

GluR4 (Ab-862) Antibody [#B0936]

Catalog Number:	B0936
Concentration:	1mg/ml
Swiss-Prot No.:	P48058
Other Names:	GLR4, GLUR4, GRIA4, GluR-4, GluR-D, GluR4, Glutamate receptor 4 precursor, Glutamate receptor ionotropic, AMPA 4
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 non-phosphopeptide derived from human GluR4 around the phosphorylation site of serine 862 (R-L-S ^P -I-T).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GluR4 (Ab-862) antibody detects endogenous levels of total GluR4 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Maxim Mokin, J Neurophysiol, Jan 2006; 95: 215 - 224. Alejandro Caicedo, Chem Senses, Jul 2004; 29: 463 - 471. LJ Martin, J. Neurosci., Feb 1993; 13: 782. Masahiro Kondo, J. Neurosci., Mar 1997; 17: 1570.