

MTR1L Antibody [#G400]

Catalog Number:	G400
Concentration:	1mg/ml
Swiss-Prot No.:	Q13585
Other Names:	Melatonin-related receptor; G protein-coupled receptor 50; H9; GPR50
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 MTR1L.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MTR1L antibody detects endogenous levels of total MTR1L protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:10000
References:	Reppert S.M., FEBS Lett. 386:219-224(1996).