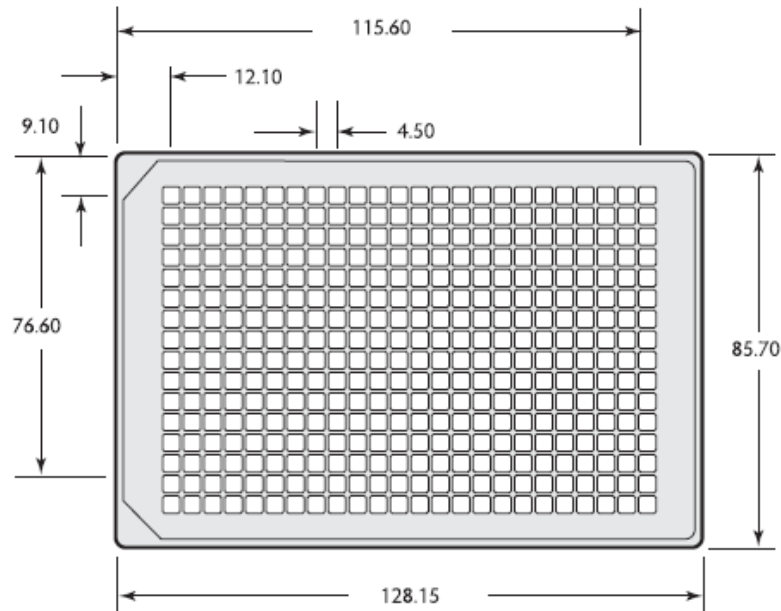
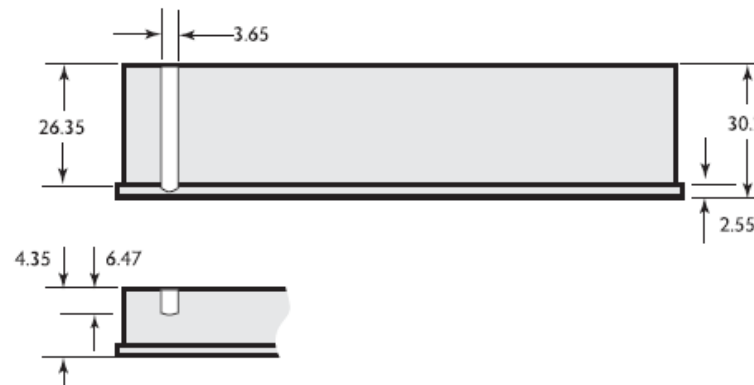


**General Arrangement Drawings  
Microplates for Chromatography-Certified**

Cat.No.	Description	A	B	F	G1	H	J	K	L	M	N
		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	FLANGE SHORT SIDE
60180-P203 60180-P213	384-Well MicroWell Microplate	128.15	85.70	30.20	3.65	12.10	9.10	4.50	6.47	3.85	2.55
60180-P204 60180-P214	384-Well Deep Well Microplate	128.15	85.70	30.20	3.65	12.10	9.10	4.50	26.25	3.85	2.55
60180-P205 60180-P215	384-Well MicroWell Microplate-V bottom	128.15	85.70	30.20	3.65	12.10	9.10	4.50	10.50	3.85	2.55



**60180-P204/60180-P214**



**60180-P203/60180-P213 60180-P205/60180-P215**

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