

MKP-1/2 (Ab-296/318) Antibody [#B8372]

Catalog Number:	B8372
Concentration:	1mg/ml
Swiss-Prot No.:	P28562/Q13115
Other Names:	DUSP1; CL100; MKP1; PTPN10; VH1; MAP kinase phosphatase 1
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 MKP-1/2 around the phosphorylation site of serine 296/318 (I-I-S ^P -P-N).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MKP-1/2 (Ab-296/318) antibody detects endogenous levels of total MKP-1/2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
References:	Keyes S.M., Nature 359:644-647(1992). Kwak S.P., J. Biol. Chem. 269:3596-3604(1994). The MGC Project Team; Genome Res. 14:2121-2127(2004).