

MASTL Antibody [#C11491]

Catalog Number:	C11491
Concentration:	1mg/ml
Swiss-Prot No.:	Q96GX5
Other Names:	Microtubule-associated serine/threonine-protein kinase-like, EC 2.7.11.1, THC2
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 MASTL.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MASTL antibody detects endogenous levels of total MASTL protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Ota T., Nat. Genet. 36:40-45(2004). Deloukas P., Nature 429:375-381(2004). Beausoleil S.A., Nat. Biotechnol. 24:1285-1292(2006).