

OR2A25 Antibody [#G430]

Catalog Number:	G430
Concentration:	1mg/ml
Swiss-Prot No.:	A4D2G3
Other Names:	O2A25; olfactory receptor 2A25; olfactory receptor, family 2, subfamily A, member 25; OR2A24P; OR2A25P; OR2A27
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 OR2A25.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OR2A25 antibody detects endogenous levels of total OR2A25 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
References:	Scherer S.W., Science 300:767-772(2003).