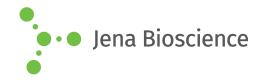
DATA SHEET





■ West Nile Virus Pre-M^{His}

WNV Pre-M recombinant, E. coli

Cat. No.	Amount
PR-1274	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:** > 95 % (SDS-PAGE)

Form: liquid (Supplied in 20 mM phosphate buffer pH 7.5)

Applications:

Antigen in ELISA and Western Blots, excellent antigen for detection of West-Nile virus with minimal specificity problems.

Description:

The $\it E.~coli$ derived recombinant protein contains the West-Nile N-Terminal Pre-M Virus immunodominant regions. The protein is fused with $\it 6xHistag.$

West Nile virus (WNV) is a virus of the family Flaviviridae part of the Japanese encephalitis (JE) antigenic complex of viruses. Image reconstructions and cryoelectron microscopyreveal a 45-50 nm virion covered with a relatively smooth proteinsurface. This structure is similar to virus, both belong to the genus flavivirus within the family Flaviviridae. WNV is a positive-sense, single strand of RNA, it is between 11,000 and 12,000 nucleotides long which encode seven non-structural proteins and three structural proteins. The RNA strand is held within a nucleocapsid formed from 12 kDa protein blocks, the capsid is contained within a host-derived membrane altered by two viral glycoproteins.

The protein is purified by proprietary chromatographic technique.

Selected References:

Wengler (1989) Cell-associated West Nile flavivirus is covered with E+pre-M protein heterodimers which are destroyed and reorganized by proteolytic cleavage during virus release. *J Virol.* **63**:2521.