
Product Manual

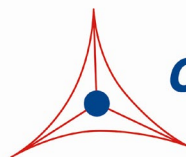
West Nile Virus Envelope Protein ELISA Kit

Catalog Number

VPK-5154

96 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

West Nile virus (WNV) is a single-stranded RNA virus that causes West Nile fever and is a member of the genus *Flavivirus*, which includes Zika virus, dengue virus, and yellow fever virus. WNV is transmitted mostly by mosquitoes, but the primary hosts of WNV are birds and the transmission cycle occurs from bird to mosquito to bird.

The WNV genome is a positive single stranded RNA encased by a nucleocapsid which is contained in a lipid bi-layered 50 nm envelope. A single polyprotein is expressed and cleaved by host and viral proteases into three structural (C, prM/M, and envelope, or E protein) and seven nonstructural (NS1, NS2A, NS2B, NS3, NS4A, NS4B, and NS5) proteins. The WNV E protein is responsible for cellular entry of WNV. The crystal structure of the WNV E protein shows that there are three domains: a β -barrel-shaped domain I, an elongated finger-like domain II, and a C-terminal immunoglobulin-like domain III⁸. Domain I is glycosylated at amino acid position 154, which is critical for WNV infection. The internal fusion peptide loop in domain II allows for the trimerization of the WNV E protein as well as initiation of virus entry into cells. Domain III controls binding of WNV to host cells.

Cell Biolabs' West Nile Virus Envelope Protein ELISA Kit is an enzyme immunoassay designed to measure WNV E Protein from cell or tissue samples. It provides sufficient reagents for up to 96 tests in a 96-well plate including standard curve and unknown samples. Detection sensitivity is 31.3 ng/mL.

Related Products

1. VPK-5145: SARS-Cov-2 Nucleocapsid ELISA Kit
2. VPK-107: Quick Titer™ Lentivirus Titer Kit (Lentivirus-associated p24 ELISA)
3. VPK-109: QuickTiter™ Adenovirus Titer Immunoassay Kit
4. VPK-112: QuickTiter™ Lentivirus Quantitation Kit
5. VPK-120: QuickTiter™ Retrovirus Quantitation Kit
6. VPK-145: QuickTiter™ AAV Quantitation Kit

Kit Components

Box 1 (shipped at room temperature)

1. Anti-WNV E Protein Antibody Coated Plate (Part No. 51541B): One 96-well strip plate (8 x 12).
2. Biotinylated Anti-WNV E Protein Antibody (1000X) (Part No. 51542C): One 10 μ L vial.
3. Streptavidin-Enzyme Conjugate (Part No. 310803): One 20 μ L vial.
4. Assay Diluent (Part No. 310804): One 50 mL bottle.

5. Lysis Solution (Part No. 51543A): One 15 mL bottle containing 0.5% Tween 20 and 1% Triton X-100 in 10X PBS. The lysis solution was developed according to Mayo and Beckwith, 2002 and Colavita et al, 2017 (See refs 7 and 8).
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. WNV E Protein Standard (Part No. 51544D): One 50 μ L vial of 100 μ g/mL WNV E Protein.

Materials Not Supplied

1. Recombinant Purified WNV Samples
2. Microcentrifuge
3. 10 μ L to 1000 μ L adjustable single channel micropipettes with disposable tips
4. 50 μ L to 300 μ 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 receipt, aliquot and store the WNV E Protein Standard at -80°C and the Biotinylated Anti-WNV E Protein Antibody at -20°C . Avoid multiple freeze/thaw cycles. Store all other components at 4°C .

Preparation of Reagents

- 1X Wash Buffer: Dilute the 10X Wash Buffer to 1X with deionized water. Stir to homogeneity.
- Biotinylated Anti-WNV E Protein Antibody and Streptavidin Enzyme Conjugate: Immediately before use dilute the Anti-WNV E Protein Antibody and the Streptavidin Enzyme Conjugate 1:1000 with Assay Diluent. Do not store diluted solutions.

Preparation of Standard Curve

1. Prepare a dilution series of WNV E Protein standards in the concentration range of 0 to 2000 ng/mL into Assay Diluent (Table 1).

Standard Tubes	WNV E Protein Standard (μL)	Assay Diluent (μL)	WNV E Protein (ng/mL)
1	10	490	2000
2	250 of Tube #1	250	1000
3	250 of Tube #2	250	500
4	250 of Tube #3	250	250
5	250 of Tube #4	250	125
6	250 of Tube #5	250	62.5
7	250 of Tube #6	250	31.3
8	0	250	0

Table 1. Preparation of WNV E Protein Standards

- Transfer 225 μL of each dilution (Standard Tubes 1-8) to a microcentrifuge tube containing 25 μL of Lysis Solution. Perform the assay as described in Assay Instructions.

Preparation and Inactivation of Samples

- (Optional) Dilute viral supernatant in culture medium as needed. For unknown samples we recommend several dilutions for each sample. Include culture medium as a negative control.
- Transfer 225 μL of each sample to a microcentrifuge tube containing 25 μL of Lysis Solution (see refs 7 and 8). Vortex well.
- Incubate 60 minutes at room temperature.

