

---

Product Manual

# Herpes Simplex Virus Glycoprotein D (gD) ELISA Kit

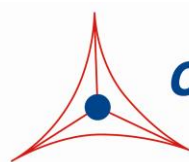
Catalog Numbers

VPK-5178

96 assays

**FOR RESEARCH USE ONLY**  
**Not for use in diagnostic procedures**

---



**CELL BIOLABS, INC.**  
*Creating Solutions for Life Science Research*

## **Introduction**

Herpes simplex virus (HSV) is a relatively large enveloped virus with a DNA linear double-stranded genome. There are two main HSV types, HSV-1 and HSV-2. HSV produce a wide variety of illnesses, including mucocutaneous infections, central nervous system infections, and occasionally, infections of the visceral organs.

Four viral glycoproteins (gD, gB, gH, and gL) are required for the entry of HSV into cells. Glycoprotein D (gD) is the receptor binding protein, and its binding to the receptor is proposed to activate the heterodimer gH/gL to trigger gB to mediate the fusion of viral and cellular membranes. Receptors for HSV gD include nectin-1, herpes virus entry mediator 8, and a modified form of heparan sulfate.

Cell Biolabs' Herpes Simplex Virus Glycoprotein D (gD) ELISA Kit is an enzyme immunoassay developed for detection and quantitation of the HSV glycoprotein D. The ELISA antibodies recognize the glycoprotein D from both HSV-1 and HSV-2. The kit has a detection sensitivity limit of 313 pg/mL HSV gD. Each kit provides sufficient reagents to perform up to 96 assays including standard curve and HSV lysate samples.

## **Assay Principle**

An anti-HSV gD monoclonal antibody is adsorbed onto a microtiter plate. HSV gD present in the sample or standard binds to the antibody adsorbed on the plate; an FITC-conjugated anti-HSV gD monoclonal antibody is added and binds to the HSV gD captured by the first antibody. Following incubation and wash steps, an HRP-conjugated mouse anti-FITC antibody is added and binds to the FITC conjugated anti-HSV gD monoclonal antibody. Unbound HRP-conjugated mouse anti-FITC antibody is removed during a wash step, and Substrate Solution reactive with HRP is added to the wells. A colored product is formed in proportion to the amount of HSV gD present in the sample. The reaction is terminated by addition of Stop Solution and absorbance is measured at 450 nm. A standard curve is prepared from the provided recombinant HSV gD standard, and the sample HSV gD concentration is then determined.

## **Related Products**

1. VPK-108-H: QuickTiter™ Lentivirus Quantitation Kit (HIV-1 p24 ELISA)
2. VPK-150: QuickTiter™ Hepatitis B Core Antigen (HBcAg) ELISA Kit
3. VPK-151: QuickTiter™ Hepatitis C Core Antigen (HCcAg) ELISA Kit
4. VPK-156: QuickTiter™ MuLV Core Antigen (MuLV p30) ELISA Kit
5. VPK-5169: Adenovirus Hexon ELISA Kit
6. VPK-5170: RSV Fusion Protein ELISA Kit
7. VPK-5171: RSV Nucleoprotein ELISA Kit
8. VPK-5172: Human Cytomegalovirus Glycoprotein B ELISA Kit
9. VPK-5174: Influenza A Nucleoprotein ELISA Kit
10. VPK-5175: Influenza B Nucleoprotein ELISA Kit

## **Kit Components**

### **Box 1 (shipped at room temperature)**

1. Anti-HSV gD Antibody Coated Plate (Part No. 51781B): One strip well 96-well plate.
2. FITC-Conjugated Anti-HSV gD Monoclonal Antibody (Part No. 51782C): One 20  $\mu$ L vial.
3. HRP-Conjugated Anti-FITC Monoclonal Antibody (Part No. 310811): One 20  $\mu$ L vial.
4. Assay Diluent (Part No. 310804): One 50 mL bottle.
5. 10X Viral Lysis Buffer (Part No. 51693B): One 15 mL bottle containing 200 mM Tris, pH 7.5, 1500 mM NaCl, 10% Triton X-100, 1% SDS.
6. 10X Wash Buffer (Part No. 310806): One 100 mL bottle.
7. Substrate Solution (Part No. 310807): One 12 mL amber bottle.
8. Stop Solution (Part. No. 310808): One 12 mL bottle.

