

APPLICATION NOTE

The Naveni® Technology

A trustworthy neuroscience tool

Naveni® Proximity Ligation Assay in neuroscience research

The world of neuroscience is intricate and requires cutting-edge tools for the discovery of synaptic interactions and protein dynamics holding the key to the secrets of neurodegenerative diseases, cognitive functions, and even behavioral disorders. The Naveni® Technology, based on proximity ligation, is designed to explore protein-protein interactions with unparalleled precision and has been instrumental in several neuroscience publications.

Use case 1.

In a study by Vitet et al., the Naveni® Proximity Ligation method demonstrated the interaction between HTT and KIF1A, a protein interaction previously unconfirmed. Researchers found that neurons expressing constitutive phospho-HTT exhibited a significantly higher number of HTT/KIF1A interactions compared to wild type (WT). This discovery unveiled a critical pathway regulating synaptic transmission and motor skill learning. See figure 1.

Use case 2.

In another study led by Ross et al. the Naveni® Proximity Ligation Technology was used to investigate the dynamic interaction between AMPA receptors and the synaptic scaffolding protein Shank3. Shank3 is a master regulator

of synaptic architecture, intricately involved in regulating synaptic receptor density and, consequently, synaptic strength. Genetic associations extend its relevance to autism spectrum disorders.

Ross et al. explored the stability of the GluA1-Shank3 interaction under conditions of prolonged neuronal depolarization induced by elevated potassium levels. The results were intriguing: chronic depolarization led to a reduction in GluA1-Shank3 interactions, revealing the modulatory role of neuronal activity on this crucial synaptic interaction.

The benefits of the Naveni® Proximity Ligation Assay for neuroscience

Findings like these highlight the potential of proximity ligation assays in neuroscience research and the power of spatial interactomics. The sensitivity and specificity of the Naveni® Technology enables a deep dive into the molecular intricacies of synaptic connections, uncovering signaling pathways and protein interplay. Visualizing function is pivotal in understanding how proteins orchestrate synaptic plasticity, and how disturbances in the formation of these dynamic protein complexes contribute to neurological disorders.

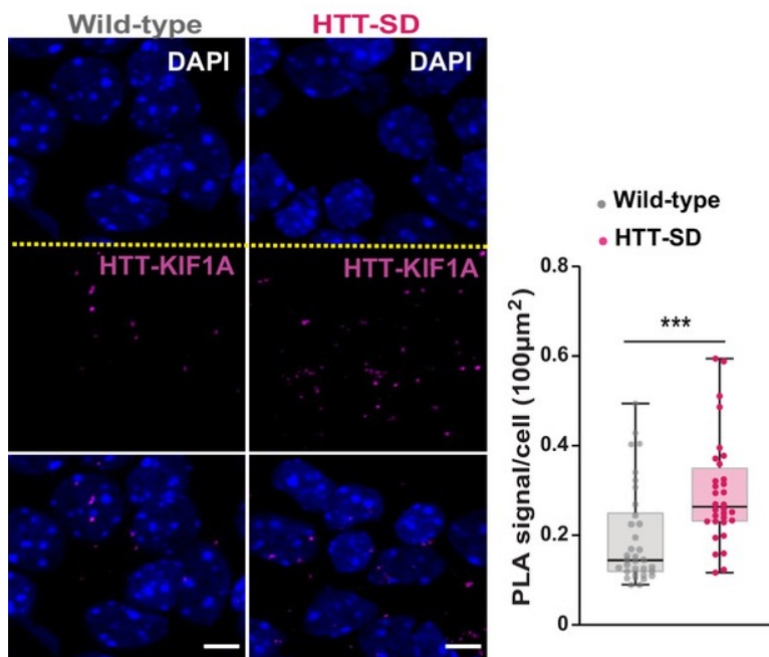
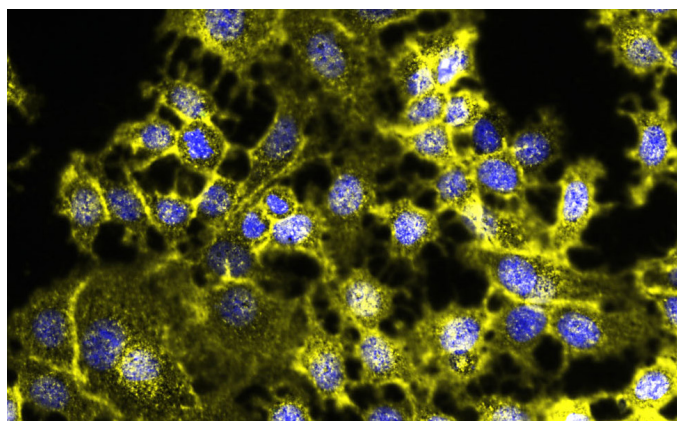


