

HIV-2 Envelope

Human Immunodeficiency Virus 2 Antigen recombinant, *E. coli*

Cat. No.	Amount
PR-1218	100 μ g

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 6 M urea.

Endotoxin

Less than 0.1 ng/ μ g (IEU/ μ g) of HIV-2 Envelope.

Purity

>97% by SDS-PAGE and RP-HPLC

Description

This antigen is an *E. coli*-derived recombinant protein that composes all of the reported immunogenic determinants found in gp36. The gene encoding this fusion protein was synthesized using codons optimized for *E. coli* expression and doesn't represent a linear HIV-2 envelope sequence. rHIV-2 is a non-glycosylated, 135 amino acid polypeptide chain, having a molecular mass of 16.1 kDa.

Recombinant HIV-2 Envelope is purified by proprietary chromatographic technique.

Background

HIV belongs to the retrovirus family, distinguished by possession of a viral reverse transcriptase that transcribes viral RNA into DNA which is integrated into the host-cell genome.

gp36 is the transmembrane envelope protein from HIV-2 comparable to gp41 from HIV-1.

Selected References:

Chen *et al.* (1995) HIV-2 transmembrane protein gp36 like HIV-1 gp41 binds to human lymphocytes and monocytes. *AIDS*. **9**:1193.