

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

Cat. No.	Amount
PR-1290	100 µg

**For general laboratory use.****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 100 mM Tris-HCl pH 8.0, 60 mM NaCl and 5 % Glycerol)**Description:**

The *E. coli* derived recombinant protein contains the NS1 Dengue Virus n-end Type 2 immunodominant regions. The Recombinant Dengue Virus NS1n is fused to a GST tag.

The dengue protein is purified by proprietary chromatographic technique.