

OR10G9 Antibody [#G420]

Catalog Number:	G420
Concentration:	1mg/ml
Swiss-Prot No.:	Q8NGN4
Other Names:	Olfactory receptor 10G9; OR10G10P
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 OR10G9.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OR10G9 antibody detects endogenous levels of total OR10G9 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:40000
References:	Suwa M., Submitted (JUL-2001) to the EMBL/GenBank/DDBJ databases.