

## OR13C4 Antibody [#G426]

<b>Catalog Number:</b>	G426
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q8NGS5
<b>Other Names:</b>	Olfactory receptor 13C4; Olfactory receptor OR9-7
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human OR13C4.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	OR13C4 antibody detects endogenous levels of total OR13C4 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Humphray S.J., Nature 429:369-374(2004). Fuchs T., Genomics 80:295-302(2002). Malnic B., Proc. Natl. Acad. Sci. U.S.A. 101:2584-2589(2004).