

GSK-J4

JMJ H3K27 Demethylase inhibitor (Cell Permeable)

Histone demethylase JMJD3/UTX inhibitor. Inhibits demethylation of histone H3K27. Reduces LPS-induced proinflammatory cytokine production in primary human macrophages. Rescues newborn pups from embryonic lethality in BRAF knockin mice which recapitulate major features of RASopathies.

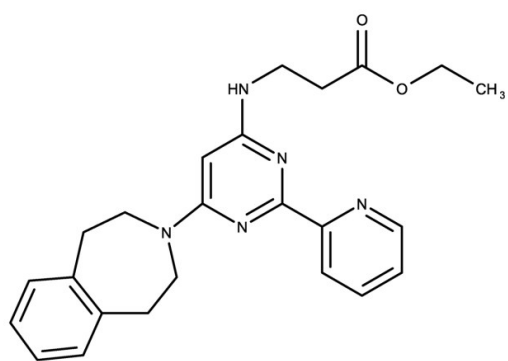
Ordering Information

[Order Online »](#)

ENZ-CHM482-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-[2-(2-Pyridinyl)-6-(1,2,4,5-tetrahydro-3H-3-benzazepin-3-yl)-4-pyrimidinyl]-β-alanine ethyl ester

Appearance Yellow waxy solid.

CAS 1373423-53-0

Couple Target Demethylase

Couple Type Inhibitor

Formula $C_{24}H_{27}N_5O_2$

Identity Determined by NMR.

MW 417.48

Purity ≥98% (TLC)

Solubility Soluble in DMSO (40 mg/ml) or ethanol (40 mg/ml).

Last modified: February 27, 2025



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)