

MNK1 (Ab-255) Antibody [#B8143]

Catalog Number:	B8143
Concentration:	1mg/ml
Swiss-Prot No.:	Q9BUB5
Other Names:	EC 2.7.11.1; kinase Mnk1; Map kinase interacting kinase; MAP kinase signal-integrating kinase 1; MAP kinase-interacting serine/threonine kinase 1; MKNK1
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 MNK1 around the phosphorylation site of threonine 255 (L-T-T ^P -P-C).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MNK1 (Ab-255) antibody detects endogenous levels of total MNK1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Fukunaga R., EMBO J. 16:1921-1933(1997). Knauf U., Mol. Cell. Biol. 21:5500-5511(2001). O'Loghlen A., Exp. Cell Res. 299-343-355(2004).