

CD3 ζ (Ab-42) Antibody [#B0468]

Catalog Number:	B0468
Concentration:	1mg/ml
Swiss-Prot No.:	P20963
Other Names:	CD3-zeta, T-cell receptor T3 zeta chain, T-cell surface glycoprotein CD3 zeta chain precursor, T3Z, TCRZ
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 non-phosphopeptide derived from human CD3 ζ around the phosphorylation site of tyrosine 42 (G-L-Y ^P -Q-G).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CD3 ζ (Ab-42) antibody detects endogenous levels of total CD3 ζ protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Linda A. Trimble, Blood, Aug 2000; 96: 1021 - 1029. Sandeep Krishnan, J. Immunol., Sep 2005; 175: 3417 - 3423. Linda A. Trimble, J. Virol., Aug 2000; 74: 7320 - 7330. Hisashi Arase, J. Immunol., Jan 2001; 166: 21.