

GPR176 Antibody [#G115]

Catalog Number:	G115
Concentration:	1mg/ml
Swiss-Prot No.:	Q14439
Other Names:	Probable G-protein coupled receptor 176; HB-954; GPR176
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 C-terminal of human GPR176.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GPR176 antibody detects endogenous levels of total GPR176 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:10000
References:	Hata S., Biochim. Biophys. Acta 1261:121-125(1995). The MGC Project Team; Genome Res. 14:2121-2127(2004).