

SPR1 Antibody [#C11103]

Catalog Number:	C11103
Concentration:	1mg/ml
Swiss-Prot No.:	Q15743
Other Names:	Sphingosylphosphorylcholine receptor, Ovarian cancer G-protein coupled receptor 1, OGR-1, G-protein coupled receptor 68, GPR12A, OGR1
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 SPR1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SPR1 antibody detects endogenous levels of total SPR1 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	An S., FEBS Lett. 375:121-124(1995). Xu Y., Genomics 35:397-402(1996). King M.M., Submitted (APR-2004) to the EMBL/GenBank/DDBJ databases.