

MRGRE Antibody [#G391]

Catalog Number:	G391
Concentration:	1mg/ml
Swiss-Prot No.:	Q86SM8
Other Names:	Mas-related G-protein coupled receptor member E; G-protein coupled receptor 167; MRGPRES; GPR167, MRGE
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 MRGRE.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRGRE antibody detects endogenous levels of total MRGRE protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	The MGC Project Team; Genome Res. 14:2121-2127(2004). Vassilatis D.K., Proc. Natl. Acad. Sci. U.S.A. 100:4903-4908(2003).