

## TNF Receptor II Antibody [#C10436]

<b>Catalog Number:</b>	C10436
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P20333
<b>Other Names:</b>	Tumor necrosis factor receptor superfamily member 1B [Precursor], Tumor necrosis factor receptor 2, TNF-R2, Tumor necrosis factor receptor type II, p75, p80 TNF-alpha receptor, CD120b antigen, Etanercept, Tumor necrosis factor receptor superfamily member 1b, membrane form, Tumor necrosis factor-binding protein 2, TNFRSF1B, TNFBR, TNFR2
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from C-terminal of human TNF Receptor II.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	TNF Receptor II antibody detects endogenous levels of total TNF Receptor II protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Smith C.A., Science 248:1019-1023(1990). Kohno T., Proc. Natl. Acad. Sci. U.S.A. 87:8331-8335(1990). Beltinger C.P., Genomics 35:94-100(1996).