

OR2AE1 Antibody [#G537]

Catalog Number:	G537
Concentration:	1mg/ml
Swiss-Prot No.:	Q8NHA4
Other Names:	Olfactory receptor 2AE1; OR2AE1; OR2AE2
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 OR2AE1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OR2AE1 antibody detects endogenous levels of total OR2AE1 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. The MGC Project Team, Genome Res. 14:2121-2127(2004).