

NF-κB p105/p50 (Ab-927) Antibody [#B0021]

Catalog Number:	B0021
Concentration:	1mg/ml
Swiss-Prot No.:	P19838
Other Names:	DNA-binding factor KBF1, EBP-1, NF-kappa-B1 p84/NF-kappa-B1 p98, NFKB1, NFκB-p50, Nuclear factor NF-kappa-B p105 subunit, nuclear factor NF-kappa-B p50 subunit
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 NF-κB p105/p50 around the phosphorylation site of serine 927 (C-D-S ^P -G-V).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NF-κB p105/p50 (Ab-927) antibody detects endogenous levels of total NF-κB p105/p50 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	John J. Haddad, J. Pharmacol. Exp. Ther., Feb 2002; 300: 567. Sabine Kirchhoff, Eur. J. Biochem., Apr 1999; 261: 546. Gerald Thiel, Eur. J. Biochem., Jul 2004; 271: 2855 – 2862.