# **PRODUCT** INFORMATION



**JNK Inhibitor XVI** 

Item No. 18096

CAS Registry No.:	1410880-22-6
Formal Name:	3-[[4-(dimethylamino)-1-oxo-2-
	buten-1-yl]amino]-N-[3-methyl-4-
	[[4-(3-pyridinyl)-2-pyrimidinyl]amino]
	phenyl]-benzamide
Synonyms:	JNK-IN-8,
	c-Jun N-terminal Kinase Inhibitor XVI
MF:	$C_{20}H_{20}N_7O_2$
FW:	507.6
Purity:	≥95% <sup>0</sup> <sup>1</sup>
UV/Vis.:	$\lambda_{max}$ : 277 nm
Supplied as:	A crystalline solid
Storage:	-20°C
Stability:	≥4 years
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis	

specifications. Batch specific analytical results are provided on each certificate of analysis

# Laboratory Procedures

JNK inhibitor XVI is supplied as a crystalline solid. A stock solution may be made by dissolving the JNK inhibitor XVI in the solvent of choice, which should be purged with an inert gas. JNK inhibitor XVI is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of JNK inhibitor XVI in these solvents is approximately 2, 16, and 20 mg/ml, respectively.

JNK inhibitor XVI is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JNK inhibitor XVI should first be dissolved in DMF and then diluted with the aqueous buffer of choice. JNK inhibitor XVI has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

# Description

c-Jun N-terminal kinases (JNKs) are MAP kinase family members that become highly activated after cells are exposed to stress conditions and are poorly activated by exposure to growth factors or mitogens. JNK inhibitor XVI is a selective, irreversible JNK inhibitor (IC<sub>50</sub>s = 4.67, 18.7, and 0.98 nM for JNK1, 2, and 3, respectively) that prevents phosphorylation of c-Jun in A375 and HeLa cells with  $EC_{50}$  values of 338 and 486 nM, respectively.<sup>1</sup> It has been shown to inhibit JNK kinase activity by a mechanism that depends on covalent modification of cysteine<sup>116</sup> in the ATP-binding motif.<sup>1</sup> This compound has been used to explore the role of JNK in mediating cancer cell death.<sup>2,3</sup>

# References

- 1. 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).
- 2. Li, Q., Song, X., Ji, Y., et al. The dual mTORC1 and mTORC2 inhibitor AZD8055 inhibits head and neck squamous cell carcinoma cell growth in vivo and in vitro. Biochem. Bioph. Res. Commun. 440(4), 701-706 (2013).
- 3. Fallahi-Sichani, M., Moerke, N.J., Niepel, M., et al. Systematic analysis of BRAF<sup>V600E</sup> melanomas reveals a role for JNK/c-Jun pathway in adaptive resistance to drug-induced apoptosis. Mol. Syst. Biol. 11(3), 797 (2015).

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

## SAFFTY 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.

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, 12/01/2022

# 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