

## NF- $\kappa$ B p65 (Ab-281) Antibody [#B8049]

<b>Catalog Number:</b>	B8049
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q04206
<b>Other Names:</b>	NFKB3; nuclear factor NF-kappa-B p65 subunit; p65; RELA; TF65; transcription factor p65
<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 NF- $\kappa$ B around the phosphorylation site of serine 281 (E-L-S <sup>P</sup> -E-P).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	NF- $\kappa$ B p65 (Ab-281) antibody detects endogenous levels of total NF- $\kappa$ B protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
<b>References:</b>	Ruben S.M., Science 251:1490-1493(1991). Deloukas P., Hum. Mol. Genet. 2:1895-1900(1993). Lyle R., Gene 138:265-266(1994).