

CMV pp52 (residues 202-434)

Cytomegalo Virus Phosphoprotein 52
recombinant, *E. coli*

Cat. No.	Amount
PR-1250-1	1 mg

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied as 1 mg/ml solution containing 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

51 kDa

Purity

>95% by SDS-PAGE

Description

The protein contains the CMV pp52 immunodominant regions, amino acids 202-434.

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.

CMV pp52 (residues 202-434)

Cytomegalo Virus Phosphoprotein 52

recombinant, *E. coli*

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.