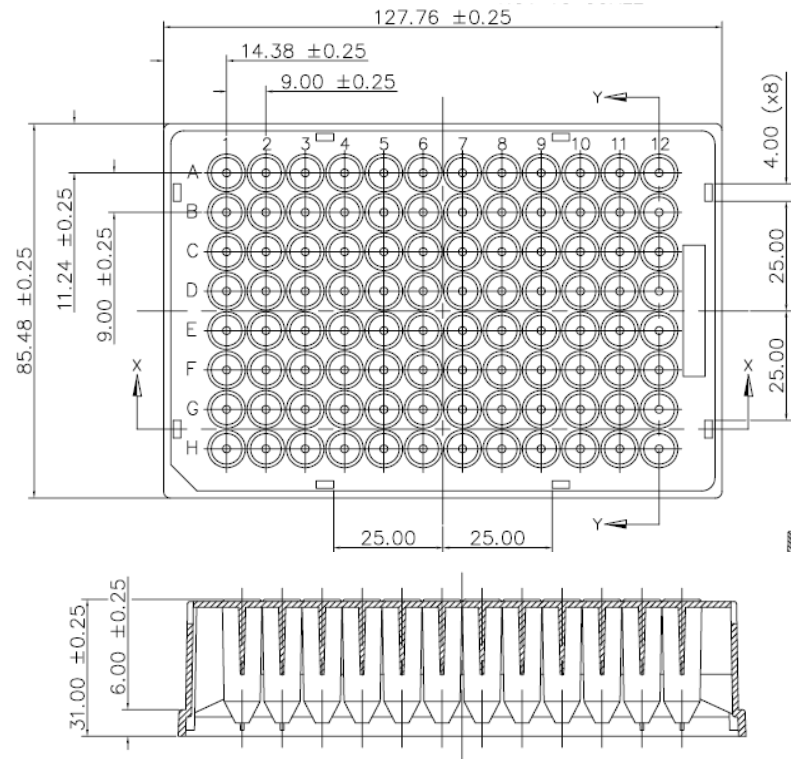


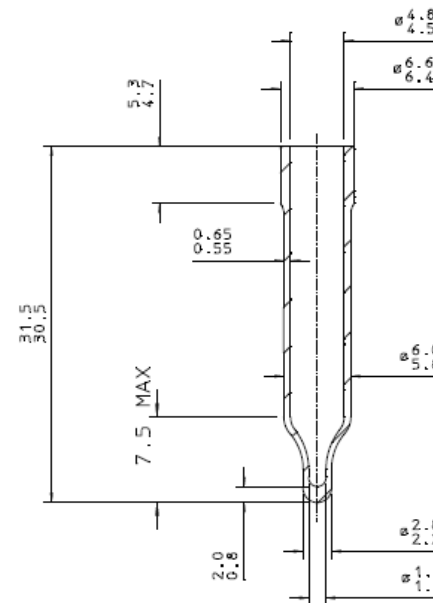
**General Arrangement Drawings
Plates for Chromatography-G3-60180-K100**

Cat.No.	Description	A	B	F	G1	H	J	K	L	M
		BASE LENGTH	BASE WIDTH	OVERALL HEIGHT	DIA TOP	A - I LOCATION (X)	A - I LOCATION (Y)	WELL TO WELL SPACING	WELL DEPTH	BOTTOM OF WELL DISTANCE
60180-K100	96 Well small Volume Microplate with Sealing Mat	127.76	85.48	35.00 (vial)	4.65 (vial)	14.38	11.24	9.00	27.00 (30.00 vial)	4.0

60180-K100



60180-V100



Notes.

Drawings are for instrument set-up purposes and are not manufacturing drawings

Thermo Fisher Scientific is providing information that has been supplied by the original manufacturer and is not responsible for subsequent changes or modifications.

All dimensions mm