

# Grik2 (C2) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Grik2 (C2)

## Instruction Manual

Catalog Number	PK-AB718-4387
Synonyms	Grik2 Antibody: Glutamate receptor ionotropic kainate 2, glutamate receptor 6, gluR6, excitatory amino acid receptor 4, eea4
Description	Glutamate receptors are the predominant excitatory neurotransmitter receptors in the mammalian brain and are activated in a variety of normal neurophysiologic processes. Grik2, also known as glutamate receptor 6, belongs to the kainate family of glutamate receptors, which are composed of four subunits and function as ligand-activated ion channels. Recent reports have suggested that defects in the Grik2 protein may be associated with autosomal recessive mental retardation and possibly other neurological disorders such as schizophrenia. Numerous isoforms of Grik2 are known to exist and may be subject to RNA editing within the second transmembrane domain, which is thought to alter the properties of ion flow. This Grik2 antibody may exhibit some cross-reactivity to Grik3.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Grik2 antibody was raised in rabbits against a 13 amino acid peptide near the carboxy terminus of the human Grik2.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB <b>Note:</b> Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Grik2 antibody can be used for detection of Grik2 by Western blot at 0.5 - 1 µg/mL.
Images	Available upon request.
References	Tanaka K. Functions of glutamate transports in the brain. <i>Neurosci. Res.</i> 2000; 37:15-9. Pinheiro P and Mulle C. Kainate receptors. <i>Cell Tissue Res.</i> 2006; 326:457-82. Paschen W, Blackstone CD, Haganir RL, et al. Human GluR6 kainate receptor (GRIK2): molecular cloning, expression, polymorphism, and chromosomal assignment. <i>Genomics</i> 1994; 20:435-40. Motazacker MM, Rost BR, Hucho T, et al. A defect in the ionotropic glutamate receptor 6 gene (GRIK2) is associated with autosomal recessive mental retardation. <i>Am. J. Hum. Genet.</i> 2007; 81:792-8.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4387P; Grik2 Peptide Cat.No. PK-AB718-1303; Human Brain Tissue Lysate

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.