

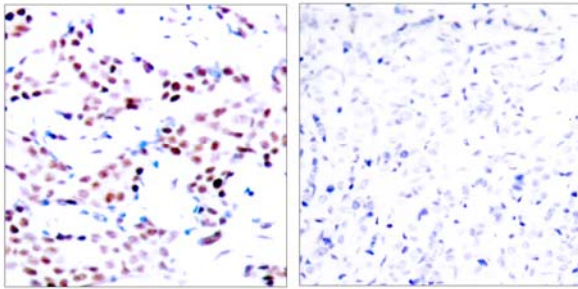


c-Jun(Phospho-Thr91) Antibody

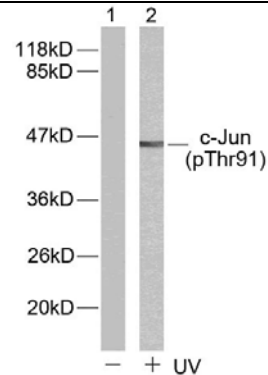
Catalog Number 11021

(For Research Use Only)

- Catalog Number:** 11021-1, 11021-2, 11021
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 phosphopeptide derived from human c-Jun around the phosphorylation site of Threonine 91.
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.
Specificity/Sensitivity: c-Jun(Phospho-Thr91) Antibody detects endogenous levels of c-Jun only when phosphorylated at Threonine 91.
Reactivity: Human, Mouse, Rat
Applications: WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:20000
Swiss-Prot No. : P05412
References: Binetruy B, et al. (1991) Nature. 351: 122-127.
Smeal T, et al. (1991) Nature. 354:494-496.
Derijard B, et al. (1994) Cell. 76:1025-1037.
Kyriakis J M, et al. (1994) Nature. 369: 156-160.



Immunohistochemical analysis of paraffin-embedded breast carcinoma. Left: Using c-Jun (Phospho-Thr91) Antibody(#11021); Right: The same antibody preincubated with synthesized phosphopeptide.



Western blot analysis of extracts using c-Jun (Phospho-Thr91) Antibody(#11021). Line1: The extracts from HeLa cells untreated; Line2: The extracts from HeLa cells treated with UV.