

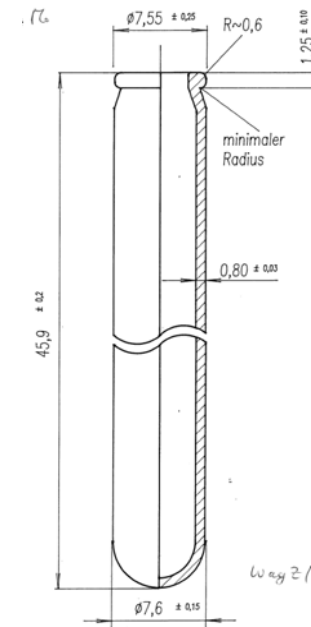
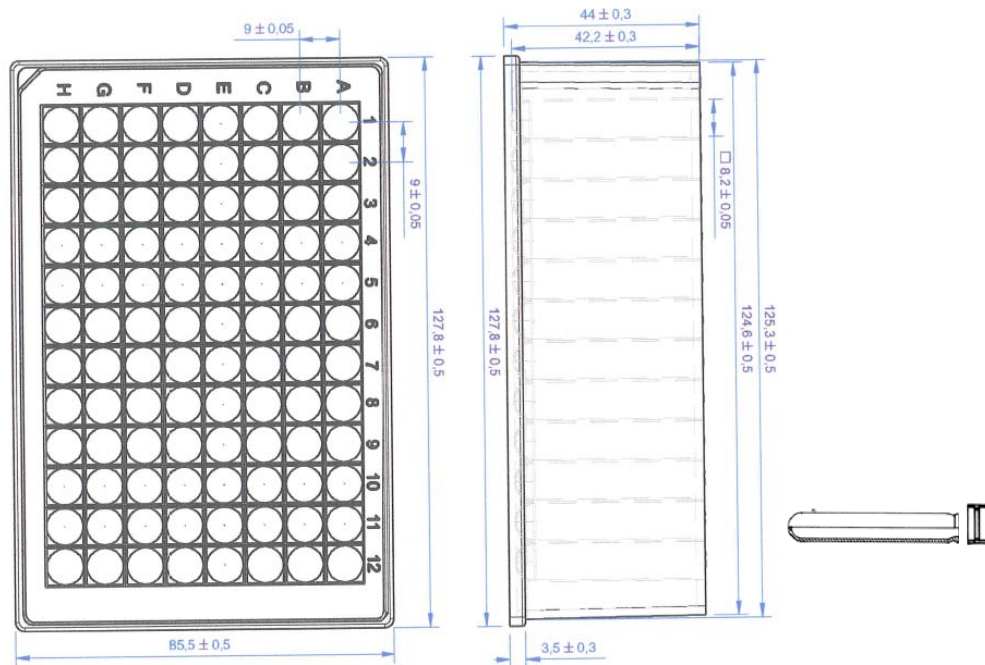
**General Arrangement Drawings
Deepwell Plates for Chromatography- Glass Vials-Pre-inserted**

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-K202	SQW-Block, PP, 96 positions, filled with 1ml glass Micro-Insert, with assembled 9mm PE Cap	127.8	85.5	48.5	7.55	14.00	11.00	9.0	45.1	3.5
60180-K203	SQW-Block, PP, 96 positions, filled with 1ml glass Micro-Insert, with assembled 9mm PE Cap with slit.	127.8	85.5	48.5	7.55	14.00	11.00	9.0	45.1	3.5

60180-P406

60180-K202/K203

60180-V110



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