

LTBR Antibody [#C10577]

Catalog Number:	C10577
Concentration:	1mg/ml
Swiss-Prot No.:	P36941
Other Names:	Tumor necrosis factor receptor superfamily member 3 [Precursor], Lymphotoxin-beta receptor, Tumor necrosis factor receptor 2-related protein, Tumor necrosis factor C receptor, LTBR, D12S370, TNFCR, TNFRSF3
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 LTBR.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	LTBR antibody detects endogenous levels of total LTBR protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:40000
References:	Baens M., Genomics 16:214-218(1993). Crowe P.D., Science 264:707-710(1994). Wu M.-Y., J. Biol. Chem. 274:11868-11873(1999).