### **Box 2 (shipped on blue ice packs)**

1. Recombinant HSV-2 gD Standard (Part No. 51783D): One 100  $\mu$ L vial of 2  $\mu$ g/mL recombinant HSV-2 Glycoprotein D (Lys26-Tyr310) in PBS containing BSA.

## **Materials Not Supplied**

1. HSV Sample: purified virus or unpurified viral supernatant
2. Microcentrifuge
3. 10  $\mu$ L to 1000  $\mu$ L adjustable single channel micropipettes with disposable tips
4. 50  $\mu$ L to 300  $\mu$ L adjustable multichannel micropipette with disposable tips
5. Multichannel micropipette reservoir
6. Microplate reader capable of reading at 450 nm (620 nm as optional reference wave length)

## **Storage**

Upon receiving, aliquot and store the HSV-2 gD Standard at -20°C and avoid freeze/thaw. Store all other components at 4°C.

## **Safety Considerations**

Remember that your HSV samples contain infectious viruses before inactivation; you must follow the recommended NIH guidelines for all materials containing infectious organisms.

## **Preparation of Reagents**

- 1X Wash Buffer: Dilute the 10X Wash Buffer Concentrate to 1X with deionized water. Stir to homogeneity.

- FITC-Conjugated Anti-HSV gD Monoclonal Antibody and HRP-Conjugated Anti-FITC Monoclonal Antibody: Immediately before use dilute the FITC-conjugated antibody 1:1000 and HRP-conjugated antibody 1:1000 with Assay Diluent. Do not store diluted solutions.

### **Preparation of Standard Curve**

1. Prepare a dilution series of HSV-2 gD Standard in the concentration range of 20 ng/mL – 0.313 ng/mL by diluting the stock solution in Assay Diluent (Table 1).

<b>Standard Tubes</b>	<b>2 µg/mL HSV-2 gD Standard (µL)</b>	<b>Assay Diluent (µL)</b>	<b>HSV-2 gD (ng/mL)</b>
1	10	990	20
2	500 of Tube #1	500	10
3	500 of Tube #2	500	5
4	500 of Tube #3	500	2.5
5	500 of Tube #4	500	1.25
6	500 of Tube #5	500	0.625
7	500 of Tube #6	500	0.313
8	0	500	0

**Table 1. Preparation of HSV gD Standard**

2. Transfer 225 µL of each dilution to a microcentrifuge tube containing 25 µL of 10X Lysis Buffer. Perform the assay as described in Assay Protocol.

### **HSV Sample Inactivation and Lysis**

1. (Optional) Dilute HSV samples in culture medium. Include culture medium as a negative control.
2. Transfer 225 µL of each sample to a microcentrifuge tube containing 25 µL of 10X Lysis Buffer, vortex well. Inactivate HSV sample at 56°C for 30 min.
3. Centrifuge at 12,000 x g for 5 minutes at 4°C. Collect the supernatant as HSV lysate.

