

Research use only. Not for use in diagnostic procedures.

## Technical datasheet: Dimensions of polypropylene Storplate-96, round well, round bottom (2.0 mL volume)

- Part numbers:** PN 6008490 Natural polypropylene, untreated
- Material and treatment:** Round well, polypropylene 96-well microplate with U-shaped bottoms. Well volume 2.0 mL. Untreated, non-sterile.
- Packaging:** Part number 6008490 is a case with 50 plates, packed as 10 bags of 5 plates each.

**Dimensions:**

