

NF- κ B p65 (Ab-529) Antibody [#B7162]

- Catalog Number:** B7162-1, B7162-2, B7162
- Amount:** 50 μ g/50 μ l, 100 μ g/100 μ l, 200 μ g/200 μ l
- Swiss-Prot No. :** Q04206
- All Names:** NF- κ B p65, NFKB3, Nuclear factor NF-kappa-B p65 subunit, RELA, TF65, Transcription factor p65, p65
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized non-phosphopeptide derived from human NF- κ B p65 around the phosphorylation site of serine 529.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** NF- κ B p65 (Ab-529) antibody detects endogenous levels of total NF- κ B p65 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:10000
- References:** Xu C, et al (2005) Oncogene:24(28): 4486-95.
McNulty SE, et al. (2004) Pigment Cell Res Apr; 17(2): 173-80.
Madrid LV, et al. (2001) J Biol Chem: 276(22): 18934-40.
Wang D, et al. (2000) J Biol Chem : 275(42): 32592-7.