

HTLV-1+2 (fusion protein) Human T-cell leukemia virus 1+2 recombinant, *E. coli*

Cat. No.	Amount
PR-1222	25 µg

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied as 1 mg/ml solution containing 25 mM Tris-HCl, pH 8.0, 250 mM glycine and 0.1% SDS.

Applications

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 and HTLV-2 infected individuals.

Purity

>95% by SDS-PAGE and RP-HPLC.

Description

Recombinant Human T Lymphotropic Virus fused envelope proteins (HTLV-1/2), having a molecular weight of 56 kDa.

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.

Selected References:

- Lee *et al.* (1994) A multicentric seroepidemiological survey of HTLV-I/II in Italy. *Clin. Diagn. Virol.* **2**:139.
Bocalandro *et al.* (2004) Antigenic mixture of synthetic peptides for the immunodiagnosis of HTLV I/II infection. *J. Immunoassay. Immunochem.* **25**:205.
Tamim *et al.* (2004) Is seroprevalence of HTLV-I/II among blood donors in Lebanon relevant? *Am. J. Infect. Control.* **32**:220.
Gastaldello *et al.* (2004) Seroepidemiology of HTLV-I/II in Argentina: an overview. *J. Acquir. Immune. Defic. Syndr.* **35**:301.