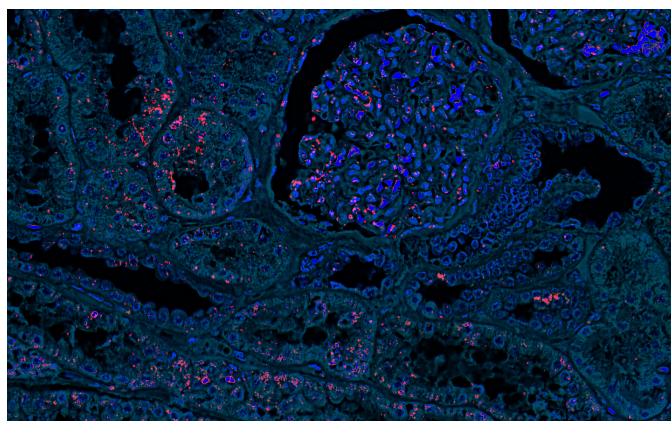
Figure 1. Naveni® Proximity Ligation Assay was used to confirm the cellular interaction between HTT and KIF1A. From Vitet et al (2023), eLife 12:e81011.

The NaveniFlex Cell and Tissue kits include two Navenibodies conjugated to proprietary oligo arms (depicted as orange antibodies in the illustration above). Only when the Navenibodies are in close proximity will they generate a rolling circle amplification reaction, resulting in a strong and distinct dot-like signal that is easily quantifiable.

NaveniFlex™ Cell



NaveniFlex™ Tissue



Selection of publications

Vitet Hélène, Bruyère Julie, Xu, Hao, Séris Claire, Brocard Jacques, et al. (2023) *loHuntingtin recruits KIF1A to transport synaptic vesicle precursors along the mouse axon to support synaptic transmission and motor skill learning.* Cambridge: eLife Sciences Publications Ltd. [DOI:10.7554/eLife.81011](https://doi.org/10.7554/eLife.81011)

Madeleine Ross, Elias Aizenman. *GluA1-Shank3 interaction decreases in response to chronic neuronal depolarization.* Neuroscience Letters, 2023, 137305, ISSN 0304-3940, [DOI:10.1016/j.neulet.2023.137305](https://doi.org/10.1016/j.neulet.2023.137305)

Estela Díaz-Sánchez a b, Alexander López-Salas a, Marina Mirchandani-Duque a, Jose Erik Alvarez-Contino a, Jose Andrés Sánchez-Pérez c, Kjell Fuxe d, Dasiel O. Borroto-Escuela a d e, Natalia García-Casares a, Manuel Narváez a b d. *Decreased medial prefrontal cortex activity related to impaired novel object preference task performance following GALR2 and Y1R agonists intranasal infusion.* Biomedicine & Pharmacotherapy. Volume 161, May 2023, 114433. [DOI: 10.1016/j.biopha.2023.114433](https://doi.org/10.1016/j.biopha.2023.114433)

Athanasios S. Alexandris, Jiwon Ryu, Labchan Rajbhandari, Robert Harlan, James McKenney, Yiqing Wang, Susan Aja, David Graham, Arun Venkatesan, Vassilis E. Koliatsos. *Protective effects of NAMPT or MAPK inhibitors and NaR on Wallerian degeneration of mammalian axons.* Neurobiology of Disease, Volume 171, 2022, [DOI:10.1016/j.nbd.2022.10580](https://doi.org/10.1016/j.nbd.2022.10580)



Distributed by:
Cayman Chemical
1180 E. Ellsworth Rd.
Ann Arbor, MI 48108 USA

Customer Service: (800) 364-9897
Technical Support: (888) 526-5351
www.caymanchem.com