

NF- κ B p65 (Ab-310) Antibody [#D0038]

Catalog Number:	D0038
Concentration:	1mg/ml
Swiss-Prot No.:	Q04206
Other Names:	NFKB3; nuclear factor NF-kappa-B p65 subunit; p65; RELA; TF65; transcription factor p65
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-acetylated peptide derived from human NF- κ B p65 around the acetylation site of lysine 310 (T-F-K ^A -S-I).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NF- κ B p65 (Ab-310) antibody detects endogenous levels of total NF- κ B p65 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Ruben S.M., Science 251:1490-1493(1991). Deloukas P., Hum. Mol. Genet. 2:1895-1900(1993). Lyle R., Gene 138:265-266(1994).