

OR2AK2 Antibody [#G434]

Catalog Number:	G434
Concentration:	1mg/ml
Swiss-Prot No.:	Q8NG84
Other Names:	Olfactory receptor 2AK2; Olfactory receptor OR1-47; OR2AK2; OR2AK1P
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 OR2AK2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OR2AK2 antibody detects endogenous levels of total OR2AK2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:5000
References:	Suwa M., Submitted (JUL-2001) to the EMBL/GenBank/DDBJ databases. Gregory S.G., Nature 441:315-321(2006). Malnic B., Proc. Natl. Acad. Sci. U.S.A. 101:2584-2589(2004)