

## NF- $\kappa$ B p105/p50 (Ab-893) Antibody [#B7167]

- Catalog Number:** B7167-1, B7167-2, B7167
- Amount:** 50 $\mu$ g/50 $\mu$ l, 100 $\mu$ g/100 $\mu$ l, 200 $\mu$ g/200 $\mu$ l
- 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- $\kappa$ B p105/p50 around the phosphorylation site of serine 893.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** NF- $\kappa$ B p105/p50 (Ab-893) antibody detects endogenous levels of NF- $\kappa$ B p105/p50 protein around serine 893.
- Reactivity:** Human
- Applications:** WB: 1:500~1:1000    IHC: 1:50~1:100  
ELISA: 1:20000
- Swiss-Prot No. :** P19838
- References:** Hou S, et al. (2003) J Biol Chem. 278(46): 45994-45998.  
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.