

NF- κ B p65 (Ab-468) Antibody [#B7170]

- Catalog Number:** B7170-1, B7170-2, B7170
- Amount:** 50 μ g/50 μ l, 100 μ g/100 μ l, 200 μ g/200 μ l
- 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:** This antiserum was produced against synthesized non-phosphopeptide derived from human NF- κ B p65 around the phosphorylation site of serine 468.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** NF- κ B p65 (Ab-468) 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: 1:20000
- Swiss-Prot No. :** Q04206
- References:** Buss H, et al. (2004) J Biol Chem.
Schmitz ML, et al. (2004) Blood.