

TNFSF15 Antibody [#C10215]

Catalog Number:	C10215
Concentration:	1mg/ml
Swiss-Prot No.:	O95150
Other Names:	Tumor necrosis factor ligand superfamily member 15, Vascular endothelial cell growth inhibitor, TNF ligand-related molecule 1, TNFSF15, TL1, VEGI
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 C-terminal of human TNFSF15.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TNFSF15 antibody detects endogenous levels of total TNFSF15 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
References:	Zhai Y., FASEB J. 13:181-189(1999).