

CMV pp38 (residues 117-373) Cytomegalo Virus Phosphoprotein 38 recombinant, *E. coli*

Cat. No.	Amount
PR-1249	100 μ 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 CMV with minimal specificity problems.

Specificity

Immunoreactive with sera of CMV-infected individuals.

Molecular Weight

52.8kDa

Purity

>95% by SDS-PAGE and RP-HPLC

Description

The protein contains the CMV pp38 immunodominant regions, amino acids 117-373.

The protein is purified by proprietary chromatographic technique.

Background

Human cytomegalovirus (HCMV), a member of the herpesvirus family, demonstrates cell specificity for virus assembly and release.

In the viral particle this antigen is bound via S-S bridge to a lower Mr compound to form a final complex of 62 kDa. p38 is present both in the nucleus and in the cytoplasm of CMV-infected cells starting from 24 h postinfection and increasingly thereafter. Immunoelectron microscopy revealed that this antigen is mainly associated with the internal portion of viral capsids both in the nucleus and in the cytoplasm.

Selected References:

Landinin *et al.* (1991) Human cytomegalovirus structural components: intracellular and intraviral localization of p38. *Virus. Res.* **19**:189.