

GPR52 Antibody [#G344]

Catalog Number:	G344
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y2T5
Other Names:	Probable G-protein coupled receptor 52
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 C-terminal of human GPR52.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GPR52 antibody detects endogenous levels of total GPR52 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Sawzdargo M., Brain Res. Mol. Brain Res. 64:193-198(1999). Gregory S.G., Nature 441:315-321(2006). The MGC Project Team; Genome Res. 14:2121-2127(2004).