

**VZV-gE (aa 48-135)**

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

Cat. No.	Amount
PR-1253	100 µg

For in vitro use only!

Shipping: shipped on blue ice

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 60 mM NaCl, 50 mM Tris-HCl pH 8.0, 0.25% Sarkosil, 10 mM Glutathione 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-gE immunodominant regions, amino acids 48-135. 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. VZV-gE is a glycoprotein that plays an active or supportive role in VZV cell membrane fusion. VZV-gE was found to enhance the fusogenic potential of VZV gB.

Specificity: Immunoreactive with sera of VZV-infected individuals.

Selected References:

Maresova *et al.* (2005) Incorporation of three endocytosed varicellazoster virus glycoproteins, gE, gH, and gB, into the virion envelope. *J. Virol.* **79**:997.

Taha *et al.* (2004) Are false negative direct immunofluorescence assays caused by varicella zoster virus gE mutant strains? *J. Med. Virol.* **73**:631.

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.

Mo *et al.* (2002) The requirement of varicella zoster virus glycoprotein E (gE) for viral replication and effects of glycoprotein I on gE in melanoma cells. *Virology.* **304**:176.

Kenyon *et al.* (2002) Phosphorylation by the varicella-zoster virus ORF47 protein serine kinase determines whether endocytosed viral gE traffics to the trans-Golgi network or recycles to the cell membrane. *J. Virol.* **76**:10980.

Jaquet *et al.* (2002) Immunogenicity of a recombinant varicellazoster virus gE-IE63 fusion protein, a putative vaccine candidate against primary infection and zoster reactivation. *Vaccine.* **20**:1593.