

Research use only. Not for use in diagnostic procedures.

# Technical datasheet: Dimensions of polypropylene microplate-384 U-bottom, 35µL volume

**Part numbers:** PN 6008890 Natural polypropylene, untreated

**Material and treatment:** Low volume (35 µL), U-bottom, untreated, clear polypropylene microplate.

**Packaging:** Part number 6008890 is a case with 50 plates, packed as 5 bags of 10 plates each.

**Dimensions -- Customer Drawing:**

