

GALR2 Antibody [#G268]

Catalog Number:	G268
Concentration:	1mg/ml
Swiss-Prot No.:	O43603
Other Names:	GAL2-R; galanin receptor 2; GALNR2
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 GALR2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GALR2 antibody detects endogenous levels of total GALR2 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Bloomquist B.T., Biochem. Biophys. Res. Commun. 243:474-479(1998). Borowsky B., Peptides 19:1771-1781(1998). Fathi Z., Brain Res. Mol. Brain Res. 58:156-169(1998).