



AnaSpec is pleased to introduce the industry's first commercially available assay kits for the detection of WNV protease NS3 inhibitors – the SensoLyte[®] series of WNV Protease Assay Kits.

West Nile virus (WNV), from the family Flaviviridae,¹ was first identified in the West Nile district of Uganda in 1937.² WNV outbreaks have been reported in Israel in the 50's, France in the 60's and South Africa in the 70's.³ In 1999, the first documented WNV infection in the US was reported in New York.⁴ The main route of human infection is through infected mosquito bites. WNV infection can cause severe neurological disease and fatalities in both human and animal hosts.

- ✓ Industry's First & Only
- ✓ Fluorogenic
- ✓ Sensitive
- ✓ Homogeneous
- ✓ Ideal for HTS of WNV Protease Inhibitor

WNV contains a single-stranded, positive-sense RNA genome, which encodes three structural proteins (capsid (C), membrane (M), envelope (E)), and seven non-structural proteins (NS1, NS2A, NS2B, NS3, NS4A, NS4B, NS5).^{5, 6} NS3 protease is essential (along with viral-encoded cofactor NS2B) for post-translational processing of a viral polypeptide precursor in infected host cells. This polypeptide provides the structural and functional viral proteins. Inhibition of its processing could represent a potential treatment for viral infections. With no effective vaccine or antiviral drug to protect against WNV infection,⁷ this protease represents a potentially key target for developing anti-WNV drugs.^{8, 9}

SensoLyte[®] West Nile Virus Protease Assay Kits provide a convenient, homogeneous assay for high throughput screening of West Nile Virus protease NS3 inhibitors. Utilizing a fluorogenic peptide (**SensoLyte[®] 440 West Nile Virus Protease Assay Kit**) or a FRET peptide (**SensoLyte[®] 570 West Nile Virus Protease Assay Kit**), these assays provide continuous quantification of protease activity.¹⁰ Upon NS3 protease cleavage, the fluorescence of AMC in the former and 5-TAMRA in the latter is recovered. Their fluorescence can be monitored at their characteristic emission wavelength.

SensoLyte[®] West Nile Virus Protease Assay Kits:
SensoLyte[®] 440 West Nile Virus Protease Assay Kit (Ex/Em=354/442 nm), Cat# 72079
SensoLyte[®] 570 West Nile Virus Protease Assay Kit (Ex/Em=540/575 nm), Cat# 72080

Related Product:
WNV NS3 Protease, recombinant, Cat# 72081

1. Prepare working solutions:

- Fluorogenic substrate (Component A)
- WNV protease (not provided)
- WNV inhibitor (Component D)

2. Set up enzymatic reaction (protease+ inhibitor, controls), 37C for 10 min.

3. Add diluted fluorogenic substrate, mix and start reading fluorescence signal (kinetic or end-point).

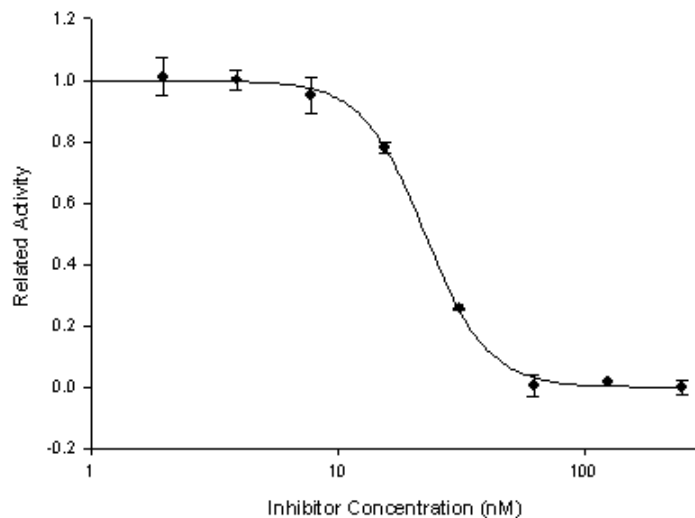


Figure 1. Inhibition of WNV by protease inhibitor, undeca-D-Arg-NH₂ using the Sensolyte[®] 440 WNV Protease Assay Kit, Cat# 72079.

Related Products: *Dengue Viral peptides*- contact service@anaspec.com

References:

1. Hayes, CG. *Ann. N.Y. Acad. Sci* 951, 25 (2001)
2. Smithburn, KC, et al. *Am J Trop Med.* 20, 471 (1940).
3. Internet. <http://www.nmnh.si.edu/BIRDNET/WNV.html>, accessed April 13, 2007.
4. Klee, AL. et al. *Emerg Infect Dis.* 10, 1405 (2004).
5. Brinton, MA. *Annu. Rev Microbiol.* 56, 371 (2002).
6. Lanciotti, RS. et al., *Science* 286, 2333 (1999).
7. Van Der Meulen, KM et al., *Arch. Virol.* 150, 637 (2005).
8. Mueller, NH. et al., *Int. J Biochem. Cell Biol.* 39(3), 606 (2007).
9. Shiryayev, SA. et al., *Biochem. J* 393, 503 (2006).
10. Chappell, KJ. et al., *J Biol. Chem.* 281, 38448 (2006).