

OR2D3 Antibody [#G546]

Catalog Number:	G546
Concentration:	1mg/ml
Swiss-Prot No.:	Q8NGH3
Other Names:	olfactory receptor 2D3; olfactory receptor OR11-89; olfactory receptor, family 2, subfamily D, member 3; OR11-89
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 OR2D3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OR2D3 antibody detects endogenous levels of total OR2D3 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. Fuchs T., Genomics 80:295-302(2002).