

HSV-1 gG (residues 84-175)

Herpes Simplex Virus-1 glycoprotein G
recombinant, *E. coli*

Cat. No.	Amount
PR-1109-1	1 mg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 25 mM Tris-HCl pH 8.0, 5 mM glutathione, 1.5 M urea and 50% glycerol.

Protein synonyms/aliases

Glycoprotein G.

Application

Recombinant rHSV-1 gG Antigen may be used in ELISA and Western blots, excellent for detection of HSV with minimal specificity problems.

Specificity

Immunoreactive with sera of HSV-infected individuals.

MW

about 40 kDa

Tag

GST

Purity

>95% by SDS-PAGE

Description

Recombinant HSV-1 gG contains the immunodominant regions of glycoprotein G, amino acids: 84-175.

HSV-1-gG is purified by proprietary chromatographic techniques.

Background

Glycoprotein G-1 (gG-1), a viral envelope glycoprotein that was suggested to contribute to viral entry through apical surfaces of polarized cells, has been utilized as a prototype antigen for HSV-1 type-specific diagnosis due to lack of cross-reactivity with its counterpart in HSV-2, i.e., gG-2.

Herpes simplex virus type 1 (HSV-1) encodes 11 envelope glycoproteins, of which glycoprotein G-1 (gG-1) induces a type-specific antibody response.

HSV-1 gG (residues 84-175)

Herpes Simplex Virus-1 glycoprotein G
recombinant, *E. coli*

Selected References:

- Rekabdar *et al.* (2002) Dichotomy of glycoprotein g gene in herpes simplex virus type 1 isolates. *J. Clin. Microbiol.* **40**:3245.
- Tunback *et al.* (2000) Glycoprotein G of herpes simplex virus type 1: identification of type-specific epitopes by human antibodies. *J. Gen. Virol.* **81**:1033.
- Anderson *et al.* (2000) Pseudotyping of glycoprotein D-deficient herpes simplex virus type 1 with vesicular stomatitis virus glycoprotein G enables mutant virus attachment and entry. *J. Virol.* **74**:2481.
- Rekabdar *et al.* (1999) Variability of the glycoprotein G gene in clinical isolates of herpes simplex virus type 1. *Clin. Diagn. Lab. Immunol.* **6**:826.
- Sullivan *et al.* (1987) Expression and characterization of herpes simplex virus type 1 (HSV-1) glycoprotein G (gG) by recombinant vaccinia virus: neutralization of HSV-1 infectivity with anti-gG antibody. *J. Gen. Virol.* **68**:2587.