

## SAPK/JNK (Ab-183) Antibody [#B7128]

<b>Catalog Number:</b>	B7128
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P45983/P45984/P53779
<b>Other Names:</b>	c-Jun N-terminal kinase 1; EC 2.7.11.24; JNK-46; JNK1-alpha-2; kinase JNK1; MAPK8; Mitogen-activated protein kinase 8; MK08; p54 gamma; PRKM8; SAPK gamma; Stress-activated protein kinase JNK1
<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 non-phosphopeptide derived from human SAPK/JNK around the phosphorylation site of threonine 183 (M-M-T <sup>P</sup> -P-Y).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	SAPK/JNK (Ab-183) antibody detects endogenous levels of total SAPK/JNK protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Derijard B., Cell 76:1025-1037(1994). Gupta S., EMBO J. 15:2760-2770(1996). Derijard B., Science 267:682-685(1995).