

GRIA3 Polyclonal Antibody

(Catalog # A70499)

Background

Glutamate receptors are the predominant excitatory neurotransmitter receptors in the mammalian brain and are activated in a variety of normal neurophysiologic processes. These receptors are heteromeric protein complexes composed of multiple subunits, arranged to form ligand-gated ion channels. The classification of glutamate receptors is based on their activation by different pharmacologic agonists. The subunit encoded by this gene belongs to a family of AMPA (alpha-amino-3-hydroxy-5-methyl-4-isoxazole propionate)-sensitive glutamate receptors, and is subject to RNA editing (AGA->GGA; R->G). Alternative splicing at this locus results in different isoforms, which may vary in their signal transduction properties.

Description

GRIA3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P42263

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 200-490 of human GRIA3 (NP_000819.3).

Storage

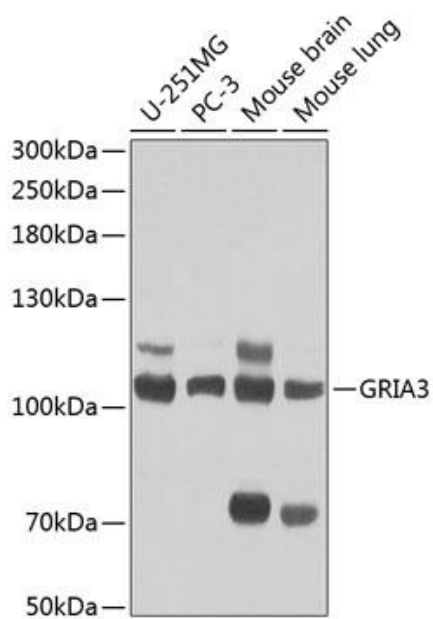
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

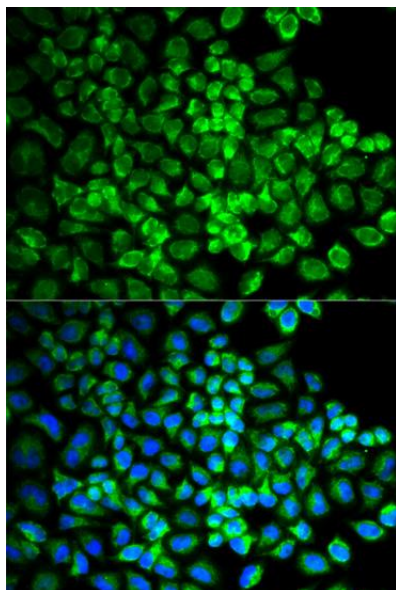
GRIA3; GLUR-C; GLUR-K3; GLUR3; GLURC; GluA3; MRX94; glutamate receptor 3

Application

WB, IHC, IF; Recommended dilution: WB: 1:500-1:2000, IHC: 1:50-1:200, IF:1:50-1:200



Western blot analysis of extracts of various cell lines, using GRIA3 Polyclonal Antibody at 1:1000 dilution. Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug per lane. Blocking buffer: 3% nonfat dry milk in TBST.



Immunofluorescence analysis of U2OS cells using GRIA3 Polyclonal Antibody. Blue: DAPI for nuclear staining.