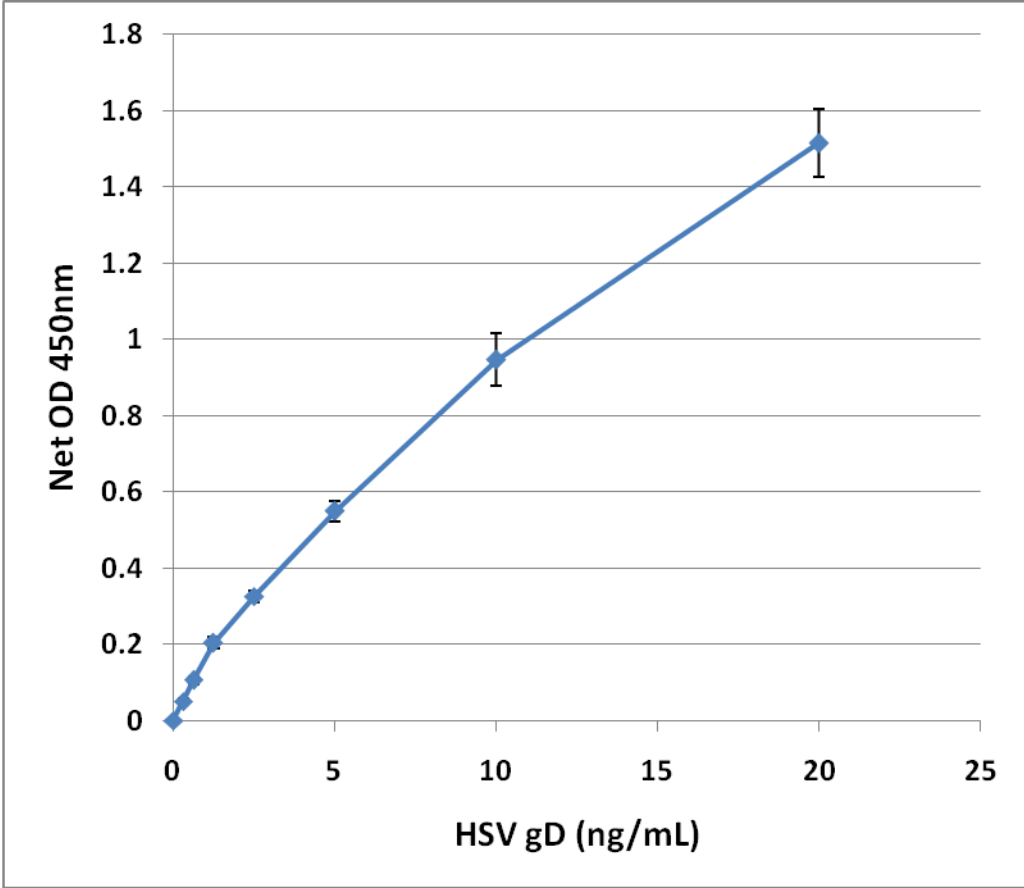
### **Assay Protocol**

1. Prepare and mix all reagents thoroughly before use.
2. Each HSV lysate sample, HSV gD standard, blank, and control medium should be assayed in duplicate.
3. Add 100 µL of HSV lysate or HSV gD standard to the Anti- HSV gD Antibody Coated Plate.
4. Cover with a plate cover and incubate at 37°C for 2 hours.

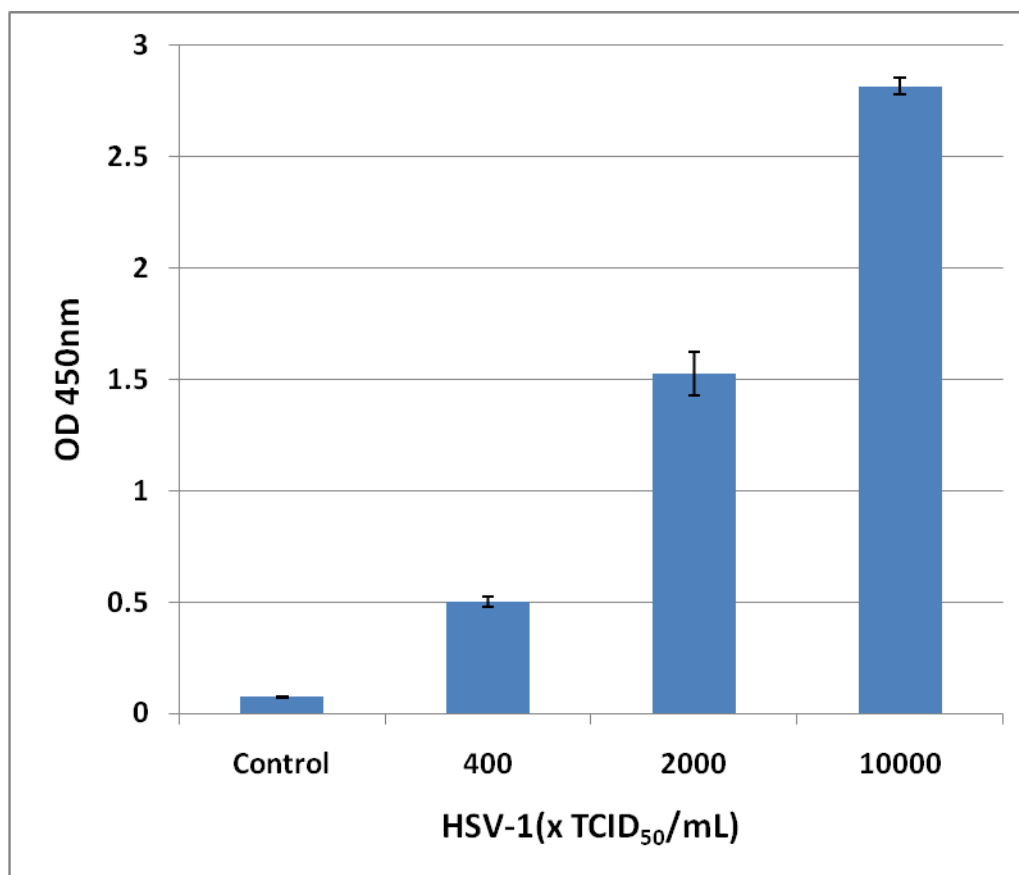
5. Remove plate cover and empty wells. Wash microwell strips 5 times with 250  $\mu$ L 1X Wash Buffer per well with thorough aspiration between each wash. After the last wash, empty wells and tap microwell strips on absorbent pad or paper towel to remove excess 1X Wash Buffer.
6. Add 100  $\mu$ L of the diluted FITC-Conjugated Anti-HSV gD Monoclonal Antibody to each well.
7. Cover with a plate cover and incubate at room temperature for 1 hour on an orbital shaker.
8. Remove plate cover and empty wells. Wash the strip wells 5 times according to step 5 above.
9. Add 100  $\mu$ L of the diluted HRP-Conjugated Anti-FITC Monoclonal Antibody to all wells.
10. Cover with a plate cover and incubate at room temperature for 1 hour on an orbital shaker.
11. Remove plate cover and empty wells. Wash microwell strips 5 times according to step 5 above. Proceed immediately to the next step.
12. Warm Substrate Solution to room temperature. Add 100  $\mu$ L of Substrate Solution to each well, including the blank wells. Incubate at room temperature on an orbital shaker. Actual incubation time may vary from 2-30 minutes.  
  
