

TNFSF9 Antibody [#C10600]

Catalog Number:	C10600
Concentration:	1mg/ml
Swiss-Prot No.:	P41273
Other Names:	Tumor necrosis factor ligand superfamily member 9, 4-1BB ligand, 4-1BBL, TNFSF9
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 TNFSF9.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TNFSF9 antibody detects endogenous levels of total TNFSF9 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
References:	Alderson M.R., Eur. J. Immunol. 24:2219-2227(1994). The MGC Project Team, Genome Res. 14:2121-2127(2004).