

c-Rel(Ab-503) Antibody [#B7211]

- Catalog Number:** B7211-1, B7211-2, B7211
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- 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 c-Rel around the phosphorylation site of Serine 503.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** c-Rel(Ab-503) Antibody detects endogenous levels of c-Rel protein around Serine 503.
- Reactivity:** Human
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
- Swiss-Prot No. :** Q04864
- References:** Martin A G, et al. (2000) J Biol Chem. 275(32): 24383-24391.