

LAT (Ab-191) Antibody [#B0504]

Catalog Number:	B0504
Concentration:	1mg/ml
Swiss-Prot No.:	O43561
Other Names:	36 kDa phospho-tyrosine adaptor protein, Linker for activation of T cells, p36-38, pp36
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 LAT around the phosphorylation site of tyrosine 191 (R-E-Y ^P -V-N).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	LAT (Ab-191) antibody detects endogenous levels of total LAT protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	M Nicosia, J. Virol., Dec 1993; 67: 7276 - 7283. Marta Pineda, J. Biol. Chem., Jul 1999; 274: 19738. Shin-ichiroh Saitoh, J. Exp. Med., Sep 2003; 198: 831. DA Garber, J. Virol., Aug 1997; 71: 5885 - 5893.