

**HIV-1 M gp160 50 Kda**

Human immunodeficiency virus type 1 group M gp160 protein recombinant, *E. coli*

Cat. No.	Amount
PR-1434-S	200 µg
PR-1434-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 50 kDa

Purity: ≥ 85 %

Form: liquid (Supplied in Sodium Phosphate 100 mM pH 4.5, Tris-HCl 10 mM, Imidazole 300 mM, Urea 8 M)

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-1 group M gp160, amino acids 283-674. The protein is fused to a 6 His TAG.

Specificity:

Immunoreactive with sera of HIV-infected individuals