

AS601245

Jnk inhibitor

Potent and cell permeable ATP competitive JNK inhibitor (hJNK1: IC_{50} =150nM, hJNK2: IC_{50} =220nM and hJNK3: IC_{50} =70 nM). Displays anti-inflammatory properties and has been shown to reduce TNF- α plasma levels induced by LPS in mice.

Citations: 8

[View Online »](#)

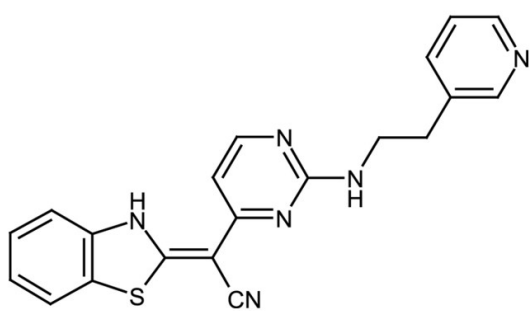
Ordering Information

[Order Online »](#)

ALX-270-443-M001	1mg
ALX-270-443-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 3 months when stored at -20°C.
Handling	Hygroscopic. Packaged under inert gas. Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1,3-Benzothiazol-2-yl-(2-[[2-(3-pyridinyl)ethyl]amino]-4-pyrimidinyl) acetonitrile, JNK Inhibitor V
Appearance	White to yellow solid.
CAS	345987-15-7
Couple Target	Jnk
Couple Type	Inhibitor
Formula	$C_{20}H_{16}N_6S$
Identity	Determined by 1H -NMR.
MW	372.5
Purity	≥95% (HPLC)
Purity Detail	Sum of two isomers.
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-
be@enzolifesciences.com](mailto:info-be@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)