

CMV pp52 (residues 202-434) Cytomegalo Virus Phosphoprotein 52 recombinant, *E. coli*

Cat. No.	Amount
PR-1250	100 µg

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 50mM Tris-HCl pH 7.2, 60mM NaCl, 10mM glutathione 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

51 kDa +26 kDa GST

Purity

>95% by SDS-PAGE

Description

The protein contains the CMV pp52 immunodominant regions, amino acids 202-434. It is fused to a 26 kDa GST-tag.

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) DNA polymerase is composed of a catalytic subunit, Pol, or UL54, which possesses a basal DNA polymerase activity, and an accessory protein, UL44 (pp52). UL44 can form homodimers in both its crystal structure and in solution.

Selected References:

- Nevels *et al.* (2004) Human cytomegalovirus immediate-early 1 protein facilitates viral replication by antagonizing histone deacetylation. *Proc. Natl. Acad. Sci. U.S.A.* **101**:17234.
- Lerner *et al.* (2004) Prevalence of abnormal cardiac wall motion in the cardiomyopathy associated with incomplete multiplication of Epstein-barr Virus and/or cytomegalovirus in patients with chronic fatigue syndrome. *In Vivo.* **18**:417.
- Xu *et al.* (2004) Human cytomegalovirus UL84 insertion mutant defective for viral DNA synthesis and growth. *J. Virol.* **78**:10360.
- Loregian *et al.* (2004) Specific residues in the connector loop of the human cytomegalovirus DNA polymerase accessory protein UL44 are crucial for interaction with the UL54 catalytic subunit. *J. Virol.* **78**:9084.
- Lee *et al.* (2004) Sumoylation of the major immediateearly IE2 protein of human cytomegalovirus Towne strain is not required for virus growth in cultured human fibroblasts. *J. Gen. Virol.* **85**:2149.
- Appelton *et al.* (2004) The cytomegalovirus DNA polymerase subunit UL44 forms a C clamp-shaped dimer. *Mol. Cell.* **15**:233.