

MYLIP Antibody [#C15558]

Catalog Number:	C15558
Concentration:	1mg/ml
Swiss-Prot No.:	Q8WY64
Other Names:	E3 ubiquitin-protein ligase MYLIP; EC 6.3.2.-; Myosin regulatory light chain-interacting protein; MIR; MYLIP; BZF1; BM-023, PP5242
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 internal of human MYLIP.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MYLIP antibody detects endogenous levels of total MYLIP protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
References:	Olsson P.-A., J. Biol. Chem. 274:36288-36292(1999). Shi W., Submitted (MAY-1997) to the EMBL/GenBank/DDBJ databases. Zhang Q.-H., Genome Res. 10:1546-1560(2000).