

TR10A Antibody [#C10014]

Catalog Number:	C10014
Concentration:	1mg/ml
Swiss-Prot No.:	O00220
Other Names:	Tumor necrosis factor receptor superfamily member 10A [Precursor], Death receptor 4, TNF-related apoptosis-inducing ligand receptor 1, TRAIL receptor 1, TRAIL-R1, CD261 antigen, TNFRSF10A, APO2, DR4, TRAILR1
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 TR10A.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TR10A antibody detects endogenous levels of total TR10A protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Pan G., Science 276:111-113(1997). The MGC Project Team, Genome Res. 14:2121-2127(2004). Chaudhary P.M., Immunity 7:821-830(1997).