

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

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

Cat. No.	Amount
PR-1255	100 $\mu$ g

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

### **Avoid freeze / thaw cycles**

### **Form**

Liquid. Supplied in 25 mM Tris-HCl pH 8.0, 1 mM EDTA and 50% glycerol.

### **Application**

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

### **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.

### **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.