

## TNAP3 Antibody [#C10444]

<b>Catalog Number:</b>	C10444
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P21580
<b>Other Names:</b>	A20, DNA binding protein A20, EC 3.-.-., TNAP3, TNFA1P2, tumor necrosis factor, alpha-induced protein 3, zinc finger protein A20
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human TNAP3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	TNAP3 antibody detects endogenous levels of total TNAP3 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Opipari A.W. Jr., J. Biol. Chem. 265:14705-14708(1990). Mungall A.J., Nature 425:805-811(2003). Eliopoulos A.G., J. Virol. 73:1023-1035(1999).