



## c-Jun(Ab-63) Antibody

Catalog Number 21001

(For Research Use Only)

- Catalog Number:** 21001-1, 21001-2, 21001  
**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 c-Jun around the phosphorylation site of Serine 63.  
**Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.  
**Specificity/Sensitivity:** c-Jun( Ab-63) Antibody detects endogenous levels of c-Jun protein around Serine 63.  
**Reactivity:** Human, Mouse, Rat  
**Applications:** WB: 1:500~1:1000      ELISA: 1:4000  
**Swiss-Prot No. :** P05412  
**References:** Wei W, et al. (2005). *Cancer Cell*. 8(1): 25-33.  
Sevilla A, et al. (2004). *Oncogene*.23(55): 8950-8958.  
Li L, et al. (2004). *J Biol Chem*.279(6): 4058-4065.  
Hurd C, et al. (2002). *Oncogene*.21(14): 2154-2160.

