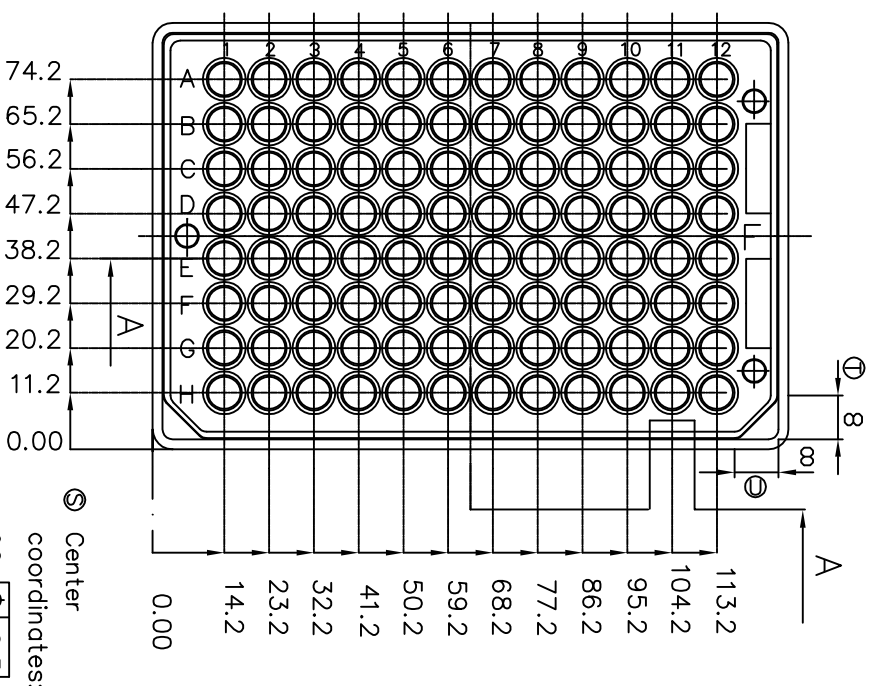


www.nunckbrand.com

DRYING BY DATE:	CHANGE ORDER NO.:	VALID FROM:	APPROVED BY DATE:	FORMAT:
TBS/30.11.99	CC-1763	2007.06.04	AMU 2007.05.30	A3
ITEM NUMBERS:	MATERIAL:	DRAWING NO.:	VERSION:	PAGE TYPE:
-	PS	2238	10	-C
				SCALE:
				1:1 (2:1)
				METRIC