

PRODUCT INFORMATION



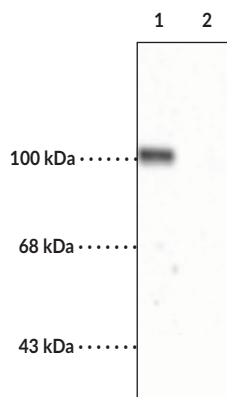
GluR1 (Phospho-Ser845) Polyclonal Antibody

Item No. 10601

Overview and Properties

Contents:	This vial contains affinity-purified IgG polyclonal antibody from pooled serum.
Synonyms:	AMPA-selective Glutamate Receptor 1, GluR-1, GluR-A, GluA1
Immunogen:	Phosphopeptide corresponding to amino residues surrounding the phospho-Ser ⁸⁴⁵ of rat GluR1
Molecular Weight:	~100 kDa
Cross Reactivity:	(+) GluR1 (phospho-Ser ⁸⁴⁵); (-) Non-phosphorylated GluR1
Species Reactivity:	(+) Mouse, rat, turtle
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	10 mM HEPES, pH 7.5, with 150 mM sodium chloride, 100 µg/ml BSA, and 50% glycerol
Host:	Rabbit
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution for IHC and WB is 1:1,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: - λ-Ptase
Lane 2: + λ-Ptase

WB of rat hippocampal lysate showing specific immunolabeling of the ~100 kDa GluR1 protein phosphorylated at Ser⁸⁴⁵ in the first lane (-). Phosphospecificity is shown in the second lane (+) where immunolabeling is completely eliminated by blot treatment with lambda phosphatase (λ-Ptase, 1200 units for 30 min).

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, 11/16/2020

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

GluR1 is a subunit of the AMPA ionotropic glutamate receptor, which is responsible for fast excitatory synaptic transmission in the CNS.¹ AMPA receptors are composed of four subunits, GluR1, GluR2, GluR3, and GluR4, which combine into heterotetramers to form a cation-permeable pore in the plasma membrane. Each subunit has two isoforms with the primary isoform designated as flip and a second isoform generated through alternative splicing designated as flop.^{2,3} The GluR1 flip and flop isoforms do not affect desensitization or channel opening and closing kinetics.³ GluR1 can be phosphorylated by PKA at serine 845 (Ser⁸⁴⁵), which increases the peak open probability of the ion channel, and dephosphorylation is required for AMPA receptor endocytosis.¹ GluR1 (phospho-Ser⁸⁴⁵) levels increase in the ventromedial prefrontal cortex (vmPFC) and nucleus accumbens (NAc) following cocaine-cue extinction.⁴ Levels of GluR1 (phospho-Ser⁸⁴⁵) are reduced following NMDA receptor activation in rat hippocampal slices and by amyloid- β (β) oligomers in primary mouse neurons.^{5,6} Cayman's GluR1 (Phospho-Ser⁸⁴⁵) Polyclonal Antibody can be used for immunohistochemistry (IHC) and Western blot (WB) applications. The antibody recognizes GluR1 (phospho-Ser⁸⁴⁵) at approximately 100 kDa from mouse, rat, and turtle samples.

References

1. Banke, T.G., Bowie, D., Lee, H.-K., *et al.* Control of GluR1 AMPA receptor function by cAMP-dependent protein kinase. *J. Neurosci.* **20**(1), 89-102 (2000).
2. Sommer, B., Keinänen, K., Verdoorn, T.A., *et al.* Flip and flop: A cell-specific functional switch in glutamate-operated channels of the CNS. *Science* **249**(4976), 1580-1585 (1990).
3. Pei, W., Huang, Z., Wang, C., *et al.* Flip and flop: A molecular determinant for AMPA receptor channel opening. *Biochemistry* **48**(17), 3767-3777 (2009).
4. Dhonnchadha, B.Á.N., Lin, A., Leite-Morris, K.A., *et al.* Alterations in expression and phosphorylation of GluA1 receptors following cocaine-cue extinction learning. *Behav. Brain Res.* **238**, 119-123 (2013).
5. Ai, H., Yang, W., Ye, M., *et al.* Differential regulation of AMPA receptor GluA1 phosphorylation at serine 831 and 845 associated with activation of NMDA receptor subpopulations. *Neurosci. Lett.* **497**(2), 94-98 (2011).
6. Miñano-Molina, A.J., España, J., Martín, E., *et al.* Soluble oligomers of amyloid- β peptide disrupt membrane trafficking of α -amino-3-hydroxy-5-methylisoxazole-4-propionic acid receptor contributing to early synapse dysfunction. *J. Biol. Chem.* **286**(31), 27311-27321 (2011).

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