

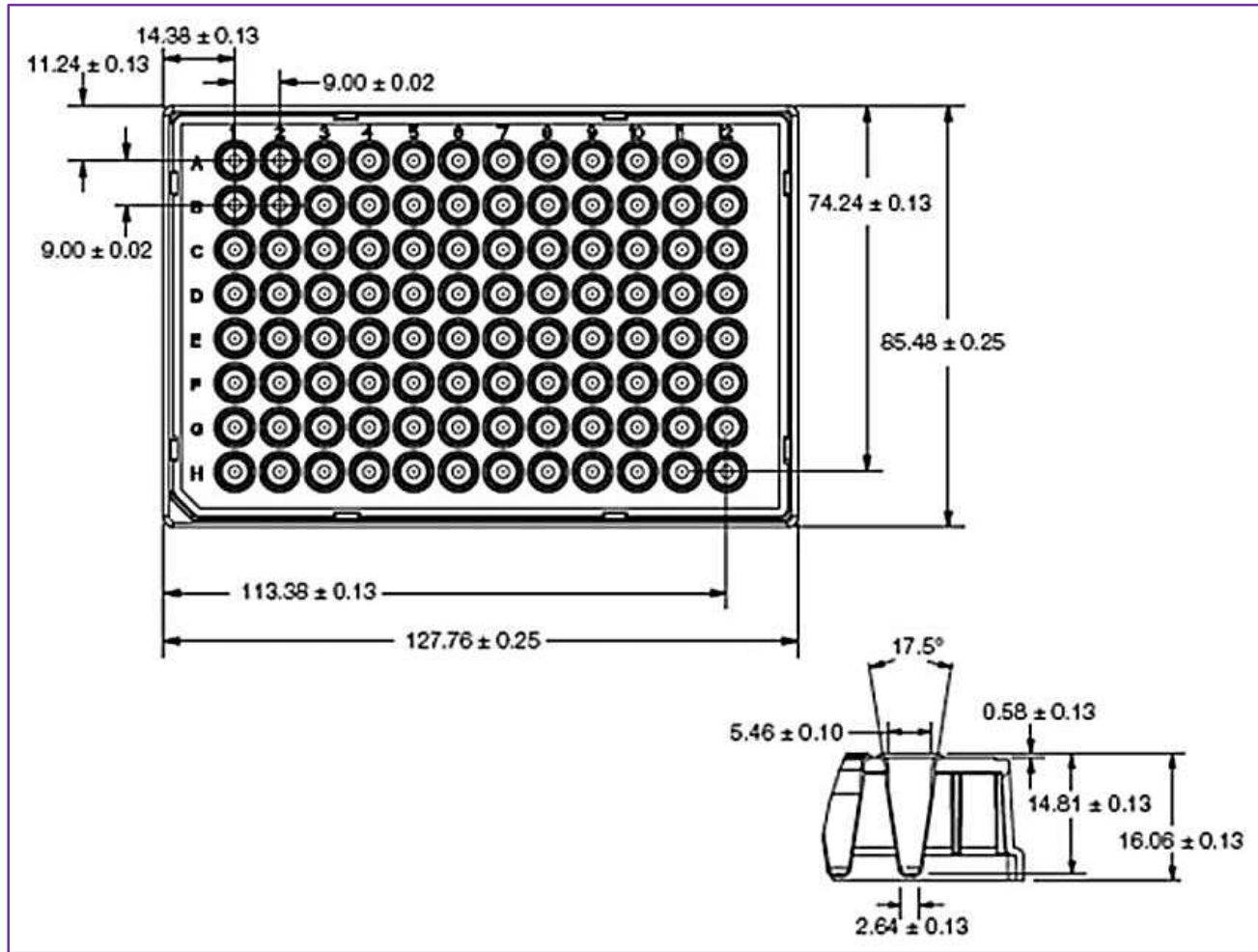
Research use only. Not for use in diagnostic procedures.

Technical datasheet:

Dimensions of Hardshell PCR Plate-96 Blue

- Part numbers:** PN 6008870 Blue rigid frame with clear wells
PN 6008610 Blue rigid frame with white wells
- Material and treatment:** Patented, rigid two-component design 96-well PCR microplate. Blue frame & clear wells (6008870) or blue frame & white wells (6008610). Low-profile, thin-wall (polypropylene tubes). Free of DNase, RNase and human genomic DNA. Well volume 200 µL.
- Packaging:** Part numbers 6008870 and 6008610 are cases with 50 plates, packed as 5 bags of 10 plates each.

Dimensions -- Customer Drawing:



Hard-Shell plates are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. patent Nos. 7,347,977, and 6,528,302