

GPR75 Antibody [#G351]

Catalog Number:	G351
Concentration:	1mg/ml
Swiss-Prot No.:	O95800
Other Names:	Probable G-protein coupled receptor 75; GPR75
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 GPR75.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GPR75 antibody detects endogenous levels of total GPR75 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
References:	Tarttelin E.E., Biochem. Biophys. Res. Commun. 260:174-180(1999). Hillier L.W., Nature 434:724-731(2005). The MGC Project Team; Genome Res. 14:2121-2127(2004).