

**HIV-1 O env mosaic 9 kDa**

Human immunodeficiency virus type 1 group O envelope protein, mosaic recombinant, *E. coli*

Cat. No.	Amount
PR-1437-S	200 µg
PR-1437-L	1 mg

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at 4 °C

Shelf Life: 6 months

Molecular Weight: about 9 kDa

Purity: ≥ 85 %

Form: liquid (Supplied in Sodium Phosphate 100 mM pH 8.0, Tris-HCl 10 mM, Imidazole 300 mM, Urea 8 M)

Applications:

Suitable for use in ELISA and LIA. Each laboratory should determine an optimum working titer for use in its particular application. Other applications have not been tested but use in such assays should not necessarily be excluded.

Description:

The protein contains immunodominant regions of HIV-1 group O gp160, amino acids (297-327)(581-610). The protein is fused to a 6 His TAG.

Specificity:

Immunoreactive with sera of HIV-infected individuals