

OR4C13 Antibody [#G588]

Catalog Number:	G588
Concentration:	1mg/ml
Swiss-Prot No.:	Q8NGP0
Other Names:	Olfactory receptor 4C13; Olfactory receptor OR11-260
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 C-terminal of human OR4C13.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OR4C13 antibody detects endogenous levels of total OR4C13 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
References:	Suwa M., Submitted (JUL-2001) to the EMBL/GenBank/DDBJ databases. Fuchs T., Genomics 80:295-302(2002). Malnic B., Proc. Natl. Acad. Sci. U.S.A. 101:2584-2589(2004).