

OR51A7 Antibody [#G619]

Catalog Number:	G619
Concentration:	1mg/ml
Swiss-Prot No.:	Q8NH64
Other Names:	Olfactory receptor 51A7
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 OR51A7.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OR51A7 antibody detects endogenous levels of total OR51A7 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:10000
References:	Suwa M., Submitted (JUL-2001) to the EMBL/GenBank/DDBJ databases.