

EDG3 Antibody [#G235]

Catalog Number:	G235
Concentration:	1mg/ml
Swiss-Prot No.:	Q99500
Other Names:	Sphingosine 1-phosphate receptor 3; S1P receptor 3; S1P3; Sphingosine 1-phosphate receptor Edg-3; S1P receptor Edg-3; Endothelial differentiation G-protein coupled receptor 3; S1PR3; EDG3
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 EDG3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	EDG3 antibody detects endogenous levels of total EDG3 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Yamaguchi F., Biochem. Biophys. Res. Commun. 227:608-614(1996). An S., FEBS Lett. 417:279-282(1997). Humphray S.J., Nature 429:369-374(2004).