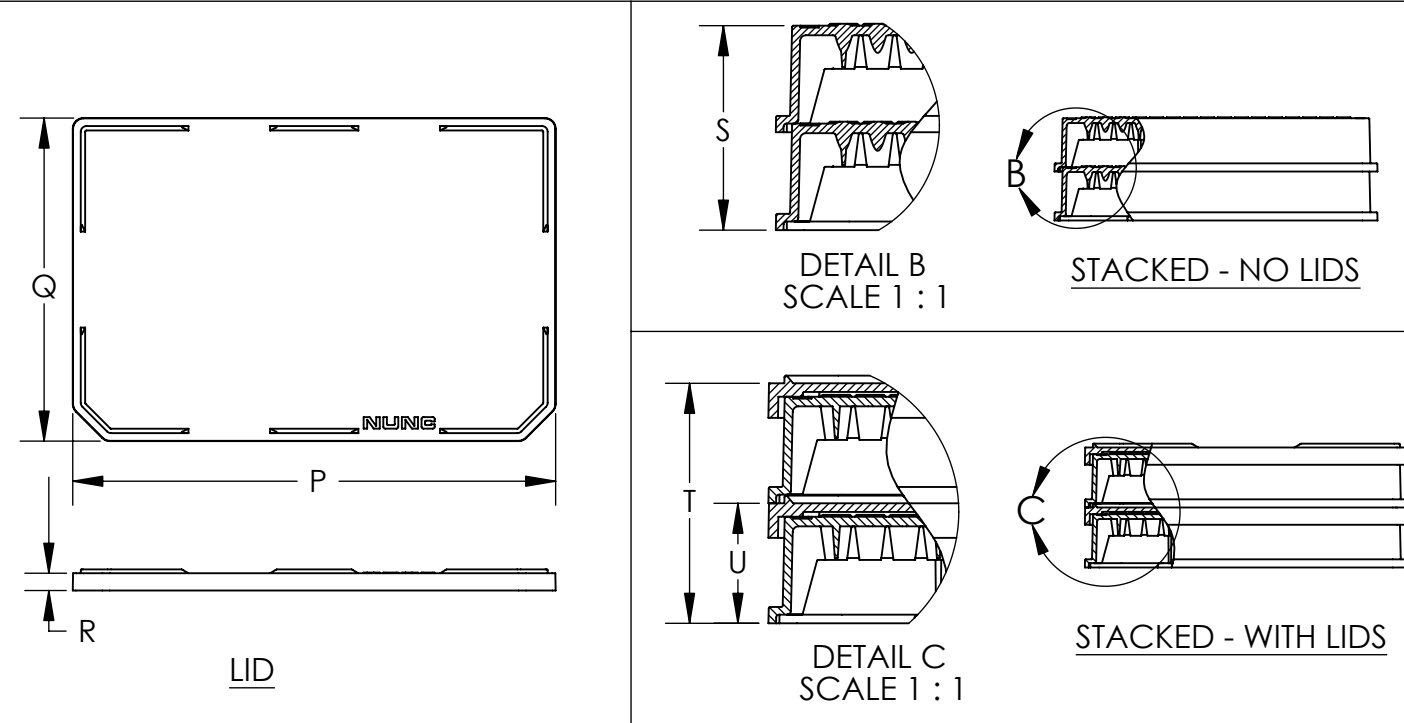
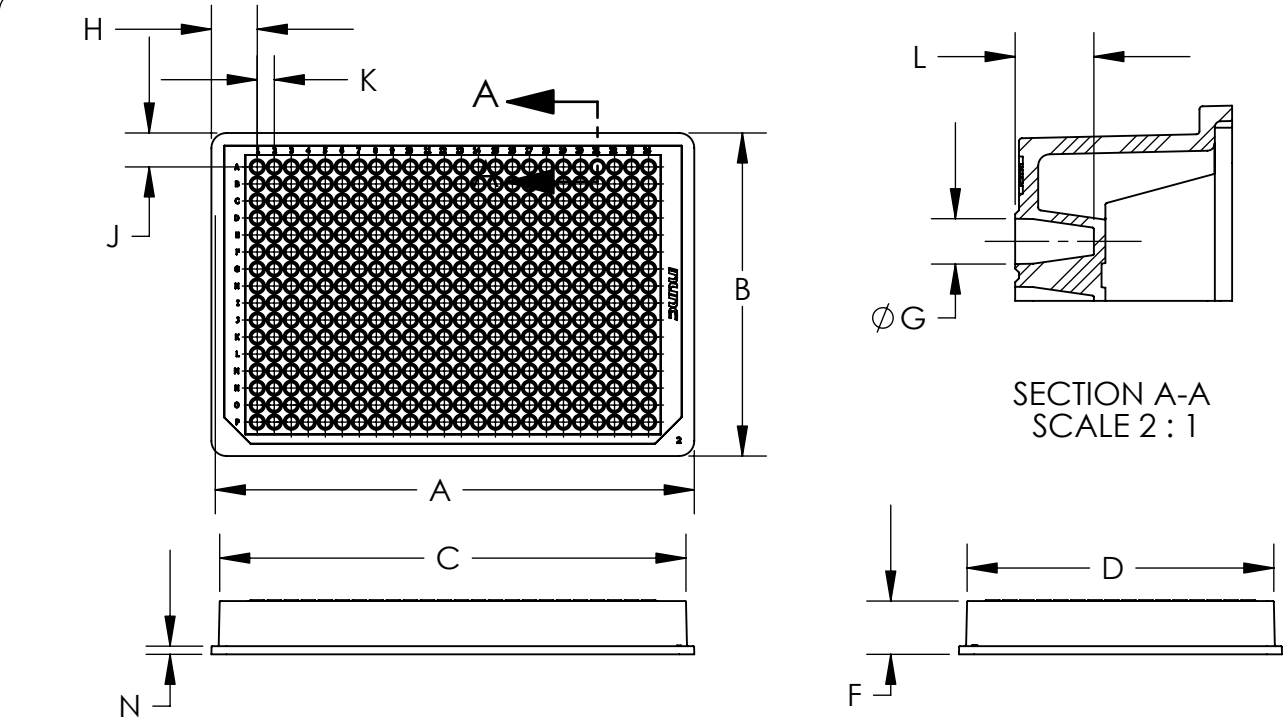


