

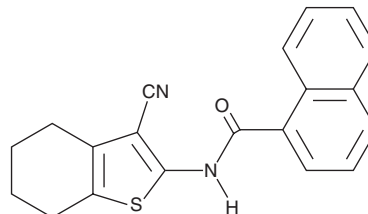
PRODUCT INFORMATION



JNK Inhibitor IX

Item No. 15624

CAS Registry No.: 312917-14-9
Formal Name: N-(3-cyano-4,5,6,7-tetrahydrobenzo[b]thien-2-yl)-1-naphthalenecarboxamide
Synonyms: c-Jun N-terminal Kinase Inhibitor IX, TCS JNK 5a
MF: C₂₀H₁₆N₂OS
FW: 332.4
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 316 nm
Supplied as: A crystalline solid
Storage: Room temperature
Stability: ≥4 years



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

Laboratory Procedures

JNK inhibitor IX is supplied as a crystalline solid. A stock solution may be made by dissolving the JNK inhibitor IX in the solvent of choice, which should be purged with an inert gas. JNK inhibitor IX is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of JNK inhibitor IX in these solvents is approximately 30 mg/ml.

JNK inhibitor IX is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JNK inhibitor IX should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. JNK inhibitor IX has a solubility of approximately 0.25 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 IX is a thienynaphthamide compound that targets the ATP binding site of JNK2 and JNK3, disrupting activity with pIC₅₀ values of 6.5 and 6.7, respectively.¹ It demonstrates little activity against JNK1, p38α, and a panel of more than 30 other kinases (pIC₅₀ < 5.0).¹

Reference

1. Angell, R.M., Atkinson, F.L., Brown, M.J., *et al.* N-(3-Cyano-4,5,6,7-tetrahydro-1-benzothien-2-yl)amides as potent, selective, inhibitors of JNK2 and JNK3. *Bioorg. Med. Chem. Lett.* **17**(5), 1296-1301 (2007).

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, 10/13/2022

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