DATA SHEET





Dengue Virus NS1 Type 2

recombinant, E. coli

Cat. No.	Amount
PR-1288	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: > 85 % (SDS-PAGE, 10%, commassie staining) **Form:** liquid (supplied in PBS and 25 mM K_2CO_3)

Description:

The *E. coli* derived recombinant protein contains the NS1 Dengue Virus full length Type 2 immunodominant regions, having an MW of 45 kDa.

The dengue protein is fused to 6 x His tag at C-terminus and purified by proprietary chromatographic technique.