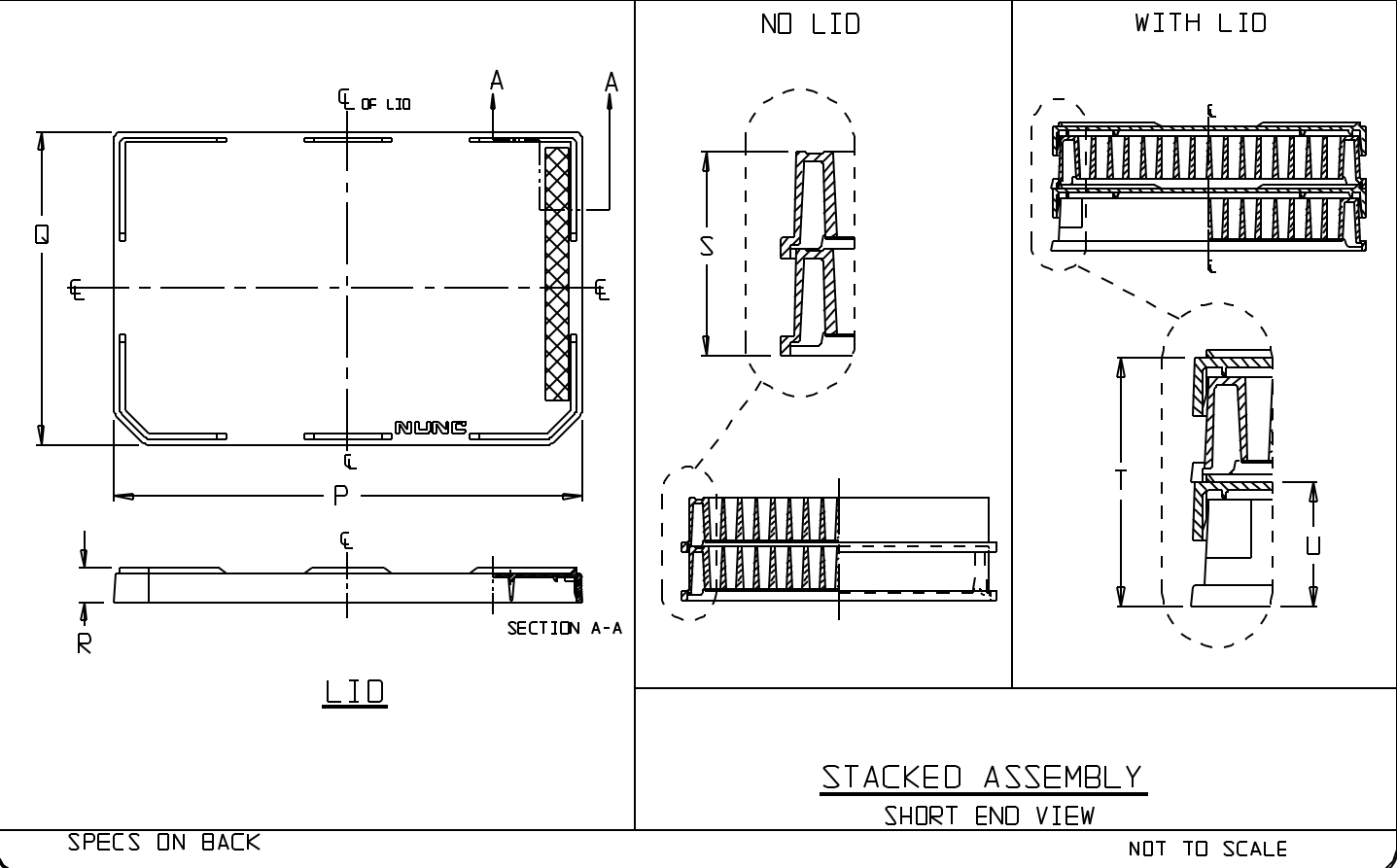
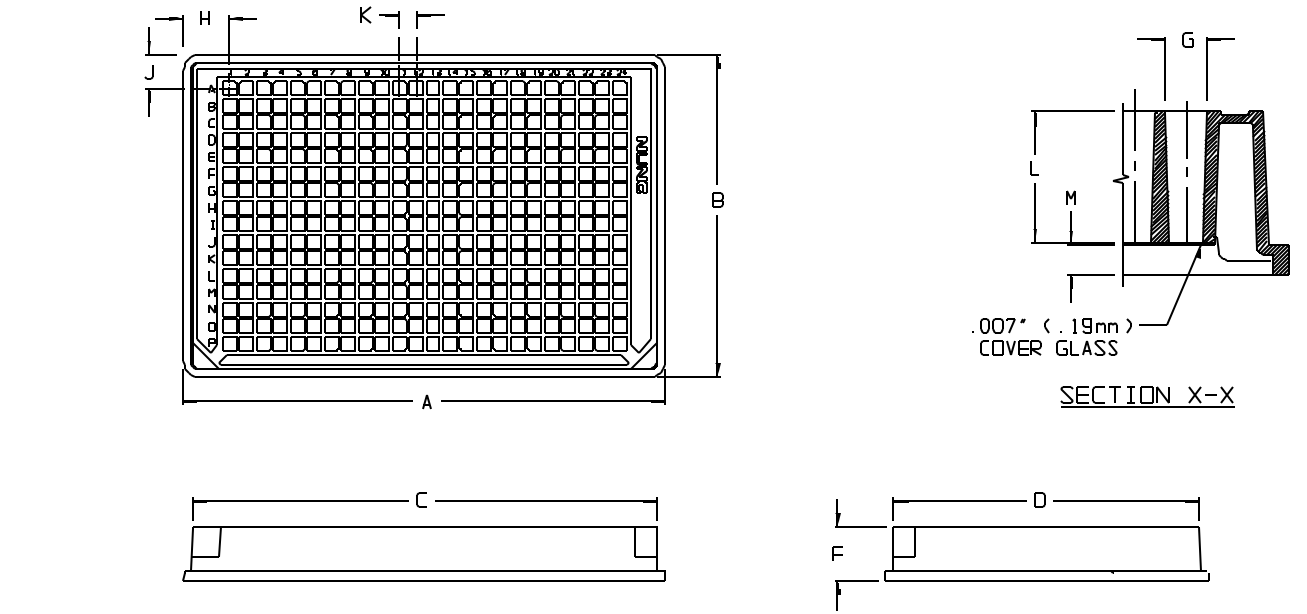


