

OR2Y1 Antibody [#G574]

Catalog Number:	G574
Concentration:	1mg/ml
Swiss-Prot No.:	Q8NGV0
Other Names:	Olfactory receptor 2Y1; Olfactory receptor OR5-2
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 OR2Y1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OR2Y1 antibody detects endogenous levels of total OR2Y1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
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).