

TAL-1 (Ab-122) Antibody [#B8431]

Catalog Number:	B8431
Concentration:	1mg/ml
Swiss-Prot No.:	P17542
Other Names:	T-cell acute lymphocytic leukemia protein 1; SCL, TCL5
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 TAL-1 around the phosphorylation site of serine 122 (Q-L-S ^P -P-P).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TAL-1 (Ab-122) antibody detects endogenous levels of total TAL-1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Aplan P.D., Mol. Cell. Biol. 10:6426-6435(1990). Gregory S.G., Nature 441:315-321(2006). Chen Q., J. Exp. Med. 172:1403-1408(1990).