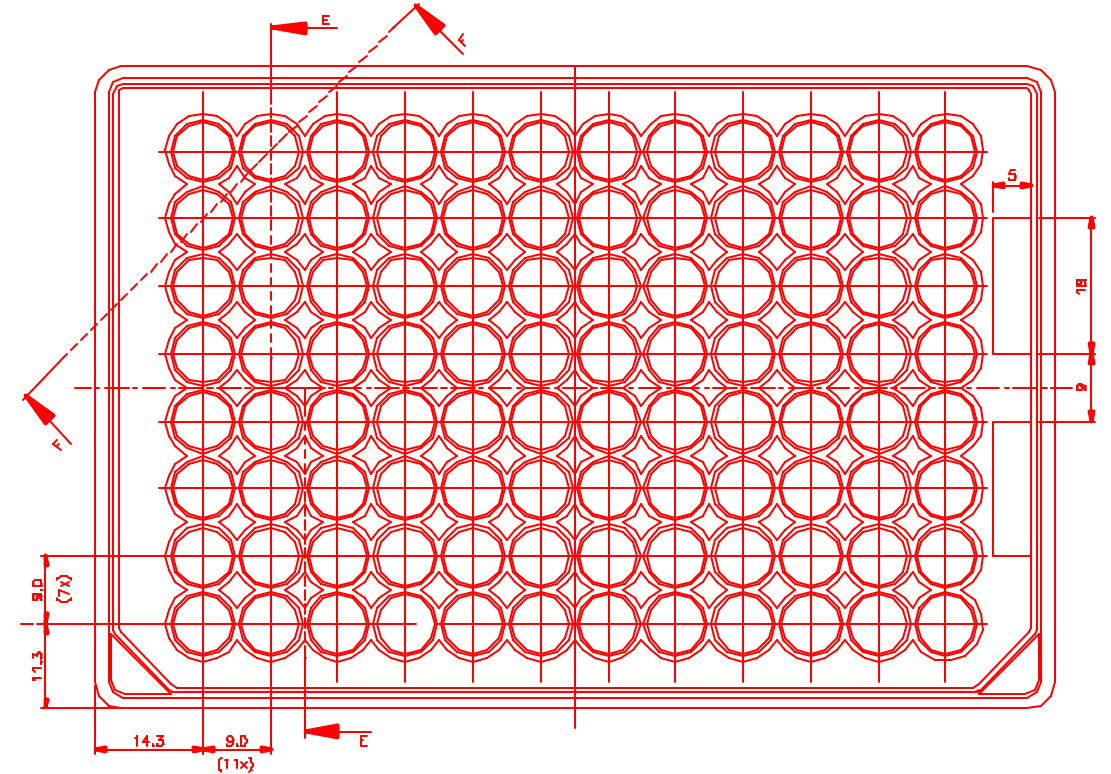


SECTION EE
SCALE 2:1



Tolerances unless otherwise specified:
 Without Decimal: $\pm 0.5\text{mm}$
 1 Place Decimal: $\pm 0.2\text{mm}$
 Material: Polypropylene Homopolymer PF511 or PD702 or Equivalent
 PLEASE NOTE that dimns. 127.7x85.8 (footprint) is as measured near to the corners.

UNLESS OTHERWISE SPECIFIED DIMENSIONS ARE IN MM TOLERANCES ARE: ANGLES: \pm DECIMAL: \pm FRACTIONS: \pm		DWG TITLE DEEP WELL PLATE 2mil	
APPROVALS	DATE		
DRAWN: CRN	07/12/00		
CHECKED: D.TAUNK	07/13/00		
BY: [Signature]		SIZE: D	DWG. NO.: SLD079
file: e:\drgs\app\pdrge\sl0078AA		SCALE: 1:1, 2:1, 8:1	REV: AA
		SHEET 1 OF 1	