

OR2D2 Antibody [#G545]

Catalog Number:	G545
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H210
Other Names:	Olfactory receptor 2D2; Olfactory receptor OR11-88; Olfactory receptor 11-610; OR11-610; HB2; OR2D2; OR2D1
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 OR2D2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OR2D2 antibody detects endogenous levels of total OR2D2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:5000
References:	Lane R.P., Proc. Natl. Acad. Sci. U.S.A. 98:7390-7395(2001). Suwa M., Submitted (JUL-2001) to the EMBL/GenBank/DDBJ databases. Buettner J.A., Genomics 53:56-68(1998).