

## GPR157 Antibody [#G306]

<b>Catalog Number:</b>	G306
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q5UAW9
<b>Other Names:</b>	Probable G-protein coupled receptor 157
<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 GPR157.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	GPR157 antibody detects endogenous levels of total GPR157 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Bonner T.I., Submitted (SEP-2004) to the EMBL/GenBank/DDBJ databases. Ota T., Nat. Genet. 36:40-45(2004). Gregory S.G., Nature 441:315-321(2006).