

## OR4F4/4F5/4F17 Antibody [#G444]

<b>Catalog Number:</b>	G444
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q8NGA8
<b>Other Names:</b>	Olfactory receptor 4F17; OR4F17; OR4F11P, OR4F18, OR4F19
<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 OR4F4/4F5/4F17.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	OR4F4/4F5/4F17 antibody detects endogenous levels of total OR4F4/4F5/4F17 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000   ELISA: 1:5000
<b>References:</b>	Suwa M., Submitted (JUL-2001) to the EMBL/GenBank/DDBJ databases.