

## JNKK Antibody [#C10624]

<b>Catalog Number:</b>	C10624
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P45985
<b>Other Names:</b>	C-JUN N-terminal kinase kinase 1, Dual specificity mitogen-activated protein kinase kinase 4, EC 2.7.12.2, JNK activating kinase 1, JNK kinase 1, JNKK, JNKK 1, JNKK1, MAP kinase kinase 4, MAP2K4, MAPK/ERK kinase 4, MAPKK 4, MKK4, MP2K4, PRKMK4, SAPK/ERK kinase 1, SEK1, SERK1, c-Jun N- terminal kinase kinase 1, kinase SEK1
<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 JNKK.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	JNKK antibody detects endogenous levels of total JNKK protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Lin A., Science 268:286-290(1995). Derijard B., Science 267:682-685(1995). Su G.H., Cancer Res. 58:2339-2342(1998).