*Note: Watch plate carefully; if color changes rapidly, the reaction may need to be stopped sooner to prevent saturation.*
13. Stop the enzyme reaction by adding 100  $\mu$ L of Stop Solution into each well, including the blank wells. Results should be read immediately (color will fade over time).
14. Read absorbance of each microwell on a spectrophotometer using 450 nm as the primary wave length.

**Example of Results**

The following figures demonstrate typical Herpes Simplex Virus Glycoprotein D (gD) ELISA results. One should use the data below for reference only. This data should not be used to interpret actual results.



**Figure 1: HSV gD ELISA Standard Curve.**



**Figure 2: HSV gD in HSV-1 Culture Fluid.** HSV-1 culture fluid ( $1.15 \times 10^6$  TCID<sub>50</sub>/mL) was first diluted with culture medium, then heat inactivated and lysed in Viral Lysis Buffer. HSV-1 lysate was subjected to Herpes Simplex Virus Glycoprotein D (gD) ELISA Kit according to the Assay Protocol.

## References

1. Heldwein EE, Krummenacher C. (2008) Entry of herpesviruses into mammalian cells. *Cell Mol Life Sci.* **65**, 1653–1668.
2. Akhtar J, Shukla D. (2009) Viral entry mechanisms: cellular and viral mediators of herpes simplex virus entry. *FEBS J.* **276**, 7228–7236.
3. Connolly, S. A., Jackson, J. O., Jardetzky, T. S. & Longnecker, R. (2011) Fusing structure and function: a structural view of the herpesvirus entry machinery. *Nature reviews. Microbiology* **9**, 369–381.
4. Eisenberg, R. J. et al. (2012) Herpes virus fusion and entry: a story with many characters. *Viruses* **4**, 800–832.
5. Menotti, L. et al. (2000) The murine homolog of human Nectin1 $\delta$  serves as a species nonspecific mediator for entry of human and animal  $\alpha$ herpesviruses in a pathway independent of a detectable binding to gD. *Proc. Nat. Acad. Sci. USA* **97**, 4867–4872.
6. Montgomery, R. I., Warner, M. S., Lum, B. J. & Spear, P. G. (1996) Herpes simplex virus-1 entry into cells mediated by a novel member of the TNF/NGF receptor family. *Cell* **87**, 427–436.
7. Shukla, D. et al. (1999) A novel role for 3-O-sulfated heparan sulfate in herpes simplex virus 1 entry. *Cell* **99**, 13–22.

## **Warranty**

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS's sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

## **Contact Information**

Cell Biolabs, Inc.  
7758 Arjons Drive  
San Diego, CA 92126  
Worldwide: +1 858-271-6500  
USA Toll-Free: 1-888-CBL-0505  
E-mail: [tech@cellbiolabs.com](mailto:tech@cellbiolabs.com)  
[www.cellbiolabs.com](http://www.cellbiolabs.com)

©2023: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.