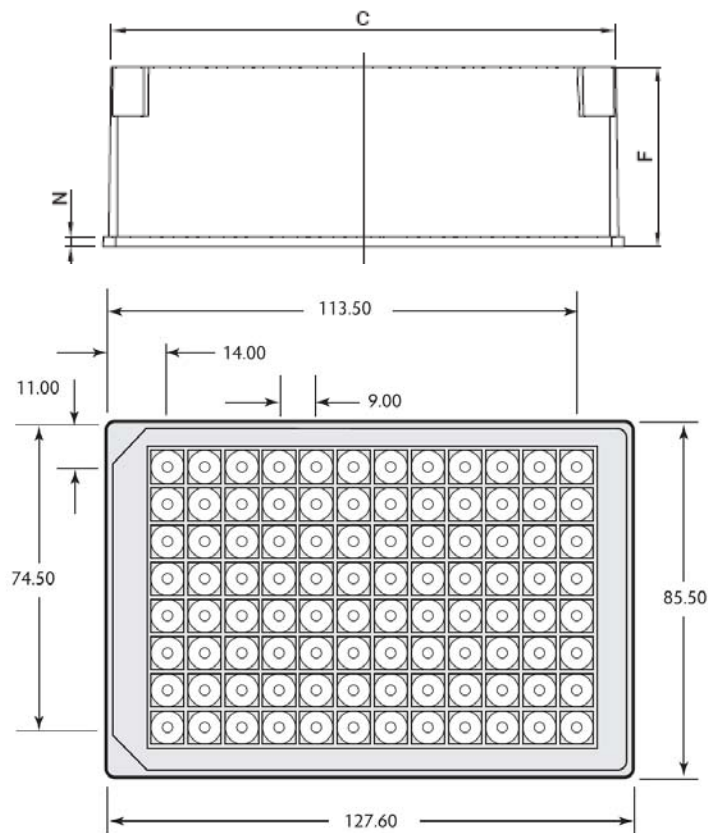


**General Arrangement Drawings  
Deepwell Plates for Chromatography-Square Well-Certified**

Cat.No.	Description	A	B	C	D	F	G1	H	J	K	L	M	N
		BASE LENGTH	BASE WIDTH	TOP LENGTH	TOP WIDTH	OVERALL HEIGHT	DIA TOP	A - I LOCATION (X)	A - I LOCATION (Y)	WELL TO WELL SPACING	WELL DEPTH	BOTTOM OF WELL DISTANCE	FLANGE SHORT SIDE
60180-P202 60180-P232	96-Well MicroWell Microplate, square well-2.0mL	127.60	85.50	125.4	84.3	44.40	8.30	14.00	11.00	9	39.90	4.5	4.0

**60180-P202/60180-P232**



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