

TNAP1 Antibody [#C11006]

Catalog Number:	C11006
Concentration:	1mg/ml
Swiss-Prot No.:	Q13829
Other Names:	B12, B61, EDP1, MGC2317, TNAP1, tumor necrosis factor, alpha-induced protein 1 (endothelial)
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 TNAP1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TNAP1 antibody detects endogenous levels of total TNAP1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Wolf F.W., J. Biol. Chem. 267:1317-1326(1992). The MGC Project Team, Genome Res. 14:2121-2127(2004). Nousiainen M., Proc. Natl. Acad. Sci. U.S.A. 103:5391-5396(2006).