

# LGI3 (IN) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat LGI3 (IN)

## Instruction Manual

Catalog Number	PK-AB718-4493
Synonyms	LGI3 Antibody: Leucine-rich, glioma inactivated 3
Description	The leucine-rich, glioma inactivated gene 3 (LGI3) is a member of the LGI family in which LGI1 is the exemplar. The LGI family consists of four of highly related proteins containing leucine-rich repeats (LRRs) which are highly similar to other transmembrane signaling molecules and receptors. LGI1 has been identified as a candidate tumor suppressor gene for glioma and plays a role in autodominate lateral temporal epilepsy (ADTLE), an epileptic syndrome characterized by focal seizures with predominant auditory symptoms. Despite its high homology with LGI1 and similar pattern of expression, mutations in LGI3 have not been found to be associated with ADTLE. LGI3 expression is induced in rat astrocyte cultures by the amyloid beta (Ab) peptide and accumulated on neuronal plasma membranes of aged monkey brains and co-localized with Ab. Two isoforms of LGI3 are known to exist; this LGI3 antibody will recognize both. This LGI3 antibody is predicted to be specific to LGI3 and not recognize other LGI proteins. The observed higher molecular weight band may represent a post-translationally modified form of LGI3.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	LGI3 antibody was raised in rabbits against a 12 amino acid peptide from near the center of human LGI3.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	Two isoforms of LGI3 are known to exist; this LGI3 antibody will recognize both. This LGI3 antibody is predicted to be specific to LGI3 and not recognize other LGI proteins. The observed higher molecular weight band may represent a post-translationally modified form of LGI3.
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 [Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.]  LGI3 antibody can be used for the detection of LGI3 by Western blot at 1 µg/mL.
Images	Available upon request.
References	Gu W, Gibert Y, Wirth T, et al. Using gene-history and expression analysis to assess the involvement of LGI genes in human disorders. <i>Mol. Biol. Evol.</i> 2005; 22:2209-16. Chernova OB, Somerville RP and Cowell JK. A novel gene, LGI1, from 10q24 is rearranged and downregulated in malignant brain tumors. <i>Oncogene</i> 1998; 17:2873-81. Berkovic SF, Izzillo P, McMahon JM, et al. LGI1 mutations in temporal lobe epilepsies. <i>Neurology</i> 2004; 62:1115-9. Kimura N, Ishii Y, Suzuki S, et al. Abeta upregulates and colocalizes with LGI3 in cultured rat astrocytes. <i>Cell Mol. Neurobiol.</i> 2007; 27:335-50.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4493P; LGI3 Peptide Cat.No. PK-AB718-1303; Human Brain Tissue Lysate

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