





VZV ORF9 (residues 6-28/76-100)

Varicella-zoster Virus Protein recombinant, E. coli

Cat. No.	Amount	Applications:
PR-1254	100 µg	Antigen in ELIS VZV with mini

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)

ISA and Western blots, excellent antigen for detection of imal specificity problems.

Description:

The protein contains the VZV ORF9 immunodominant regions, amino acids 6-28 and 76-100. The protein is purified by proprietary chromatographic technique.

Background: Varicella-zoster virus (VZV) is an extremely cellassociated 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 ORFIO 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.

Specificity: Immunoreactive with sera of VZV-infected individuals.

Selected References:

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

