

GRM1 Antibody [#G363]

Catalog Number:	G363
Concentration:	1mg/ml
Swiss-Prot No.:	Q13255
Other Names:	Metabotropic glutamate receptor 1; mGluR1; GRM1; GPRC1A; MGLUR1
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 GRM1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GRM1 antibody detects endogenous levels of total GRM1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:5000
References:	Desai M.A., Mol. Pharmacol. 48:648-657(1995). Stephan D., Neuropharmacology 35:1649-1660(1996). Mungall A.J., Nature 425:805-811(2003).