

N-(3-Cyano-4,5,6,7-tetrahydro-1-benzothien-2-yl)-1-naphthylamide

Thienyl naphthylamide compound. Potent inhibitor of JNK2 ($\text{pIC}_{50}=6.5$) and JNK3 ($\text{pIC}_{50}=6.7$) targeting the ATP-binding site. Has no activity against JNK1.

Citations: 1

[View Online »](#)

Ordering Information

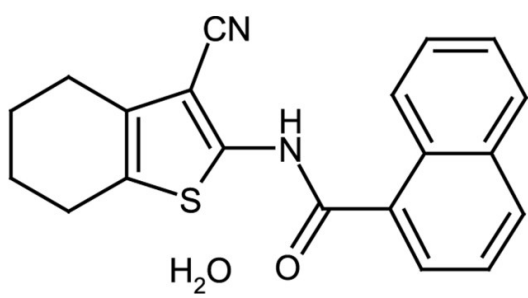
[Order Online »](#)

ALX-270-490-M005

5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 6 months when stored at -20°C.
Handling	Protect from light. Packaged under inert gas. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Tan solid.
CAS	312917-14-9
Couple Target	Jnk
Couple Type	Inhibitor
Formula	$C_{20}H_{16}N_2OS \cdot H_2O$
Identity	Determined by NMR.
MW	332.4 . 18.0
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (10mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)