

**HSV-1 gG 36 kDa**

Herpes simplex virus type 1 gG protein
recombinant, *E. coli*

Cat. No.	Amount
PR-1441-S	200 µg
PR-1441-L	1 mg

For in vitro use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: about 36 kDa

Purity: ≥ 85 %

Form: liquid (Supplied in Tris-HCl 25 mM pH 8.0, Glutathione 5 mM, Urea 1.5 M, 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 immunodominant region of HSV 1 glycoprotein G, amino acids 84-175. The protein is fused to a GST TAG.

Specificity:

Immunoreactive with sera of HSV 1-infected individuals