

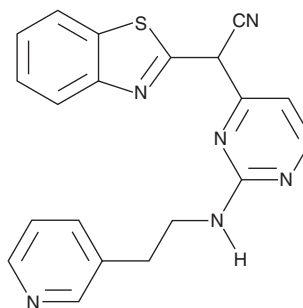
PRODUCT INFORMATION



JNK Inhibitor V

Item No. 17542

CAS Registry No.: 345987-15-7
Formal Name: a-[2-[[2-(3-pyridinyl)ethyl]amino]-4-pyrimidinyl]-2-benzothiazoleacetonitrile
Synonyms: AS-601245,
c-Jun N-terminal Kinase Inhibitor V
MF: C₂₀H₁₆N₆S
FW: 372.5
Purity: ≥98% (sum of isomers)
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

JNK inhibitor V is supplied as a crystalline solid. A stock solution may be made by dissolving the JNK inhibitor V in the solvent of choice, which should be purged with an inert gas. JNK inhibitor V is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of JNK inhibitor V in these solvents is approximately 2 mg/ml.

Description

c-Jun N-terminal kinases (JNKs) phosphorylate c-Jun and regulate transcription in response to an array of inflammatory signals.¹ JNK inhibitor V is an ATP-competitive inhibitor of JNK1, JNK2, and JNK3 (IC₅₀s = 150, 220, and 70 nM, respectively).^{2,3} While initial studies demonstrated 10- to 100-fold selectivity for JNK isoforms over a panel of 25 other kinases, strong interactions of JNK inhibitor V with GSK3B, Pim-1, Pim-3, and other kinases, evaluated as a shift in thermal melting point, suggest additional targets exist.^{2,4} *In vivo* efficacy of JNK inhibitor V against JNK isoforms has been demonstrated in gerbils, mice, and rats via oral, intravenous, or intraperitoneal administration.^{4,5} This compound is commonly used to investigate the role of JNK signaling in cells and animals.^{6,7}

References

1. Johnson, G.L. and Lapadat, R. Mitogen-activated protein kinase pathways mediated by ERK, JNK, and p38 protein kinases. *Science* **298**(5600), 1911-1912 (2002).
2. Gaillard, P., Jeanclaude-Etter, I., Ardisson, V., *et al.* Design and synthesis of the first generation of novel potent, selective, and *in vivo* active (benzothiazol-2-yl)acetonitrile inhibitors of the c-Jun N-terminal kinase. *J. Med. Chem.* **48**(14), 4596-4607 (2005).
3. Carboni, S., Hiver, A., Szyndralewicz, C., *et al.* AS601245 (1,3-benzothiazol-2-yl (2-[[2-(3-pyridinyl) ethyl] amino]-4 pyrimidinyl) acetonitrile): A c-Jun NH₂-terminal protein kinase inhibitor with neuroprotective properties. *J. Pharmacol. Exp. Ther.* **310**(1), 25-32 (2004).
4. Fedorov, O., Marsden, B., Pogacic, V., *et al.* A systematic interaction map of validated kinase inhibitors with Ser/Thr kinases. *Proc. Natl. Acad. Sci. USA* **104**(51), 20523-20528 (2007).
5. Ferrandi, C., Ballerio, R., Gaillard, P., *et al.* Inhibition of c-Jun N-terminal kinase decreases cardiomyocyte apoptosis and infarct size after myocardial ischemia and reperfusion in anaesthetized rats. *Br. J. Pharmacol.* **142**(6), 953-960 (2004).
6. Oktem, O., Buyuk, E., and Oktay, K. Preantral follicle growth is regulated by c-Jun-N-terminal kinase (JNK) pathway. *Reprod.Sci.* **18**(3), 269-276 (2011).
7. Tu, Y.F., Tsai, Y.S., Wang, L.W., *et al.* Overweight worsens apoptosis, neuroinflammation and blood-brain barrier damage after hypoxic ischemia in neonatal brain through JNK hyperactivation. *J. Neuroinflammation* **8**:40, (2011).

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, 03/13/2023

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