

MLTK Antibody [#C11781]

Catalog Number:	C11781
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NYL2
Other Names:	AZK, EC 2.7.11.25, MLTK, MRK-beta, mixed lineage kinase-related kinase, sterile alpha motif and leucine zipper containing kinase AZK
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 MLTK.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MLTK antibody detects endogenous levels of total MLTK protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
References:	Gotoh I., J. Biol. Chem. 276:4276-4286(2001). Bloem L.J., J. Mol. Cell. Cardiol. 33:1739-1750(2001). Gross E.A., J. Biol. Chem. 277:13873-13882(2002).