

GPR31 Antibody [#G334]

Catalog Number:	G334
Concentration:	1mg/ml
Swiss-Prot No.:	O00270
Other Names:	Probable G-protein coupled receptor 31; GPR31
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 GPR31.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GPR31 antibody detects endogenous levels of total GPR31 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Zingoni A., Genomics 42:519-523(1997). Mungall A.J., Nature 425:805-811(2003).