

## GPR133 Antibody [#G292]

<b>Catalog Number:</b>	G292
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q6QNK2
<b>Other Names:</b>	Probable G-protein coupled receptor 133 [Precursor]; G-protein coupled receptor PGR25; PGR25
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human GPR133.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	GPR133 antibody detects endogenous levels of total GPR133 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Bonner T.I., Submitted (JAN-2004) to the EMBL/GenBank/DDBJ databases. Bjarnadottir T.K., Genomics 84:23-33(2004). Ota T., Nat. Genet. 36:40-45(2004).