

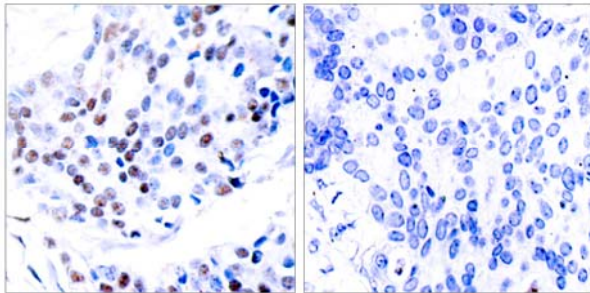


c-Jun (Ab-239) Antibody

Catalog Number 21024

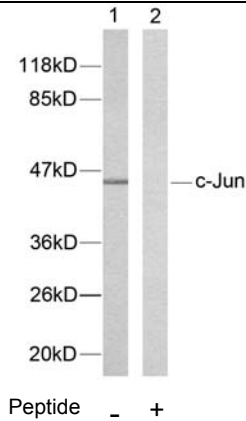
(For Research Use Only)

- Catalog Number:** 21024-1, 21024-2, 21024
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 threonine 239.
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity: c-Jun (Ab-239) antibody detects endogenous levels of total c-Jun protein.
Reactivity: Human, Mouse, Rat
Applications: WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:20000
Swiss-Prot No. : P05412
References: Boyle W J, et al. (1991) Cell. 64(3): 573-584.
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.



Peptide - +

Immunohistochemical analysis of paraffin-embedded human breast carcinoma tissue using c-Jun (Ab-239) antibody (#21024).



Western blot analysis of extracts from HeLa cells using c-Jun (Ab-239) antibody (#21024).