

GluR7 Antibody [#C0211]

Catalog Number: C0211-1, C0211-2, C0211

Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl

Swiss-Prot No.: Q14831

All Names: EAA5; GLR7; GluR7a, dJ1090M5.1 (glutamate receptor, ionotropic, kainate 3 (GLUR7)), excitatory amino acid receptor 5, glutamate receptor 7 precursor

Form of Antibody: Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.

Storage/Stability: Store at -20°C/1 year

Immunogen: The antiserum was produced against synthesized peptide derived from human GluR7.

Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.

Specificity/Sensitivity GluR7 antibody detects endogenous levels of total GluR7 protein.

Reactivity: Human, Mouse, Rat

Applications: WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000

References: Hans H. Schiffer, J. Neurosci., Dec 2000; 20: 9025.

Nathalie Strutz, J. Neurosci., Jan 2001; 21: 401.

Frédéric Jaskolski, J. Biol. Chem., Jun 2005; 280: 22968 - 22976.

DE Hamassaki-Britto, J. Neurosci., May 1993; 13: 1888.