

## CMV pp150 (residues 1011-1048) Cytomegalo Virus Phosphoprotein 150 recombinant, *E. coli*

Cat. No.	Amount
PR-1252	100 µg

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 7.2, 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.

### Purity

>95% by SDS-PAGE

### Description

The protein contains the CMV pp150 immunodominant regions, amino acids 1011-1048.

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.

The human cytomegalovirus (HCMV) basic phosphoprotein pp150, encoded by the UL32 gene, together with the two other major phosphoproteins, pp65 (ppUL83) and pp71 (ppUL82) and several minor structural proteins, form the tegument around the viral nucleocapsid.

Western blot analysis of isolated cell fractions showed that pp150 is initially (48 h post-infection) localized in the nucleus, associated either with the nuclear membrane or with viral assembly regions, and later (72 h post-infection) in the cytoplasm.

### Selected References:

- Gentile *et al.* (2000) Detection of cytomegalovirus late (pp150) antigen in peripheral blood leukocytes of patients with human immunodeficiency virus-related lymphoma. *J. Infect. Dis.* **182**:376.
- Gyulai *et al.* (2000) Cytotoxic T lymphocyte (CTL) responses to human cytomegalovirus pp65, IE1-Exon4, gB, pp150, and pp28 in healthy individuals: reevaluation of prevalence of IE1-specific CTLs. *J. Infect. Dis.* **181**:1537.
- Boivin *et al.* (1999) Expression of the late cytomegalovirus (CMV) pp150 transcript in leukocytes of AIDS patients is associated with a high viral DNA load in leukocytes and presence of CMV DNA in plasma. *J. Infect. Dis.* **179**:1101.
- Li *et al.* (1997) Human cytomegalovirus matrix protein PP150 is efficiently presented as one of target antigens for cytotoxic T lymphocyte recognition. *Chin. Med. J. (Engl)*. **110**:397.
- Hendrix *et al.* (1996) Detection and sequence analysis of the major immediate early and PP150 gene of latent human cytomegalovirus in spleen, liver, and kidney tissues of trauma victims. *J. Med. Virol.* **50**:193.
- Hensel *et al.* (1995) Nuclear localization of the human cytomegalovirus tegument protein pp150 (ppUL32). *J. Gen. Virol.* **76**:1591.