

**VZV ORF26 (residues 9-33/184-208)**

Varicella-zoster Virus Protein
recombinant, *E. coli*

Cat. No.	Amount
PR-1255	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, RP-HPLC)

Form: liquid (Supplied in 25 mM Tris-HCl pH 8.0, 1 mM EDTA and 50% glycerol)

Applications:

Antigen in ELISA and Western blots, excellent antigen for detection of VZV with minimal specificity problems.

Description:

The protein contains the VZV ORF26 immunodominant regions, amino acids 9-33 and 184-208. The protein is purified by proprietary chromatographic technique.

Background: Varicella-zoster virus (VZV) is an extremely cell-associated alpha herpesvirus. It interacts with cell surface heparan sulfate proteoglycans during virus attachment.

Specificity: Immunoreactive with sera of VZV-infected individuals.

Selected References:

Pasieka *et al.* (2004) Regulation of varicella-zoster virus-induced cell-to-cell fusion by the endocytosis-competent glycoproteins gH and gE. *J. Virol.* **78**:2884.