

**Dengue Virus NS1 c-end Type 2**recombinant, *E. coli*

Cat. No.	Amount
PR-1289	100 µg

**Description:**

The *E. coli* derived recombinant protein contains the NS1 Dengue Virus c-end Type 2 immunodominant regions.

The dengue protein is purified by proprietary chromatographic technique.

**For in vitro use only!****Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** avoid freeze/thaw cycles**Shelf Life:** 12 months**Purity:** > 90 % (SDS-PAGE)**Form:** liquid (Supplied in 20 mM Tris-HCl pH 8.0, 8 M urea and 60 mM NaCl)