



HTLV-1 Mosaic (residues 374-400/190-207)

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

| Cat. No. | Amount |
|----------|--------|
| PR-1221 | 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 20 mM Phosphate buffer pH 7.5)

Applications:

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

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

Specificity: Immuno reactive with all sera of HIV-1 infected individuals.

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.