

## MAPK10 Antibody [#C10722]

<b>Catalog Number:</b>	C10722
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P53779
<b>Other Names:</b>	EC 2.7.11.24, JNK3-alpha-2, JNK3A, MAP kinase p49 3F12, MAPK10, MK10, Mitogen-activated protein kinase 10, PRKM10, SAPK-beta, SERK2, Stress-activated protein kinase JNK3, c-Jun N-terminal kinase 3, kinase JNK3, p54- beta
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human MAPK10.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	MAPK10 antibody detects endogenous levels of total MAPK10 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Mohit A.A., Neuron 14:67-78(1995). Lisnock J., Biochemistry 39:3141-3148(2000). Jagadish N., Biochem. J. 389:73-82(2005).