

GPR114 Antibody [#G284]

Catalog Number:	G284
Concentration:	1mg/ml
Swiss-Prot No.:	Q8IZF4
Other Names:	Probable G-protein coupled receptor 114 [Precursor]; G-protein coupled receptor PGR27; GPR114; PGR27; UNQ2524/PRO6017
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 GPR114.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GPR114 antibody detects endogenous levels of total GPR114 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Clark H.F., <i>me Res.</i> 13:2265-2270(2003). Ota T., <i>Nat. Genet.</i> 36:40-45(2004). Vassilatis D.K., <i>Proc. Natl. Acad. Sci. U.S.A.</i> 100:4903-4908(2003).