

**HIV-2 Envelope**

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

Cat. No.	Amount
PR-1218	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: > 97 % (SDS-PAGE, RP-HPLC)

Form: liquid (Supplied in 50 mM Tris-HCl pH 7.2 and 6 M urea)

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.

Endotoxin: Less than 0.1 ng/µg (IEU/µg) of HIV-2 Envelope.

Selected References:

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