

Research use only. Not for use in diagnostic procedures.

Technical datasheet:

Dimensions of polypropylene pipetting reservoir with 96 wells (2.5 mL/Well)

- Part numbers:** PN 6008880 Natural polypropylene, untreated
- Material and treatment:** Pipetting reservoir with 96 wells of 2.5 mL volume each. With pyramid-shaped bottom for efficient aspiration. Made of natural, untreated polypropylene.
- Packaging:** Part number 6008880 is a case with 25 reservoirs, packed as 5 bags of 5 reservoirs each.

Dimensions -- Customer Drawing:

