

GPR115 Antibody [#G285]

Catalog Number:	G285
Concentration:	1mg/ml
Swiss-Prot No.:	Q8IZF3
Other Names:	FLJ38076; G-protein coupled receptor 115; G-protein coupled receptor PGR18; GP115; PGR18
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 GPR115.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GPR115 antibody detects endogenous levels of total GPR115 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Mungall A.J., Nature 425:805-811(2003). Vassilatis D.K., Proc. Natl. Acad. Sci. U.S.A. 100:4903-4908(2003). Fredriksson R., FEBS Lett. 531:407-414(2002).