

GPR22 Antibody [#G330]

Catalog Number:	G330
Concentration:	1mg/ml
Swiss-Prot No.:	Q99680
Other Names:	Probable G-protein coupled receptor 22
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 GPR22.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GPR22 antibody detects endogenous levels of total GPR22 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:5000
References:	O'Dowd B.F., Gene 187:75-81(1997). Hillier L.W., Nature 424:157-164(2003).