

## HTLV-1 Mosaic (residues 374-400, 190-207)

Human T-cell leukemia virus 1, Envelope mosaic protein recombinant, *E. coli*

Cat. No.	Amount
PR-1221-1	1 mg

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

### Avoid freeze / thaw cycles

### Form

Liquid. Supplied in 50 mM NaPO<sub>4</sub> pH 6.0, 1 mM DTT and 1 mM EDTA.

### Application

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

### Specificity

Immunoreactive with all sera of HTLV-1 infected individuals.

### Purity

>95% by SDS-PAGE and RP-HPLC

### Description

The protein contains the HTLV-1 gp21/gp46 immunodominant fragments, amino acids: 374-400 (gp21) and 190-207 (gp46).

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<sup>+</sup> and CD8<sup>+</sup> T-lymphocytes *in vivo*, whereas CD8<sup>+</sup> T-lymphocytes are thought to be the *in vivo* reservoir of HTLV-2.

### Selected References:

- Toedter *et al.* (1991) Development of a monoclonal antibody-based p24 capsid antigen detection assay for HTLV-I, HTLV-II, and STLV-I infection. *AIDS. Res. Hum. Retroviruses*. **8**:527.
- Segal-Eiras *et al.* (1991) HTLV-I p24 antigen in circulating immune complexes associated with acute Tlymphoblastic leukemia. *Haematologica*. **76**:441.
- Stransky *et al.* (1991) Ultrastructural localization of HTLV-I gag proteins p19 and p24 by single and double immunogold labeling. *J. Histochem. Cytochem.* **39**:185.
- Ando *et al.* (1990) p24 protein of a human T-lymphotropic virus type-I (HTLV-I) producing cell line (TCL-Kan) reacts with some normal pregnant women's sera. *Nippon. Sanka. Fujinka. Gakkai. Zasshi*. **42**:1251.
- Ziegler *et al.* (1989) Immunohistochemical localization of HTLV-I p19- and p24-related antigens in synovial joints of patients with rheumatoid arthritis. *Am. J. Pathol.* **135**:1.
- Grimaldi *et al.* (1988) HTLV-I-associated myelopathy: oligoclonal immunoglobulin G bands contain anti-HTLV-I p24 antibody. *Ann. Neurol.* **24**:727.