

MAS1L Antibody [#G383]

Catalog Number:	G383
Concentration:	1mg/ml
Swiss-Prot No.:	P35410
Other Names:	Mas-related G-protein coupled receptor MRG; MAS-R; MAS1-like; MAS1L; MRG
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 MAS1L.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MAS1L antibody detects endogenous levels of total MAS1L protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:5000
References:	Monnot C., Mol. Endocrinol. 5:1477-1487(1991). Mungall A.J., Nature 425:805-811(2003). The MGC Project Team, Genome Res. 14:2121-2127(2004).