

## OR5111 Antibody [#G449]

<b>Catalog Number:</b>	G449
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9H343
<b>Other Names:</b>	Olfactory receptor 5111; Olfactory receptor OR11-39; Odorant receptor HOR5'beta11; OR5111
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human OR5111.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	OR5111 antibody detects endogenous levels of total OR5111 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IF: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Bulger M., Proc. Natl. Acad. Sci. U.S.A. 97:14560-14565(2000). Malnic B., Proc. Natl. Acad. Sci. U.S.A. 101:2584-2589(2004).