





## **CMV pp28 (residues 130-160)**

Cytomegalo Virus Phosphoprotein 28 recombinant, *E. coli* 

Cat. No.	Amount
PR-1248	100 µg

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 50 mM Tris-HCl pH 7.2, 1 mM EDTA and 50% glycerol)

## **Applications:**

Antigen in ELISA and Western blots, excellent antigen for detection of CMV with minimal specificity problems.

## **Description:**

The protein contains the CMV pp28 immunodominant regions, amino acids 130-160. 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 pp28 is a myristylated phosphoprotein that is a constituent of the virion. The pp28 protein is positioned within the tegument of the virus particle, a protein structure that resides between the capsid and envelope. In the infected cell, pp28 is found in a cytoplasmic compartment derived from the Golgi apparatus, where the virus buds into vesicles to acquire its final membrane.

**Specificity:** Immunoreactive with sera of CMV-infected individuals.

## Selected References:

Jones *et al.* (2004) An acidic cluster of human cytomegalovirus UL99 tegument protein is required for trafficking and function. *J. Virol.* **78**:1488.

Britt *et al.* (2004) Rapid genetic engineering of human cytomegalovirus by using a lambda phage linear recombination system: demonstration that pp28 (UL99) is essential for production of infectious virus. *J. Virol.* **78**:539.

Silva *et al.* (2003) Human cytomegalovirus UL99-encoded pp28 is required for the cytoplasmic envelopment of tegument-associated capsids. *J. Virol.* **77**:10594.

Wu *et al.* (2001) Late temporal gene expression from the human cytomegalovirus pp28US (UL99) promoter when integrated into the host cell chromosome. *J. Gen. Virol.* **82**:1147.

Sanchez *et al.* (2000) Human cytomegalovirus pp28 (UL99) localizes to a cytoplasmic compartment which overlaps the endoplasmic reticulum-golgi-intermediate compartment. *J. Virol.* **74**:3842.

Kerry *et al.* (1997) Translational regulation of the human cytomegalovirus pp28 (UL99) late gene. *J. Virol.* **71**:981.

