

IKK- α / β (Ab-180/181) Antibody [#B0440]

Catalog Number:	B0440
Concentration:	1mg/ml
Swiss-Prot No.:	O15111/O14920
Other Names:	CHUK, Conserved helix-loop-helix ubiquitous kinase, I kappa-B kinase alpha, I-kappa-B kinase 1, IKK-A, IKK-alpha, IKK1, IKKA, IkbKA, IkappaB kinase, Inhibitor of nuclear factor kappa-B kinase alpha subunit, NFKBIKA, Nuclear factor NFkappaB inhibitor kinase alpha, kinase IKK-alpha, EC 2.7.11.10, I-kappa-B kinase 2, I-kappa-B-kinase beta, IKK-B, IKK-beta, IKK2, IKKB, IkbKB, NFKBIKB, Nuclear factor NF-kappa-B inhibitor kinase beta, kinase IKK-beta
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 IKK- α /IKK- β around the phosphorylation site of serine 180/181 (C-T-S ^P -F-V).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	IKK- α / β (Ab-180/181) antibody detects endogenous levels of total IKK- α / IKK- β protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
References:	Bing-Chang Chen, J. Immunol., Oct 2004; 173: 5219 - 5228. Jennifer Harris, J. Immunol., Aug 2006; 177: 2527 - 2535. Chien-Huang Lin, J. Immunol., Sep 2006; 177: 3427 - 3438.