

PRODUCT INFORMATION



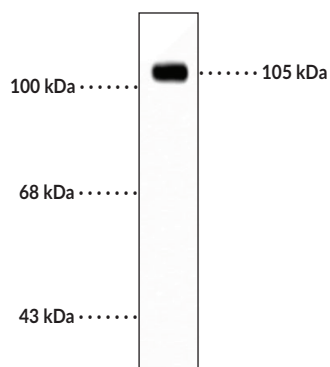
GluR1 Monoclonal Antibody (Clone RH95)

Item No. 29277

Overview and Properties

Contents:	This vial contains 100 μ l of protein G-purified mouse monoclonal antibody.
Synonyms:	AMPA 1, GluA1, Glutamate Receptor 1, GRIA1
Immunogen:	Peptide corresponding to amino acid residues from the N-terminal region of rat GluR1
Molecular Weight:	~105 kDa
Species Reactivity:	(+) Mouse, rat
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	\geq 1 year
Storage Buffer:	10 mM HEPES, pH 7.5, with 150 mM sodium chloride, 100 μ g/ml BSA, and 50% glycerol
Clone:	RH95
Host:	Mouse
Isotype:	IgG2a
Applications:	Western blot (WB); the recommended starting dilution is 1:1,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



WB of rat hippocampal lysate showing specific immunolabeling of the ~105 kDa GluR1 protein.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/30/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

AMPA receptors are ionotropic glutamate receptors that mediate excitatory synaptic transmission.^{1,2} They are tetrameric protein complexes expressed throughout the central nervous system in both neurons and glia that are assembled from combinations of GluR1, GluR2, GluR3, and GluR4, also known as GluR-A-D, subunits, each of which has extracellular N-terminal and ligand binding domains, a channel domain consisting of three membrane-spanning helices and a channel pore loop, and an intracellular C-terminus.²⁻⁵ The GluR1 subunit, which is encoded by *GRIA1* in humans, is primarily expressed in the forebrain and hippocampus. Mice lacking the GluR1 subunit exhibit normal spatial reference memory but have deficits in spatial working memory.¹ They also exhibit increased and decreased duration of nocifensive behaviors in phases I and II, respectively, of the formalin test, as well as increased duration of nocifensive behaviors in response to intraplantar injection of capsaicin (Item No. 92350), compared with wild-type mice.⁶ SNPs in *GRIA1* have been found in patients with bipolar disorder with psychotic features.⁴ Elevated levels of *GRIA1* mRNA have been found in postmortem dorsolateral prefrontal cortex from patients with schizophrenia.⁷ Cayman's GluR1 Monoclonal Antibody (Clone RH95) can be used for Western blot (WB) applications. The antibody recognizes GluR1 at approximately 105 kDa from mouse and rat samples.

References

1. Sanderson, D.J., Good, M.A., Seeburg, P.H., *et al.* The role of the GluR-A (GluR1) AMPA receptor subunit in learning and memory. *Progress in Brain Research: Essence of Memory*. Sossin, W., Lacaille, J.-C., Castellucci, V.F., *et al.*, editors, Elsevier (2008).
2. Isaac, J.T.R., Ashby, M.C., and McBain, C.J. The role of the GluR2 subunit in AMPA receptor function and synaptic plasticity. *Neuron* **54(6)**, 859-871 (2007).
3. Bassani, S., Valnegri, P., Beretta, F., *et al.* The GLUR2 subunit of AMPA receptors: Synaptic role. *Neuroscience* **158(1)**, 55-61 (2009).
4. Kerner, B., Jasinska, A.J., DeYoung, J., *et al.* Polymorphisms in the *GRIA1* gene region in psychotic bipolar disorder. *Am. J. Med. Genet. B Neuropsychiatr. Genet.* **150B(1)**, 24-32 (2009).
5. Sprengel, R. Role of AMPA receptors in synaptic plasticity. *Cell Tissue Res.* **326(2)**, 447-455 (2006).
6. Hartmann, B., Ahmadi, S., Heppenstall, P.A., *et al.* The AMPA receptor subunits GluR-A and GluR-B reciprocally modulate spinal synaptic plasticity and inflammatory pain. *Neuron* **44(4)**, 637-650 (2004).
7. O'Connor, J.A. and Hemby, S.E. Elevated *GRIA1* mRNA expression in layer II/III and V pyramidal cells of the DLPFC in schizophrenia. *Schizophr. Res.* **97(1-3)**, 277-288 (2007).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM