

GPR120 Antibody [#G127]

Catalog Number:	G127
Concentration:	1mg/ml
Swiss-Prot No.:	Q5NUL3
Other Names:	G-protein coupled receptor 120; G-protein coupled receptor PGR4; G-protein coupled receptor GT01; G-protein coupled receptor 129; GPR129, PGR4
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 GPR120.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GPR120 antibody detects endogenous levels of total GPR120 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
References:	Fredriksson R., FEBS Lett. 554:381-388(2003). Hirasawa A., Nat. Med. 11:90-94(2005). Deloukas P., Nature 429:375-381(2004).