

Research use only. Not for use in diagnostic procedures.

Dimensions of ProxiPlate-96

- Part numbers:** PN 6006290 / 6006299 White, untreated
 PN 6006270 / 6006279 Black, untreated
- Material and treatment:** Black or white polystyrene microplate with 96 flat-bottom, shallow wells. Plate is available in untreated version only.
- Packaging:** 6006290 / 6006270 is a case with 50 plates, packed as 2 bags of 25 plates each. 6006299 / 6006279 holds 200 plates as 8 bags of 25.

Dimensions -- Customer Drawing

