

**General Arrangement Drawings
Plate+ Glass Coated Microplates**

Cat.No.	Description	A	B	C	D	D'	E	F	G	H	I
60180-P302	Plate + Microplate 96 V	127.76	85.48	14.6	6.96	6.65	2.5	10.9	14.38	11.24	9
60180-P303											
60180-P300	Plate + Microplate 96 U	127.76	85.48	14.6	6.96	N/A	2.5	10.9	14.38	11.24	9
60180-P301											
60180-P310	Plate + Microplate 384	127.76	85.48	14.4	3.80	3.60	2.5	11.5	12.13	8.99	5
60180-P311											

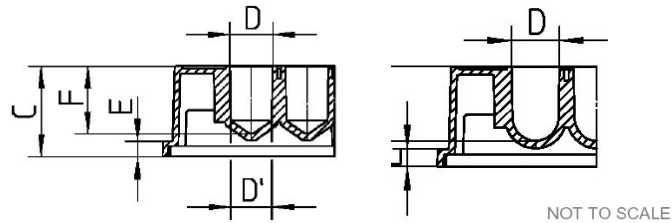
Notes.

The glass coating will have some effect on dimensions D, D' and F and these are shown as nominal values.

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.

60180-P302 / 60180-P303



60180-P300 / 60180-P301

