

OR511 Antibody [#G641]

Catalog Number:	G641
Concentration:	1mg/ml
Swiss-Prot No.:	Q13606
Other Names:	HSOLF1; OLF1; olfactory receptor 5I1; olfactory receptor OR11-159; olfactory receptor, family 5, subfamily I, member 1; olfactory receptor-like OLF1
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 OR511.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OR511 antibody detects endogenous levels of total OR511 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:10000
References:	Issel-Tarver L., Genetics 145:185-195(1997). The MGC Project Team; Genome Res. 14:2121-2127(2004). Malnic B., Proc. Natl. Acad. Sci. U.S.A. 101:2584-2589(2004).