

Product Information



GluR1 (Phospho-Ser⁸³¹) Polyclonal Antibody

Item No. 10602

Supplied as:	100 µl in 10 mM HEPES, pH 7.5, 150 mM NaCl, 100 µg/ml BSA, and 50% glycerol.
Synonym:	Glutamate Receptor Subunit 1
Host:	Rabbit
Antigen:	Phosphopeptide corresponding to amino acid residues surrounding the phospho-Ser ⁸⁴⁵ of GluR1
Cross Reactivity:	(+) rat GluR1
Stability:	≥1 year at -20°C
Application:	The recommended starting dilution for western blot is 1:1,000

The ion channels activated by glutamate are typically divided into two classes. Those that are sensitive to N-methyl-D-aspartate (NMDA) are designated NMDA receptors (NMDAR) while those activated by α -amino-3-hydroxy-5-methyl-4-isoxalone propionic acid (AMPA) are known as AMPA receptors (AMPA). The AMPAR are comprised of four distinct glutamate receptor subunits designated (GluR1-4) and they play key roles in virtually all excitatory neurotransmission in the brain. The GluR1 subunit is widely expressed throughout the nervous system. GluR1 is potentiated by phosphorylation at Ser⁸³¹ which has been shown to be mediated by either PKC or CaM kinase II. In addition, phosphorylation of this site has been linked to synaptic plasticity as well as learning and memory.

Related Products

GABAA Receptor δ subunit (N-Term) Polyclonal Antibody - Item No. 10600 • GluR1 (Phospho-Ser⁸⁴⁵) Polyclonal Antibody - Item No. 10601 • DARPP-32 (Phospho-Thr³⁴) Polyclonal Antibody - Item No. 10603 • Tyrosine Hydroxylase Polyclonal Antibody - Item No. 10604 • Synapsin I Polyclonal Antibody (neat serum) - Item No. 10605 • Synapsin I Polyclonal Antibody - Item No. 10606 • NMDA Receptor NR1 Subunit Monoclonal Antibody - Item No. 10607 • NMDA Receptor NR2A Subunit Polyclonal Antibody - Item No. 10608 • NMDA Receptor NR2B Subunit Polyclonal Antibody - Item No. 10609

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, 08/23/2010