

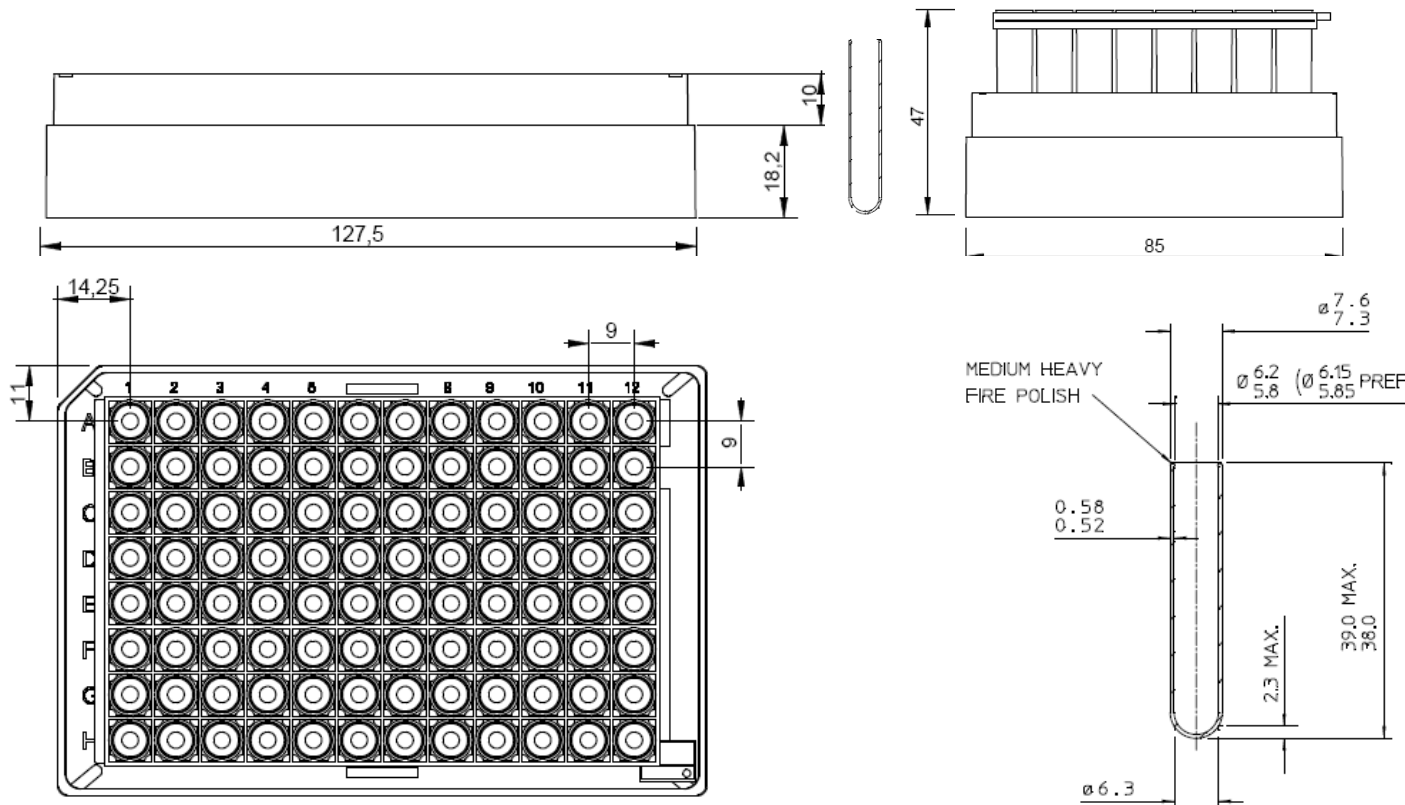
**General Arrangement Drawings
Deepwell Plates for Chromatography-G3**

Cat.No.	Description	A	B	F	G1	H	J	K	L	M
60180-K104	WEBSEAL KIT, 96 CLR, 1.1ML, ROUND, T/S MAT, 5/PK	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
		127.5	85	47	6.0	14.25	11	9	38	4

60180-P401

60180-K104

60180-V105



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