

LGI1 (IN) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat LGI1 (IN)

Instruction Manual

Catalog Number	PK-AB718-4489
Synonyms	LGI1 Antibody: Leucine-rich, glioma inactivated 1, epitempin, EPT
Description	The leucine-rich, glioma inactivated gene 1 (LGI1) was first identified as a candidate tumor suppressor gene for glioma and may play a role in other cancers. LGI1 is a member of a family of highly related proteins containing leucine-rich repeats (LRRs) which are highly similar to other transmembrane signaling molecules and receptors. LGI1 serves as a ligand to ADAM22, a metalloprotease localized at the synapse. Mutations in LGI1 account for nearly half of autodominate lateral temporal epilepsy (ADTLE), an epileptic syndrome characterized by focal seizures with predominant auditory symptoms. Two isoforms of LGI1 are known to exist; this LGI1 antibody will recognize only the longer form. This LGI1 antibody is predicted to be specific to LGI1 and not recognize other LGI proteins.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	LGI1 antibody was raised in rabbits against a 14 amino acid peptide from near the center of human LGI1.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	This LGI1 antibody is predicted to be specific to LGI1 and not recognize other LGI proteins.
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 INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. LGI1 antibody can be used for the detection of LGI1 by Western blot at 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	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. Fialka F, Gruber RM, Hitt R, et al. CPA6, FMO2, LGI1, SIAT1 and TNC are differentially expressed in early- and late-stage oral squamous cell carcinoma – A pilot study. <i>Oral Oncol.</i> 2008; 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. Fukata Y, Adesnik H, Iwanaga T, et al. Epilepsy-related ligand/receptor complex LGI1 and ADAM22 regulate synaptic transmission. <i>Science</i> 2006; 313:1792-5.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4489P; LGI1 Peptide Cat.No. PK-AB718-1201; HeLa Cell Lysate

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