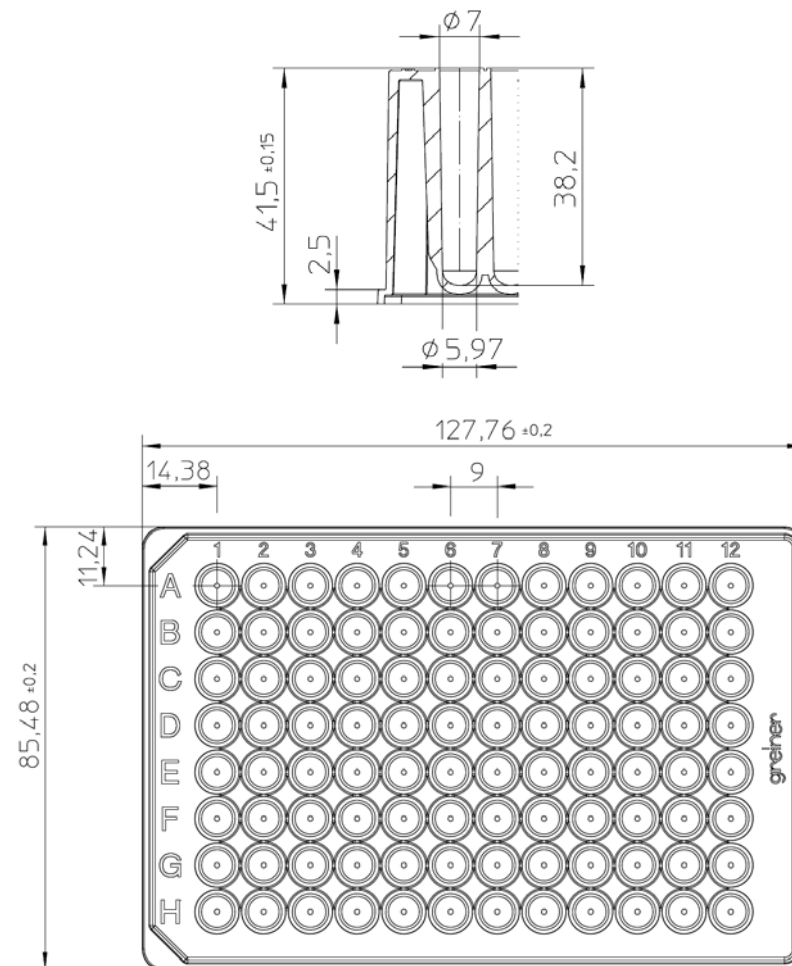


**General Arrangement Drawings**  
**Plate+ Glass Coated Microplates**

Cat.No.	Description	A	B	C	D	D'	E	F	G	H	I
60180-P306 60180-P307	Plate + Microplate 1mL	127.76	85.48	41.5	7	5.97	2.5	38.2	14.38	11.24	9

Dimensions in mm



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.