

GPR173 Antibody [#G144]

Catalog Number:	G144
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NS66
Other Names:	Probable G-protein coupled receptor 173; Super conserved receptor expressed in brain 3; GPR173; SREB3
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 GPR173.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GPR173 antibody detects endogenous levels of total GPR173 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Matsumoto M., Biochem. Biophys. Res. Commun. 272:576-582(2000). The MGC Project Team; Genome Res. 14:2121-2127(2004).