

## Recombinant c-Jun protein

**Catalog No:** 31116

**Expressed In:** Sf9 cells

**Quantity:** 5 µg

**Concentration:** 0.2 µg/µl

**Source:** Human

**Buffer Contents:** 5 µg of Recombinant c-Jun protein in Dilution Buffer AM1 (20 mM Tris-Cl (pH 8), 20% glycerol, 100 mM KCl, 1 mM DTT and 0.2 mM EDTA).

**Background:** c-Jun is a component of the heteromeric **AP-1 transcription factor complex**. AP-1 proteins play a role in the expression of many genes involved in the regulation of cellular processes such as differentiation, proliferation and apoptosis. The transcription factor AP-1 is composed of a mixture of heterodimeric protein complexes derived from the Fos and Jun families, including **c-Fos, FosB, Fra-1, c-Jun, JunB and JunD**. c-Jun can be activated by phosphorylation from the JNK pathway.

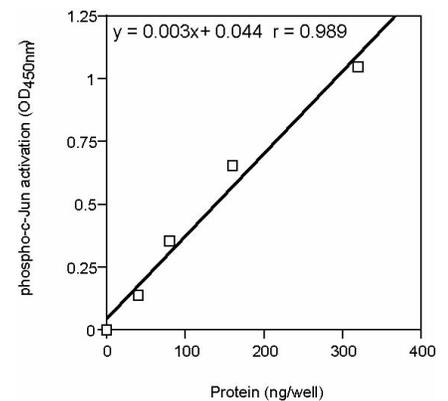
**Protein Details:** Recombinant c-Jun was expressed from full length (accession number NM 002228) with an N-terminal His-Tag in a baculovirus system and purified by an affinity column in combination with FPLC chromatography. The purified recombinant protein is greater than 95% homogeneous and contains no detectable protease, DNase and RNase activity.

**Application Notes:** Recombinant c-Jun is suitable for DNA and protein-protein interaction assays. 100 ng is sufficient for DNA-protein and protein-protein interaction studies. The molecular weight of the protein is ~43 kDa. The standard curve for TransAM® AP-1 c-Jun was generated using the range of 320-40 ng of protein.

NOTE: The presence of Poly [d(I-C)] in buffers may affect protein functionality and should be avoided.

**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



TransAM® standard curve generated using Recombinant c-Jun protein. The standard curve for TransAM® was generated using a range of 320-40 ng of protein and run on the TransAM® AP-1 c-Jun ELISA Kit.