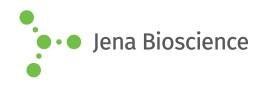
DATA SHEET





📈 🔳 HIV-1 M gp120 (v3 loop) 36 kDa

Human immunodeficiency virus type 1 group M gp120 (v3 loop regions) protein recombinant, *E. coli*

Cat. No.	Amount
PR-1436-S	200 μg
PR-1436-L	1 mg

For general laboratory use.

Shipping: shipped on gel packs **Storage Conditions:** store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 6 months

Molecular Weight: about 36 kDa

Purity: 60 - 85 %

Form: liquid (Supplied in Tris-HCl 50 mM pH 8.0, Sodium Chloride 60

mM, Glutathione 10 mM, Sarcosyl 0.25 %, Glycerol 50 %)

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 the V3 loop region of HIV-1 group M gp 120 subtype C, amino acids 301-330. The protein is fused to a GST, 6 His, S-tag.

Specificity:

Immunoreactive with sera of HIV-infected individuals