

**PCR-grade Water**

Nuclease-free Water

Cat. No.	Amount
PCR-258S	10 x 1,2 ml
PCR-258L	50 ml
PCR-258-100	100 ml
PCR-258-500	500 ml

For in vitro use only!**Shipping:** shipped at ambient temperature**Storage Conditions:** store at 4 °C or -20 °C**Shelf Life:** 12 months**Form:** liquid**Description:**

PCR-grade Water is intended for use in molecular biology applications including PCR and RT-PCR. The ultra-pure and sterile filtered water is manufactured free of detectable inhibitors, contaminants or enzymatic activity.

Quality Control:

Each lot of PCR-grade Water is functionally tested in highly sensitive Low Copy Long Range PCR (18 kb fragment from lambda DNA, template dilution series). The absence of DNA contaminations (bacterial and human DNA) and Nucleases (DNases and RNases) is certified.