

GRID1 Antibody [#C12394]

| | |
|---------------------------|--|
| Catalog Number: | C12394 |
| Concentration: | 1mg/ml |
| Swiss-Prot No.: | Q9ULK0 |
| Other Names: | GRID1, GluR delta-1, glutamate receptor delta-1 subunit |
| Storage/Stability: | Store at -20°C/1 year |
| Form of Antibody: | Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol. |
| Immunogen: | The antiserum was produced against synthesized peptide derived from internal of human GRID1. |
| Purification: | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen. |
| Specificity: | GRID1 antibody detects endogenous levels of total GRID1 protein. |
| Reactivity: | Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%) |
| Applications: | WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000 |
| References: | The MGC Project Team, Genome Res. 14:2121-2127(2004). Nagase T., DNA Res. 6:337-345(1999). |