

**EBV p18 (residues 1-119)**

Epstein-Barr Virus Capsid Antigen  
recombinant, *E. coli*

Cat. No.	Amount
PR-1225	100 µg

**For in vitro use only!**

**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)

**Form:** liquid (Supplied in 50 mM Tris-HCl pH 8.0, 60 mM NaCl and 50% glycerol)

**Applications:**

Antigen in ELISA and Western blots, excellent antigen for detection of HHV-4 (EBV) with minimal specificity problems.

**Description:**

Recombinant Epstein-Barr Virus protein contains the EBV (HHV-4) p18 fragment, amino acids 1-119. The protein is purified by proprietary chromatographic technique.

**Background:** Epstein-Barr virus, frequently referred to as EBV, is a member of the gamma herpesvirus family and one of the most common human viruses. The virus occurs worldwide, and most people become infected with EBV sometime during their lives. It persists in B lymphocytes for the life of the host. The small capsid protein p18 is highly immunogenic in humans, and the essential B-cell epitopes have been mapped to the carboxy region. It is described as a late antigen. IgG to the viral capsid antigen appears in the acute phase, peaks at 2 to 4 weeks after onset, declines slightly, and then persists for life.

**Specificity:** Immunoreactive with all sera of EBV infected individuals.

**Selected References:**

Feng *et al.* (2005) Serological diagnosis of infectious mononucleosis by chemiluminescent immunoassay using capsid antigen p18 of Epstein-Barr virus. *Clin. Chim. Acta.* **354**:77.

Faerber *et al.* (2001) Serological diagnosis of Epstein-Barr virus infection by novel ELISAs based on recombinant capsid antigens p23 and p18. *J. Med. Virol.* **63**:271.

Hinderer *et al.* (1999) Serodiagnosis of Epstein-Barr virus infection by using recombinant viral capsid antigen fragments and autologous gene fusion. *J. Clin. Microbiol.* **37**:3239.

Van Grunsven *et al.* (1993) Gene mapping and expression of two immunodominant Epstein-Barr virus capsid proteins. *J. Virol.* **67**:3908.