

## NF-κB p65 (Ab-254) Antibody [#B7172]

- Catalog Number:** B7172-1, B7172-2, B7172
- Amount:** 50μg/50μl, 100μg/100μl, 200μg/200μl
- Form of Antibody:** 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.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized non-phosphopeptide derived from human NF-κB p65 around the phosphorylation site of threonine 254.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** NF-κB p65 (Ab-254) antibody detects endogenous levels of total NF-κB p65 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000    IHC: 1:50~1:100  
ELISA: 1:20000
- Swiss-Prot No. :** Q04206
- References:** Yeh PY, et al. (2004) J Biol Chem. 279(25): 26143-26148.  
Ryo A, et al. (2003) Mol Cell.12 (6): 1413-1426.  
Baeuerle P A, et al. (1994) Annu Rev Immunol. 12:141-179.  
Baeuerle P A, et al. (1996) Cell 87:13-20.  
Haskill S, et al. (1991) Cell 65:1281-1289.