

## VZV ORF9 (residues 6-28, 76-100)

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

Cat. No.	Amount
PR-1254	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 ORF9 immunodominant regions, amino acids 6-28 and 76-100. 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. VZV tegument components include the regulatory proteins IE4, IE62, IE63 and the ORF10 protein, a protein kinase (ORF47) and an abundant protein encoded in ORF9 which is the homolog of HSV VP22. The kinase is able to phosphorylate IE62 and the ORF9 protein specifically in viral particles.

### Selected References:

Spengler *et al.* (2001) Interactions among structural proteins of varicella zoster virus. *Arch. Virol. Suppl.* **17**:71.