

c-JUN Protein

Recombinant human protein expressed in E. coli cells

Catalog # J05-55G

Lot # E287-3

Product Description

Recombinant human c-JUN (1-79) was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM_002228](#).

Gene Aliases

JUN; AP1; AP-1

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

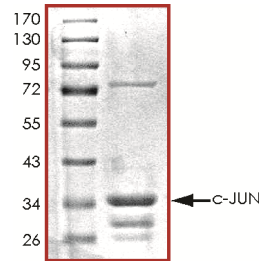
Scientific Background

c-Jun is a proto-oncogene that forms a complex with c-Fos which regulates transcription from promoters containing AP-1 activation elements (1). c-Jun has specific DNA binding activity and following in vitro translation, c-Jun binds as a homodimer to the AP-1 DNA site. The transactivating function of c-Jun is acutely regulated by a wide variety of cellular signals via modulation of phosphorylation of two serines (63 and 73). Jun N-terminal domain kinases (JNKs), are responsible for mediating S63/73 phosphorylation on c-Jun in response to a variety of cellular stimuli including TNF- α , heat stress and u.v. light (2).

References

- Halazonetis, T D. et al: c-Jun dimerizes with itself and with c-Fos, forming complexes of different DNA binding affinities. Cell. 1988 Dec 2;55(5):917-24.
- Dai, T. et al: Stress-activated protein kinases bind directly to the delta domain of c-Jun in resting cells: implications for repression of c-Jun function. Oncogene. 1995 Mar 2;10(5):849-55.

Purity



The purity was determined to be **~70%** by densitometry. Approx. MW **35kDa**.

Upstream Active Kinase

JNK1, Active	Cat # M33-10G
JNK2, Active	Cat # M34-10G
JNK3, Active	Cat # M35-10G

c-JUN Protein

Recombinant human protein expressed in E. coli cells

Catalog Number	J05-55G
Specific Lot Number	E287-3
Purity	~70%
Concentration	0.2 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.