.02	11.7 ±0.5	0.62 ±0.02	15.7 ±0.5	0.91 ±0.02	23.1 ±0.5
E (STEM HT)	2.84 ±0.04	72.1 ±1.0	3.00 ±0.04	76.2 ±1.0	4.60 ±0.04	116.8 ±1.0
H (OVERALL HT)	5.8 ±0.1	148 ±3	7.7 ±0.1	196 ±3	12.5 ±0.3	318 ±8

NOTES:

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.



NOT TO SCALE

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	PMP	APP'D	PWC	PSS 4260-XXXX-1	REV. 0516
CHK'D	JC	5/3/16			

DESCRIPTION:

HEAVY DUTY FUNNELS