CATALOG NUMBER.....	SEE BELOW	
NOM. CAP. (μL).....	120	
MATERIAL.....	PS	
	INCH	MM
A (BASE LENGTH).....	5.03	127.8
B (BASE WIDTH).....	3.37	85.5
C (TOP LENGTH).....	4.87	123.7
D (TOP WIDTH).....	3.20	81.3
F (OVERALL HEIGHT).....	0.57	14.4
G (WELL WIDTH).....	0.15	3.7
H (A-1 LOCATION) (X).....	0.48	12.1
J (A-1 LOCATION) (Y).....	0.35	9.0
K (WELL-TO-WELL SPACING)....	0.18	4.5
L (WELL DEPTH).....	0.46	11.6
M (BOTTOM OF GLASS).....	.10	2.5
*P (LID LENGTH).....	5.02	127.4
*Q (LID WIDTH).....	3.36	85.3
*R (LID HEIGHT).....	.36	9.1
S (STACKED HEIGHT - NO LID):	1.09	27.7
*T (STACKED HEIGHT W/LIDS):	1.33	33.8
*U (ASSEMBLY HEIGHT W/LID):	.665	16.9

CATALOG NUMBER	SURFACE	COLOR	STERILE	WITH LID*	UNITS/ PK/CASE
164586	CELL CULTURE	BLACK	YES	YES	6/30
240074	NON-TREATED	BLACK	NO	NO	5/30

* DIM.S APPLY ONLY WHERE "WITH LID" = YES, USING NUNC LID #8337.



- 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.

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	JDD	APP'D	PVC	PSS 16458X	REV. 0505
CHK'D	RDR		6/8/05		
DESCRIPTION: 384 WELL, PS, GLASS BOTTOM PLATE					