# **DATA SHEET**





## HIV-2 gp36 (531-851) 62 kDa

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

Cat. No.	Amount
PR-1431-S	200 μg
PR-1431-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 62 kDa

**Purity:** ≥ 85 %

Form: liquid (Supplied in Tris-HCl 20 mM pH 8.0, DTT 1 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-2 gp 36, amino acids 531-851. The protein is fused to a ß-galactosidas + 6 HIS TAG.

#### Specificity:

Immunoreactive with sera of HIV-infected individuals