



HTLV-1 Envelope

Human T cell Leukemia Virus Type-1, Envelope Protein recombinant, *E. coli*

Cat. No.	Amount
PR-1219	100 µg

For *in vitro* use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: 27 kDa

Purity: > 95 % (SDS-PAGE, RP-HPLC)

Form: liquid (Supplied in 10 mM NaPO₄ pH 6.0, 0.1% SDS and 1 mM DTT)

Applications:

Antigen in ELISA and Western Blots. Excellent reagent for correct detection of HTLV infections with minimal specificity problems.

Description:

The protein contains the C- terminus of gp46 and most of p21E of HTLV-1. This non-fusion *E. coli* derived protein starts from HTLV-1 envelope amino acids 165 and ends with 440. The protein is purified by proprietary chromatographic technique.

Background: Human T-cell leukemia virus (HTLV)-1 and -2 are deltaretroviruses that infect a wide range of cells. HTLV-1 has been found primarily in CD4⁺ and CD8⁺ T-lymphocytes *in vivo*, whereas CD8⁺ T-lymphocytes are thought to be the *in vivo* reservoir of HTLV-2.

Specificity: Immunoreactive with all sera of HTLV-1 and HTLV-2 infected individuals with antibody response to HTLV envelope.

Selected References:

Hernandez *et al.* (2001) Chimeric synthetic peptides containing two immunodominant epitopes from the envelope gp46 and the transmembrane gp21 glycoproteins of HTLV-I virus. *Biochem. Biophys. Res. Commun.* **289**:1.

Armand *et al.* (2000) Targeted expression of HTLV-I envelope proteins in muscle by DNA immunization of mice. *Vaccine.* **18**:2212.

Tallet *et al.* (2000) Expression, purification and biological properties of the carboxyl half part of the HTLV-I surface envelope glycoprotein. *J. Chromatogr. B. Biomed. Sci. Appl.* **737**:85.

Georges-Gobinet *et al.* (1998) HTLV-I associated sicca syndrome in Guadeloupe: lack of relation with a peculiar encoding sequence of surface envelope glycoprotein. *Virus. Genes.* **16**:195.

Grange *et al.* (1998) Identification of exposed epitopes on the envelope glycoproteins of human T-cell lymphotropic virus type I (HTLV-I). *Int. J. Cancer.* **75**:804.

Grange *et al.* (1997) Induction of neutralizing antibodies against HTLV-I envelope proteins after combined genetic and protein immunizations in mice. *DNA Cell. Biol.* **16**:1439.

Kim *et al.* (2004) HTLV-1 and -2 envelope SU subdomains and critical determinants in receptor binding. *Retrovirology.* **1**:41.