

# PRODUCT INFORMATION



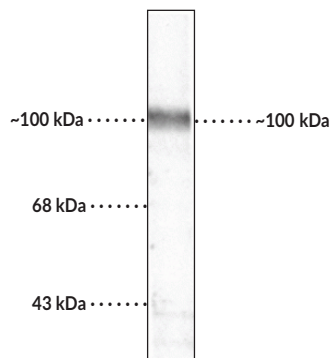
## GluR2 Polyclonal Antibody

Item No. 29278

### Overview and Properties

<b>Contents:</b>	This vial contains 100 µl of affinity-purified polyclonal antibody from pooled serum.
<b>Synonyms:</b>	AMPA 2, AMPA-selective Glutamate Receptor 2, GluA2, Glutamate Ionotropic Receptor AMPA Type Subunit 2, GRIA2
<b>Immunogen:</b>	Peptide corresponding to amino acid residues from rat GluR2, conjugated to keyhole limpet hemocyanin (KLH)
<b>Molecular Weight:</b>	~100 kDa
<b>Species Reactivity:</b>	(+) Rat
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	10 mM HEPES, pH 7.5, with 150 mM sodium chloride, 100 µg/ml BSA, and 50% glycerol
<b>Host:</b>	Rabbit
<b>Applications:</b>	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 a rat hippocampal lysate showing the specific immunolabeling of the ~100 kDa GluR2 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, 07/08/2020

**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.<sup>1,2</sup> They are tetrameric protein complexes expressed throughout the CNS 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.<sup>1,3-5</sup> GluR1 and GluR2 are predominantly expressed in the forebrain with low levels of GluR3 and GluR4, and pyramidal neurons express AMPA receptors primarily comprised of heterotetramers of GluR1 and GluR2 subunits. GluR2 is encoded by *GRIA2* and regulates AMPA receptor calcium permeability, assembly, trafficking, and single channel conductance.<sup>1,6</sup> It undergoes RNA editing at position 586, the Q/R site, to introduce an arginine residue that renders GluR2 subunit-containing AMPA receptors impermeable to calcium.<sup>1,6,7</sup> Motor neurons isolated from *Gria2*-deficient mouse embryos exhibit increased calcium influx and increased susceptibility to AMPA receptor-mediated excitotoxicity.<sup>7</sup> Heterozygous knockdown of *Gria2* accelerates motor neuron degradation and decreases lifespan in the SOD1G93A transgenic mouse model of amyotrophic lateral sclerosis. *GRIA2* expression is reduced in postmortem hippocampus from patients with chronic schizophrenia.<sup>8</sup> Cayman's GluR2 Polyclonal Antibody can be used for Western blot (WB) applications. The antibody detects GluR2 at approximately 100 kDa from rat samples.

## References

---

1. 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).
2. 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).
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. Duprat, F., Daw, M., Lim, W., et al. GluR2 protein-protein interactions and the regulation of AMPA receptors during synaptic plasticity. *Philos. Trans. R. Soc. Lond. B Biol. Sci.* **358(1432)**, 715-720 (2003).
7. Van Damme, P., Braeken, D., Callewaert, G., et al. GluR2 deficiency accelerates motor neuron degeneration in a mouse model of amyotrophic lateral sclerosis. *J. Neuropathol. Exp. Neurol.* **64(7)**, 605-612 (2005).
8. Eastwood, S.L., Burnet, P.W.J., and Harrison, P.J. GluR2 glutamate receptor subunit flip and flop isoforms are decreased in the hippocampal formation in schizophrenia: A reverse transcriptase-polymerase chain reaction (RT-PCR) study. *Brain Res. Mol. Brain Res.* **44(1)**, 92-98 (1997).

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