



Recombinant Human Active JNK2 α 1

PRODUCT ANALYSIS SHEET

Catalog Number:	PHO3034	PHO3035	PHO3031
Quantity:	10 μ g	25 μ g	100 μ g

Lot Number:	See product label
Concentration:	See product label
Molecular Weight:	~45 kDa
Purity:	>95% as determined by SDS-PAGE analysis.
Amino Acid Sequence	MSDSKCDSQF YSVQVADSTF TVLKRYQQLK PIGSGAQQIV CAAFDTVLGI NVAVKKLSRP FQNQTHAKRA YRELVLLKCV NHKNIISLLN VFPTQKTL EE FQDVYLV MEL MDANLCQVIH MELDHERMSY LLYQMLCGIK HLHSAGIIHR DLKPSNIVVK SDCTLKILDF GLARTACTNF MMTPTVVTRY YRAPEVILGM GYKENVDIWS VGCIMGELVK GCVIFQGTDH IDQWNKVIEQ LGTPSAEFMK KLQPTVRNYV ENRPKYPGIK FEELFPDWIF PSESERDKIK TSQARDLLSK MLVIDPDKRI SVDEALRHPY ITVWYDPAEA EAPPPQIYDA QLEEREHAIE EWKELIYKEV MDWEERSKNG VVKDQPSAQM QQ
Biological Activity:	~125 units/mg. One unit of JNK2 α 1 activity is equal to 1 nanomole of phosphate transferred to recombinant ATF-2 fusion protein (Catalog # PHF0041) per minute at 30°C with a final ATP concentration of 100 μ M. Recombinant active JNK2 α 1 is capable of autophosphorylation and also phosphorylates c-Jun at serine 63 and serine 73. Kinase activity may vary depending on the substrate and reaction conditions. The optimal concentration should be determined for each specific application. JNK2 α 1 is shorter than JNK2 α 2, and contains a five nucleotide insertion in the C terminal region which causes a shift in the reading frame and subsequent premature termination.
Formulation:	50 mM Tris, pH 7.5, 0.15 M NaCl, 0.27 M sucrose, 10 mM β -mercaptoethanol, 1 mM EGTA, 0.1% Triton X-100, carrier-free.
Sterility:	Filtered through a 0.22 micron sterile filter.
Production:	Recombinant human active JNK2 α 1 (amino acids 1-382) is produced in <i>E. coli</i> and purified via sequential chromatography.
Handling Recommendation:	We recommend that this vial be briefly centrifuged prior to opening to bring the contents to the bottom.
Storage:	Liquid recombinant human active JNK2 α 1 should be kept as a solution in order to maintain full activity. This stock solution should be apportioned into working aliquots and stored at $\leq -80^{\circ}\text{C}$. Keep freeze-thaw cycles to a minimum.
Expiration Date:	Expires one year from date of receipt when stored as instructed.
References:	Sluss, H.K., T. Barrett, B. Derijard, and R.J. Davis (1994) Signal transduction by tumor necrosis factor mediated by JNK protein kinases. <i>Mol. Cell. Biol.</i> 14:8376-8384.

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

www.invitrogen.com

Manufactured under ISO 13485 Quality Standard

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI Hu JNK2 α 1

(Rev 2.0) DCC-08-1232

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.



