

## OR56B4 Antibody [#G470]

<b>Catalog Number:</b>	G470
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q8NH76
<b>Other Names:</b>	O56B4; olfactory receptor 56B4; olfactory receptor OR11-67; olfactory receptor, family 56, subfamily B, 4; OR11-67
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), 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 OR56B4.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	OR56B4 antibody detects endogenous levels of total OR56B4 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Suwa M., Submitted (JUL-2001) to the EMBL/GenBank/DDBJ databases