

## GPR92 Antibody [#G360]

<b>Catalog Number:</b>	G360
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9H1C0
<b>Other Names:</b>	Probable G-protein coupled receptor 92; GPR92; GPR93
<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 GPR92.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	GPR92 antibody detects endogenous levels of total GPR92 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	White K.E., Nat. Genet. 26:345-348(2000). Lee D.K., Gene 275:83-91(2001). Takeda S., FEBS Lett. 520:97-101(2002).