



## HTLV-1 p24 Core



Human T-cell Leukemia Virus Type-1, core protein p24 recombinant, *E. coli* 

Cat. No.	Amount	<b>Applications:</b> Antigen in ELI: detection of HIV
PR-1220	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 NaPO4 pH 6.0, 1 mM DTT and 1 mM EDTA)

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

## **Description:**

The protein contains the full-length HTLV-1 major core protein p24. 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.

**Specificity:** Immuno reactive with all sera of HIV-I 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.

