

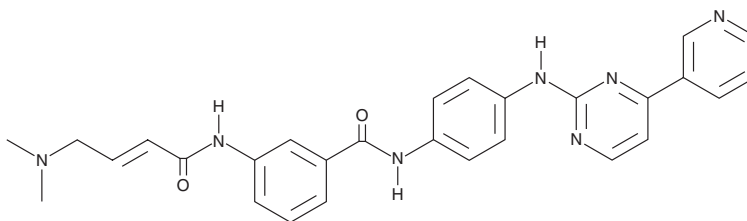
PRODUCT INFORMATION



JNK Inhibitor VII

Item No. 28831

CAS Registry No.: 1408064-71-0
Formal Name: 3-[[4-(dimethylamino)-1-oxo-2-buten-1-yl]amino]-N-[[4-[[4-(3-pyridinyl)-2-pyrimidinyl]amino]phenyl]-benzamide
Synonyms: JNK-IN-7, c-Jun N-terminal Kinase Inhibitor VII
MF: C₂₈H₂₇N₇O₂
FW: 493.6
Purity: ≥98%
UV/Vis.: λ_{max}: 284, 299 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.

Laboratory Procedures

JNK inhibitor VII is supplied as a crystalline solid. A stock solution may be made by dissolving the JNK inhibitor VII in the solvent of choice, which should be purged with an inert gas. JNK inhibitor VII is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of JNK inhibitor VII in these solvents is approximately 10 mg/ml. JNK inhibitor VII is slightly soluble in ethanol.

JNK inhibitor VII is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JNK inhibitor VII should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. JNK inhibitor VII has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

JNK inhibitor VII is a potent and non-selective inhibitor of JNKs (IC₅₀s = 1.54, 1.99, and 0.75 nM for JNK1, -2, and -3, respectively).¹ It inhibits phosphorylation of c-Jun in HeLa and A375 cells (EC₅₀s = 130 and 244 nM, respectively).

Reference

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).

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, 10/06/2022

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