



A	B	C	D	D'	E	F	G	H	I
127.76±0.2	85.48±0.2	14.4	6.96	6.58	2.5	10.9	14.38	11.4	9.0

\*all sizes in millimeters (mm)