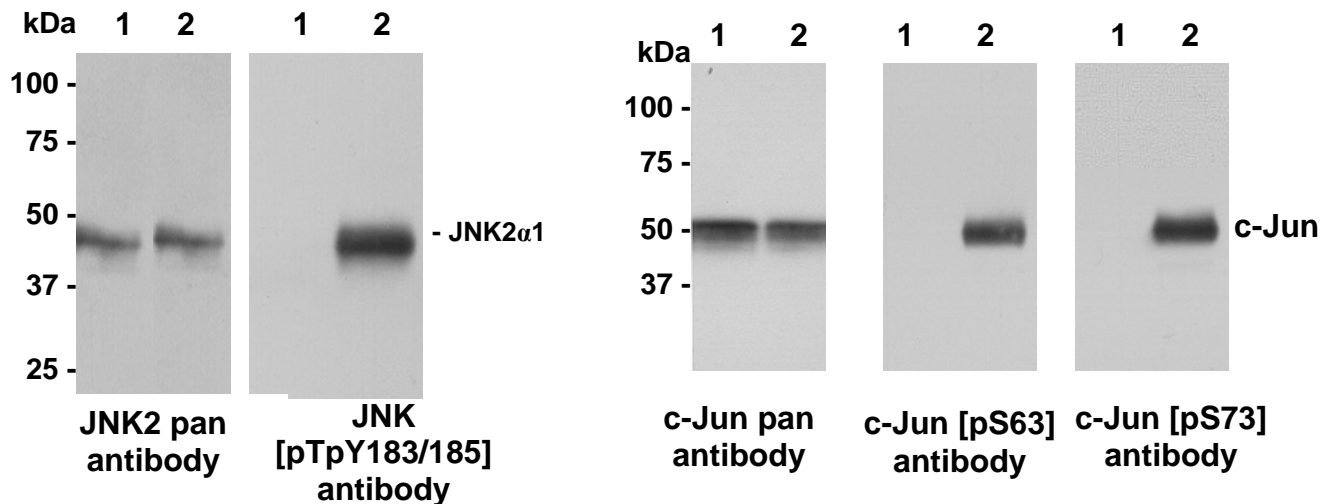
Kallunki, T., B. Su, I. Tsigelny, H.K. Sluss, B. Derijard, G. Moore, R. Davis, and M. Karin (1994) JNK2 contains a specificity-determining region responsible for efficient c-Jun binding and phosphorylation. *Genes Dev.* 8:2996-3007.

**References:
(Continued)**

Gupta, S., T. Barrett, A.J. Whitmarsh, J. Cavanagh, H.K. Sluss, B. Derijard, and R.J. Davis (1996) Selective interaction of JNK protein kinase isoforms with transcription factors. *EMBO J.* 15:2760-2770.

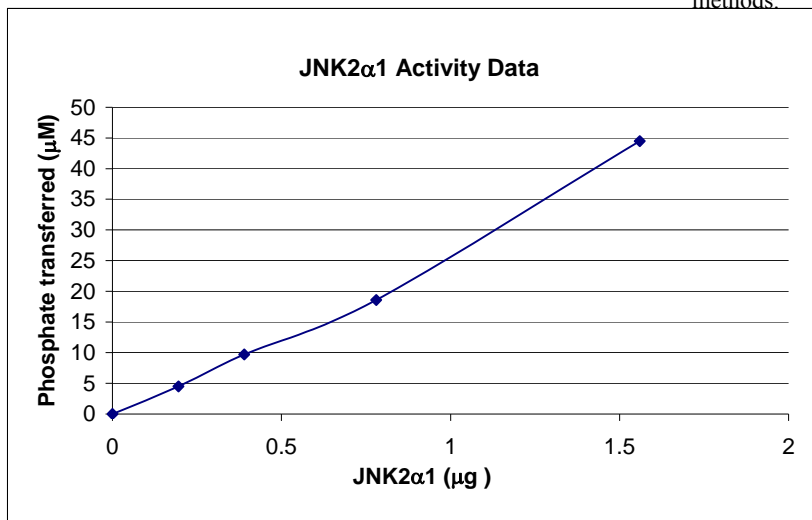
Davis, R.J. (2000) Signal transduction by the JNK group of MAP kinases. *Cell* 103:239-252.

Lin, A. (2003) Activation of the JNK signaling pathway: Breaking the brake on apoptosis. *Bioessays* 25:17-24.



Recombinant active JNK2α1 is phosphorylated at threonine 183/tyrosine 185. Recombinant human inactive (lane 1) and active (lane 2) JNK2α1 were subjected to Western blot analysis using an anti-JNK pan antibody or anti-JNK [pTpY183/185] antibody (Catalog #44-682) in conjunction with chemiluminescence detection methods.

Recombinant active JNK2α1 phosphorylates c-Jun fusion protein. Recombinant c-Jun fusion protein (2 µg; Catalog #PHF0051) was incubated without (lane 1) and with (lane 2) 0.5 µg of recombinant human active JNK2α1 in an *in vitro* kinase reaction. Kinase reactions were analyzed by Western blot analysis using an anti-c-Jun pan antibody, anti-c-Jun [pS63] antibody, or anti-c-Jun [pS73] antibody (Catalog #44-292) in conjunction with chemiluminescence detection methods.



This product is for research use only. Not for use in diagnosis.

www.invitrogen.com

Manufactured under ISO 13485 Quality Standard

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI Hu JNK2α1

(Rev 2.0) DCC-08-1232

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.

invitrogen