Assay Protocol

- Add 100 μL of WNV E Protein unknown sample or standard to the Anti-E Protein Antibody Coated Plate. Each WNV E Protein unknown sample, standard and blank should be assayed in duplicate.
- Incubate at room temperature for 1 hour on an orbital shaker.
- Wash microwell strips 3 times with 250 μL 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.
- Add 100 μL of the diluted Biotinylated Anti-WNV E Protein antibody to each well. Incubate at room temperature for 1 hour on an orbital shaker.
- Wash the strip wells 3 times according to step 3 above.
- Add 100 μL of the diluted Streptavidin-Enzyme Conjugate to each well. Incubate at room temperature for 1 hour on an orbital shaker.
- Wash the strip wells 3 times according to step 3 above. Proceed immediately to the next step.

- Warm Substrate Solution to room temperature. Add 100 μL 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.

- Stop the enzyme reaction by adding 100 μL of Stop Solution into each well, including the blank wells. Results should be read immediately (color will fade over time).
- Read absorbance of each microwell on a spectrophotometer using 450 nm as the primary wave length.

Example of Results

The following figures demonstrate typical results with the West Nile Virus Envelope Protein ELISA Kit. One should use the data below for reference only. This data should not be used to interpret actual results.

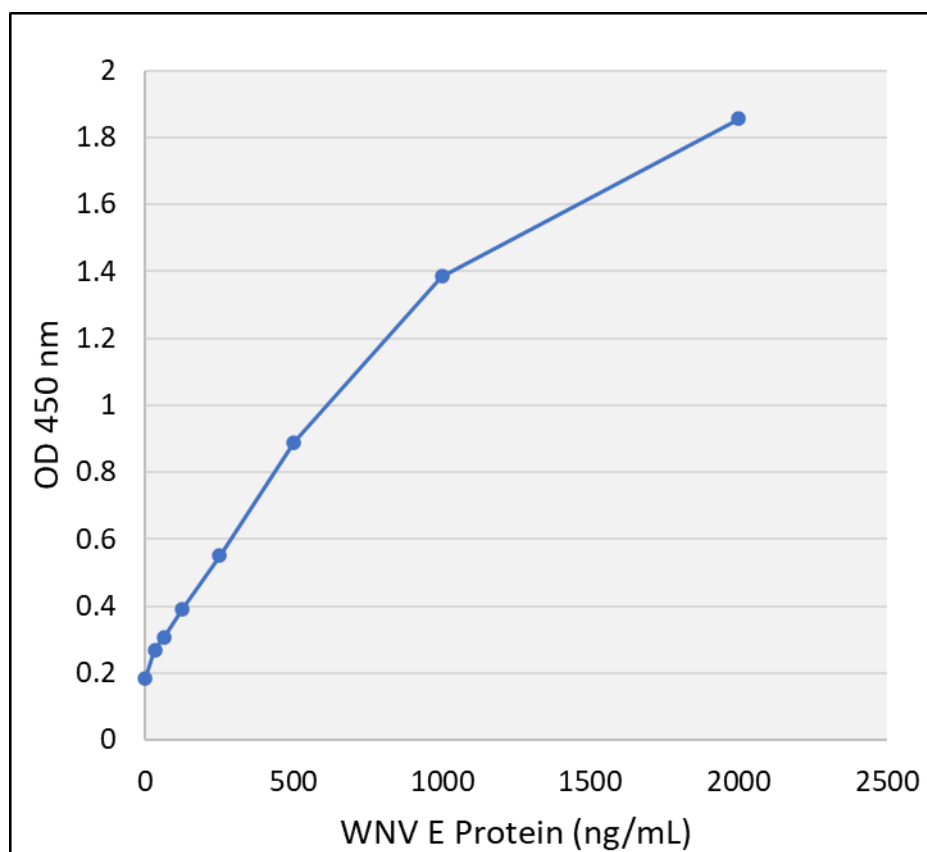


Figure 1: West Nile Virus Envelope Protein ELISA Kit Standard Curve.

References

1. Mackenzie JS, Gubler DJ, and Petersen LR (2004). *Nat. Med.* **10**: S98–S109.
2. Valiakos G, Athanasiou LV, Touloudi A, Papatsiros, Spyrou VV, Petrovska L and Billinis C (2013) *Intech*, **2**: 43-70.
3. Kobayashi S, Kaneko C, Kawakami R, Hasebe R, Sawa H, Yoshii K and Kariwa H (2020) *Sci. Reports*, **10**:7168-7177.
4. Oliphant T, Nybakken GE, Engle M, Xu Q, Nelson CA, Sukupolvi-Petty S, Marri A, Lachmi B-E, Olshevsky U, Fremont DH, Pierson TC, and Diamond MS (2006) *J Virol* **80**:12149-12159.
5. Nybakken GE, Nelson CA, Chen BR, Diamond MS, and Fremont DH (2006) *J Virol* **80**:11467-11474.
6. Maharaj PD, Langevin SA, Bolling BG, Andrade CC, Engle XA, Ramey WN, Bosco-Lauth A, Bowen RA, Sanders TA, Huang CY-H, Reisen WK, and Brault AC (2019) *PLoS Negl Trop Dis* **13**:e0007473
7. Mayo DR and Beckwith WH III (2002) *J. Clin. Microbiol.* **40**:3044-3046.
8. Colavita F, Quartu S, Lalle E, Bordi L, Lapa D, Meschi S, Vulcano A, Toffoletti A, Bordi E, Paglia MG, Di Caro A, Ippolito G, Capobianci MR, and Castilletti C. (2017) *J. Clin. Virol.* **86**: 27-30.

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' 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
www.cellbiolabs.com

©2021-2022: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.