

---

## C-jun Protein

(Catalog No. E60000)

---

### Description

C-JUN amino acids 1-81, is a non-glycosylated, polypeptide chain. C-JUN is a maltose binding protein (MBP) fusion protein with an amino-terminal polyhistidine tag and purified by proprietary chromatographic techniques. C-JUN is a gene which, in combination with c Fos, forms the AP-1 early response transcription factor. C-JUN is activated by the JNK pathway. C-JUN is the putative transforming gene of avian sarcoma virus 17. C-JUN is a protein which is highly similar to the viral protein, and which interacts directly with specific target DNA sequences to regulate gene expression. The C JUN gene is intronless and is mapped to 1p32-p31, a chromosomal region involved in both translocations and deletions in human malignancies.

### Molecular Weight (kDa)

52 kDa

### Formulation

40 mM Tris, pH 7.5, to a concentration of 0.2-1.0 mg/ml.

### Source

Expressed in E.coli expression system

### Size

5 µg

### Storage Temp & Conditions

Store at 4°C if entire vial will be used within 1-2 weeks. Store frozen at -20°C for longer periods of time. Avoid multiple freeze-thaw cycles.

### Purification Notes

Purity: Greater than 95% as determined by SDS-PAGE.

### Application Notes

Enzyme-linked Immunoabsorbent Assay; Western Blot (Recombinant protein); Antibody Production; Protein Array

### Ordering Information:

Products	Size	Cat. No.
C-jun Protein	5 µg	E60000 -1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*