

VZV ORF26 (residues 9-33, 184-208)

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

Cat. No.	Amount
PR-1255	100 µg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied as 1 mg/ml solution containing 25 mM Tris-HCl, pH 8.0, 1 mM EDTA, and 50% glycerol.

Specificity

Immunoreactive with sera of VZV-infected individuals.

Purity

>95% by SDS-PAGE, and RP-HPLC.

Description

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

Application

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

Background

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

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.