

**HIV-1 M gp41 (537-811) 58 kDa**

Human immunodeficiency virus type 1 group M gp41 protein recombinant, *E. coli*

Cat. No.	Amount
PR-1429-S	200 µg
PR-1429-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 58 kDa

Form: liquid (Supplied in Imidazole 20 mM pH 7.0, Sodium Chloride 150 mM, Urea 8M, DTT 10 mM)

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 gp41, amino acids 537-811. The protein is fused to a β-galactosidas + 6 HIS TAG.

Specificity:

Immunoreactive with sera of HIV-infected individuals