

LGI2 (IN) antibody (pAb)

Rabbit Anti-Human/Mouse LGI2 (IN)

Instruction Manual

Catalog Number	PK-AB718-4533
Synonyms	LGI2 Antibody: Leucine-rich, glioma inactivated 2, LGIL2
Description	The leucine-rich, glioma inactivated gene 2 (LGI2) 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 LGI2 have not been found to be associated with ADTLE. Three isoforms of LGI2 are known to exist; this LGI2 antibody will recognize only the longest. This LGI2 antibody is predicted to be specific to LGI2 and not recognize other LGI proteins.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	LGI2 antibody was raised in rabbits against a 10 amino acid peptide from near the center of human LGI2.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
Specificity	This LGI2 antibody is predicted to be specific to LGI2 and not recognize other LGI proteins. Three isoforms of LGI2 are known to exist; this LGI2 antibody will recognize only the longest.
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, IHC, IF I Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. LGI2 antibody can be used for the detection of LGI2 by Western blot at 1 - 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL. For immunofluorescence start at 5 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.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4533P; LGI2 Peptide Cat.No. PK-AB718-1404; Mouse Liver Tissue Lysate

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