

# PRODUCT INFORMATION



## JNK1 $\alpha$ 2 Isoform (human, recombinant)

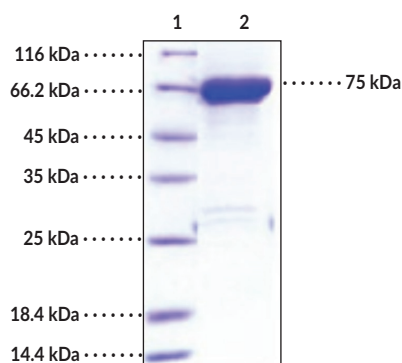
Item No. 43698

### Overview and Properties

<b>Synonyms:</b>	c-Jun N-terminal Kinase 1, JNK-46, MAPK8, MAP Kinase 8, Mitogen-activated Protein Kinase 8, SAPK1C, Stress-activated Protein Kinase 1C
<b>Source:</b>	Recombinant human N-terminal GST-tagged JNK1 $\alpha$ 2 isoform expressed in insect cells
<b>Amino Acids:</b>	1-427 (full length)
<b>Uniprot No.:</b>	P45983
<b>Molecular Weight:</b>	75 kDa
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	$\geq$ 1 year
<b>Purity:</b>	$\geq$ 90% estimated by SDS-PAGE
<b>Supplied in:</b>	Sterile 50 mM Tris, with 100 mM sodium chloride, pH 8.0, and 25% glycerol
<b>Endotoxin Testing:</b>	<1.0 EU/ $\mu$ g, determined by the LAL endotoxin assay
<b>Bioactivity:</b>	See figure for details

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Image



Lane 1: MW Markers  
Lane 2: JNK1  $\alpha$ 2 Isoform

**SDS-PAGE Analysis of JNK1  $\alpha$ 2 Isoform.** This protein has a calculated molecular weight of 75 kDa. It has an apparent molecular weight of approximately 65 kDa by SDS-PAGE under reducing conditions.

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/24/2025

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

JNK1 is a stress-activated serine/threonine protein kinase and member of the MAPK family.<sup>1</sup> It is composed of a short N-terminal helical domain, a flexible connector that contains the active site, and a large helical C-terminal domain.<sup>2</sup> JNK1 is encoded by *MAPK8*, which produces four isoforms via alternative splicing.<sup>3</sup> JNK1 is ubiquitously expressed and is found in the cytoplasm.<sup>1,3,4</sup> JNKs are activated by phosphorylation at threonine 183 (Thr183) and tyrosine 185 (Tyr185) by MAP kinase kinase 7 (MKK7) and MKK4, respectively, in response to stress stimuli, such as cellular stress, radiation, inflammation, or oxidative stress.<sup>1,5</sup> After activation, JNKs localize to the nucleus where they phosphorylate transcription factors, nuclear receptors, and adaptor proteins that broadly regulate cell proliferation, cell survival, and apoptosis.<sup>4</sup> Increased intratumoral levels of JNK1 are associated with decreased survival in patients with hepatocellular carcinoma (HCC).<sup>6</sup> Cayman's JNK1  $\alpha 2$  Isoform (human, recombinant) protein consists of 652 amino acids and has a calculated molecular weight of 75 kDa. This protein migrates at approximately 65 kDa by SDS-PAGE under reducing conditions.

## References

---

1. Kumar, A., Singh, U.K., Kini, S.G., *et al.* JNK pathway signaling: A novel and smarter therapeutic target for various biological diseases. *Future Med. Chem.* **7(15)**, 2065-2086 (2015).
2. Heo, Y.S., Kim, S.K., Seo, C.I., *et al.* Structural basis for the selective inhibition of JNK1 by the scaffolding protein JIP1 and SP600125. *EMBO J.* **23(11)**, 2185-2195 (2004).
3. Zhang, T., Inesta-Vaquero, F., Niepel, M., *et al.* Discovery of potent and selective covalent inhibitors of JNK. *Chem. Biol.* **19(1)**, 140-154 (2012).
4. Bogoyevitch, M.A. and Kobe, B. Uses for JNK: The many and varied substrates of the c-Jun N-terminal kinases. *Microbiol. Mol. Biol. Rev.* **70(4)**, 1061-1095 (2006).
5. Thévenin, A.F., Zony, C.L., Bahnson, B.J., *et al.* Activation by phosphorylation and purification of human c-Jun N-terminal kinase (JNK) isoforms in milligram amounts. *Protein Expr. Purif.* **75(2)**, 138-146 (2011).
6. Chang, Q., Chen, J., Beezhold, K.J., *et al.* JNK1 activation predicts the prognostic outcome of the human hepatocellular carcinoma. *Mol. Cancer* **8**, 64 (2009).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM