

Research use only. Not for use in diagnostic procedures.

Technical datasheet:

Dimensions of polypropylene pipetting reservoir with 4 columns (73 mL/Well)

- Part numbers:** PN 6008710 Natural polypropylene, untreated
- Material and treatment:** Pipetting reservoir with 4 column-wise wells of 73 mL volume each. Made of natural, untreated polypropylene.
- Packaging:** Part number 6008710 is a case with 25 reservoirs, packed as 5 bags of 5 reservoirs each.

Dimensions -- Customer Drawing:

