

Product Information



NMDA Receptor NR2B Subunit (Phospho-Tyr¹³³⁶) Polyclonal Antibody

Item No. 10009760

Supplied as:	100 µl of affinity-purified antibody in 10 mM HEPES, pH 7.5, containing 150 mM NaCl, 100 µg/ml BSA, and 50% glycerol.
Host:	Rabbit
Antigen:	Phosphopeptide corresponding to amino acid residues surrounding Phospho-Tyr ¹³³⁶ of the human NMDA Receptor NR2B Subunit
Cross Reactivity:	(+) Human and rat NMDA Receptor; expected to react with bovine, canine, chicken, mouse, non-human primates and zebrafish NMDA Receptor based on 100% homology with the amino acid sequence used as the antigen
Stability:	≥1 year at -20°C
Application:	The recommended starting dilution for western blot is 1:1,000.

The NMDA receptor (NMDAR) plays an essential role in memory, neuronal development, and it has also been implicated in several disorders of the central nervous system including Alzheimer's disease, epilepsy, and ischemic neuronal death.¹⁻³ The NR1 protein can form NMDA activated channels when expressed in *Xenopus* oocytes but the currents in such channels are much smaller than those seen *in situ*. Channels with more physiological characteristics are produced when the NR1 subunit is combined with one or more of the NMDAR2 (NR2 A-D) subunits.⁴ Phosphorylation of Tyr¹³³⁶ in NR2B is thought to potentiate NMDA receptor-dependent influx of calcium.⁵ Ischemia may also increase the phosphorylation of this site.

References

1. Grosshans, D.R., Clayton, D.A., Coultrap, S.J., *et al.* LTP leads to rapid surface expression of NDMA but not AMPA receptors in adult rat CA1. *Nat. Neurosci.* **5**, 27-33 (2002).
2. Wenthold R.J., Prybylowski, K., Standley, S., *et al.* Trafficking of NMDA receptors. *Annu. Rev. Pharmacol Toxicol* **43**, 335-358 (2003).
3. Carroll, R.C., Zukin, R.S. NMDA-receptor trafficking and targeting: implications for synaptic transmission and plasticity. *Trends Neurosci.* **25**, 571-577 (2002).
4. Ishii, T., Moriyoshi, K., Sugihara, H., *et al.* Molecular characterization of the family of the N-methyl-D-aspartate receptor subunits. *J. Biol. Chem.* **268**, 2836-2843 (1993).
5. Takasu, M.A., Dalva, M.B., Zigmond, R.E., *et al.*, Modulation of NMDA receptor -dependent calcium influx and gene expression through EphB receptors. *Science* **295**, 491-495 (2002).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10009760

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 12/09/2011