

EDG8 Antibody [#G089]

Catalog Number:	G089
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H228
Other Names:	Sphingosine 1-phosphate receptor Edg-8; Endothelial differentiation sphingolipid G-protein-coupled receptor 8; Sphingosine 1-phosphate receptor 5; S1P5; EDG8
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 EDG8.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	EDG8 antibody detects endogenous levels of total EDG8 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	Kothapalli R., <i>Biochim. Biophys. Acta</i> 1579:117-123(2002). Ota T., <i>Nat. Genet.</i> 36:40-45(2004). Im D.-S., <i>Biochemistry</i> 40:14053-14060(2001).