

OR13C3 Antibody [#G425]

Catalog Number:	G425
Concentration:	1mg/ml
Swiss-Prot No.:	Q8NGS6
Other Names:	O13C3; olfactory receptor 13C3; olfactory receptor, family 13, subfamily C, member 3; OR37G; OR9-8
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 OR13C3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OR13C3 antibody detects endogenous levels of total OR13C3 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:10000
References:	Suwa M., Submitted (JUL-2001) to the EMBL/GenBank/DDBJ databases. Humphray S.J., Nature 429:369-374(2004).