CATALOG NUMBER.....:	SEE CHART BELOW	
NOM. CAP. (uL).....:	25uL	
MATERIAL (BASE).....:	PS	
MATERIAL (LID*).....:	PS	
	INCH	MM
A (BASE LENGTH).....:	5.03	127.8
B (BASE WIDTH).....:	3.36	85.3
C (TOP LENGTH).....:	4.86	123.4
D (TOP WIDTH).....:	3.19	81.0
F (OVERALL HEIGHT).....:	0.57	14.4
G DIA.....:	0.12	3.0
H (A-1 LOCATION)(X).....:	0.47	11.9
J (A-1 LOCATION)(Y).....:	0.35	9.0
K (WELL-TO-WELL SPACING).....:	0.18	4.5
L (WELL DEPTH).....:	0.20	5.1
N (FLANGE SHORT SIDE).....:	0.09	2.2
*P (LID LENGTH).....:	5.03	127.8
*Q (LID WIDTH).....:	3.37	85.5
*R (LID HEIGHT).....:	0.19	4.8
S (STACKED HEIGHT - NO LID).....:	1.08	27.4
*T (STACKED HEIGHT W/LIDS).....:	1.27	32.3
*U (ASSEMBLY HEIGHT W/LID).....:	0.63	15.9

CATALOG NO.	SURFACE	COLOR	STERILE	W/LID*	UNITS/(PK/CASE)
164701	CELL CULTURE	CLEAR	YES	YES	100
164702	CELL CULTURE	BLACK	YES	YES	100
164703	CELL CULTURE	WHITE	YES	YES	100
264704	NON-TREATED	CLEAR	NO	NO	100
264705	NON-TREATED	BLACK	NO	NO	100
264706	NON-TREATED	WHITE	NO	NO	100

* DIM.S APPLY ONLY WHERE "W/LID" = "YES" USING NNI LID. #1002093

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	JDD	APP'D	PVC	PSS 1647XX	REV.	0906
CK'D	MP	11/3/06				
DESCRIPTION:						
384-SHALLOW WELL, STANDARD HEIGHT						