

OR1N1 Antibody [#G532]

Catalog Number:	G532
Concentration:	1mg/ml
Swiss-Prot No.:	Q8NGS0
Other Names:	Olfactory receptor 1N1; Olfactory receptor 1-26; OR1-26; Olfactory receptor 1N3; Olfactory receptor OR9-22; OR1N1; OR1N3
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 OR1N1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OR1N1 antibody detects endogenous levels of total OR1N1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:10000
References:	Suwa M., Submitted (JUL-2001) to the EMBL/GenBank/DDBJ databases. Rouquier S., Nat. Genet. 18:243-250(1998). Fuchs T., Genomics 80:295-302(2002).