

OR6C70 Antibody [#G481]

Catalog Number:	G481
Concentration:	1mg/ml
Swiss-Prot No.:	A6NIJ9
Other Names:	Olfactory receptor 6C70
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 OR6C70.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OR6C70 antibody detects endogenous levels of total OR6C70 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Scherer S.E., Nature 440:346-351(2006). Mural R.J., Submitted (JUL-2005) to the EMBL/GenBank/DDBJ databases.