

OR5AK3P Antibody [#G471]

Catalog Number:	G471
Concentration:	1mg/ml
Swiss-Prot No.:	Q8NH89
Other Names:	Putative olfactory receptor 5AK3; OR5AK3
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 OR5AK3P.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OR5AK3P antibody detects endogenous levels of total OR5AK3P protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Suwa M., Submitted (JUL-2001) to the EMBL/GenBank/DDBJ databases.