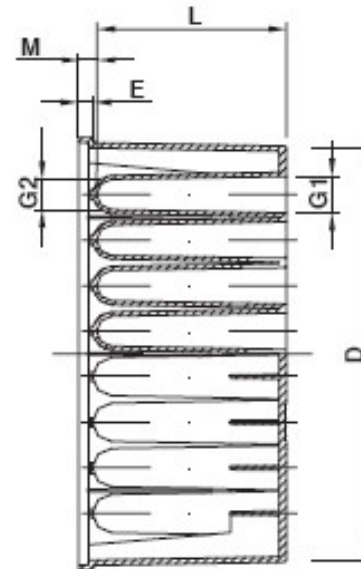
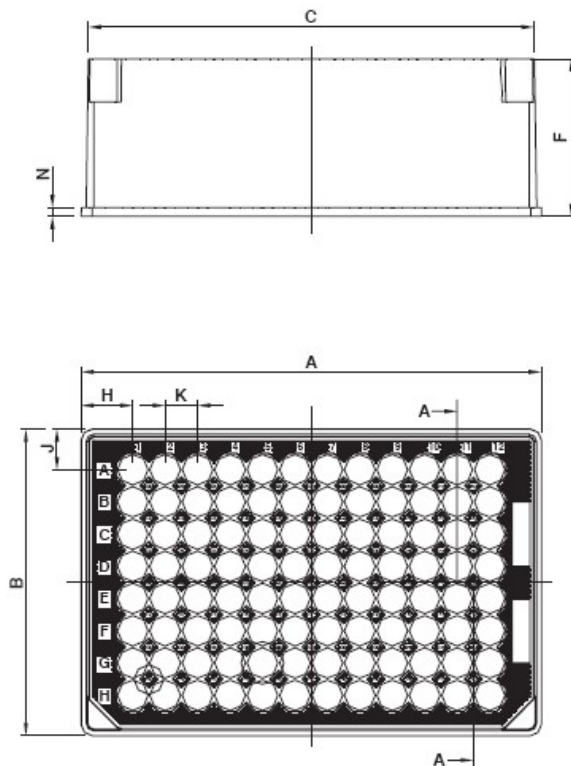


**General Arrangement Drawings
Deepwell Plates for Chromatography-Round Well**

Cat.No.	Description	A	B	C	D	E	F	G1	G2	H	J	K	L	M	N
60180-P103 60180-P133	96-Well MicroWell Microplate, round well- 1.3mL	127.8	85.5	125.4	84.3	2.5	31.6	Ø8.4	Ø7.2	14.4	11.2	9	29.1	1.4	2.5
60180-P104 60180-P134	96-Well MicroWell Microplate, round well- 2.0mL	127.8	85.5	123.6	81.6	2.5	43.6	Ø8.5	Ø7.2	14.3	11.3	9	41.5	1.4	2.5

60180-P103/60180-P133/60180-P104/60180-P134



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.