

**HIV-2 gp36 (67-98) 36 kDa**

Human immunodeficiency virus type 2 gp36 protein recombinant, *E. coli*

Cat. No.	Amount
PR-1433-S	200 µg
PR-1433-L	1 mg

For general laboratory use.

Shipping: shipped at 4 °C

Storage Conditions: store at 4 °C

Shelf Life: 6 months

Molecular Weight: about 36 kDa

Purity: ≥ 85 %

Form: liquid (Supplied in PBS pH 7.4, Sarkosyl 0.2 %, ProClin 300 0.1 %)

Applications:

Suitable for use in ELISA. 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 region of HIV-2 gp36, amino acids 67-98. The protein is fused to a GST, 6 His, S-tag.

Specificity:

Immunoreactive with sera of HIV 2-infected individuals