JNK1/2/3 (Phospho-Thr183+Tyr185) Antibody [#8A0500]

Catalog Number:	8A0500
Concentration:	1mg/ml
Origin of Antibody:	Rabbit
Swiss-Prot No.:	P45983/P45984/P53779
Other Names:	EC 2.7.11.24, JNK-46, JNK1-alpha-2, MAPK8, MK08, Mitogen-activated protein kinase 8, PRKM8, SAPK gamma, Stress-activated protein kinase JNK1, c-Jun N-terminal kinase 1, kinase JNK1, p54 gamma
All Sites:	Human: Thr183/Tyr185; Mouse: Thr183/Tyr185; Rat: Thr183/Tyr185
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 phosphopeptide derived from human JNK1/2/3 around the phosphorylation site of threonine183/tyrosine 185 (M-M- $T^{P}-P-Y^{P}-V-V$).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope- specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.
Specificity:	JNK1/2/3 (Phospho-Thr183+Tyr185) antibody detects endogenous levels of JNK1/2/3 only when phosphorylated at threonine183/tyrosine 185.
Reactivity: Applications:	Human, Mouse, Rat WB: 1:500~1:1000 ELISA: 1:10000
References:	Robert A. Kaiser, J. Biol. Chem., Sep 2005; 280: 32602 - 32608. Eleanor T. Coffey, J. Neurosci., Jun 2002; 22: 4335. Nanyue Chen, Cancer Res., Mar 2002; 62: 1300 - 1304. Nan-Shan Chang, J. Biol. Chem., Mar 2003; 278: 9195 - 9202.