

c-Jun(Ab-63) Antibody [#B7045]

Catalog Number: B7045-1, B7045-2, B7045

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-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.