

GPR158 Antibody [#G307]

Catalog Number:	G307
Concentration:	1mg/ml
Swiss-Prot No.:	Q5T848
Other Names:	G-protein coupled receptor 158; GP158
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 GPR158.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GPR158 antibody detects endogenous levels of total GPR158 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	Bonner T.I., Submitted (JAN-2004) to the EMBL/GenBank/DDBJ databases. Deloukas P., Nature 429:375-381(2004). Hirosawa M., DNA Res. 6:329-336(1999).