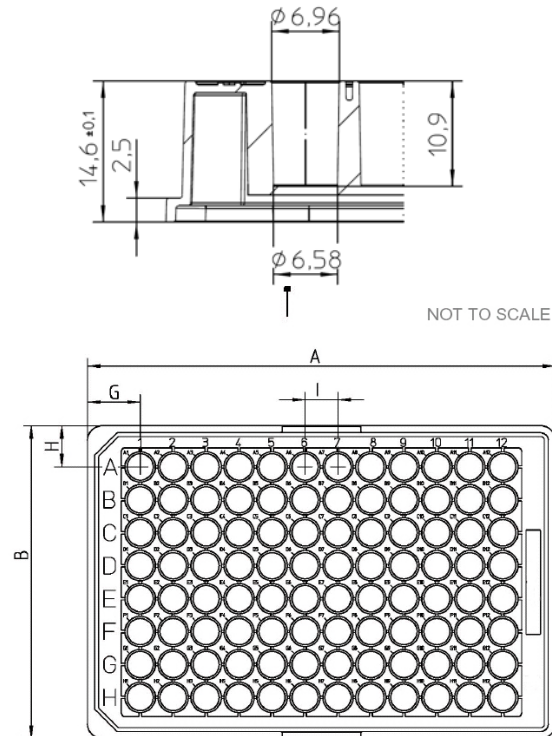


**General Arrangement Drawings**  
**Plate+ Glass Coated Microplates**

Cat.No.	Description	A	B	C	D	D'	E	F	G	H	I
60180-P304	Plate + Microplate 96-										
60180-P305	370 $\mu$ L	127.76	85.48	14.6	6.96	6.58	2.5	10.9	14.38	11.24	9

Dimensions in mm

**60180-P304 / 60180-P305**



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.