

GPR110 Antibody [#G122]

Catalog Number:	G122
Concentration:	1mg/ml
Swiss-Prot No.:	Q5T601
Other Names:	Probable G-protein coupled receptor 110 [Precursor]; G-protein coupled receptor PGR19; G-protein coupled receptor KPG_012; PGR19
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 GPR110.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GPR110 antibody detects endogenous levels of total GPR110 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:40000
References:	Mungall A.J., Nature 425:805-811(2003). Fredriksson R., FEBS Lett. 531:407-414(2002). Ota T., Nat. Genet. 36:40-45(2004).