

GPR12 Antibody [#G323]

Catalog Number:	G323
Concentration:	1mg/ml
Swiss-Prot No.:	P47775
Other Names:	Probable G-protein coupled receptor 12
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 GPR12.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GPR12 antibody detects endogenous levels of total GPR12 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:5000
References:	Song Z.-H., Genomics 28:347-349(1995). Dunham A., Nature 428:522-528(2004). The MGC Project Team, Genome Res. 14:2121-2127(2004).