

GSK-J4 (ethylester)

Epigenetic Enzyme Inhibitor

Catalog # J48-904C

Lot # P3320-20

CAS # 1373423-53-0

Product Description

Molecular Formula: C₂₄H₂₇N₅O₂

Appearance: Yellow wax

Molecular Weight: 417.5

Purity: >99%; NMR (Conforms)

Solubilization: May be dissolved in DMSO (40 mg/ml); or ethanol (40 mg/ml)

Alias

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

Specific Activity

JMJ H3K27 Demethylase Inhibitor (Cell Perm)

Storage and Stability

Store, as supplied, at -20°C. Store solutions at -20°C for up to 3 months.

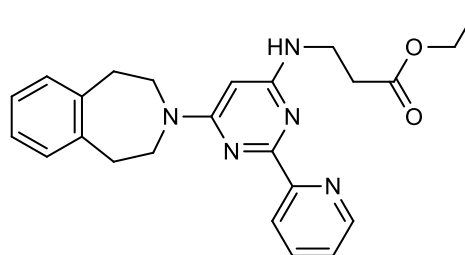
Scientific Background

Histone demethylase JMJD3/UTX inhibitor. Inhibits demethylation of histone H3K27. Reduces LPS-induced proinflammatory cytokine production in primary human macrophages (IC₅₀ = 9 μM for the inhibition of TNFα release). Cell permeable, ethyl ester of GSK-J1¹. Rescues newborn pups from embryonic lethality in BRAF knockin mice which recapitulate major features of RASopathies².

References

1. L Kruidenier et al. Nature 2012 488:404
2. Si Inoue et al. Hum. Mol. Genet. 2014 July 17, Epub

Molecular Structure



GSK-J4 (ethylester)

Epigenetic Enzyme Inhibitor

Catalog #	J48-904C
CAS #	1373423-53-0
Lot #	P3320-20
Purity	>99%
Format	Yellow wax
Solubilization	May be dissolved in DMSO (40 mg/ml); or ethanol (40 mg/ml)
Storage & Shipping	Store, as supplied, at -20°C. Store solutions at -20°C for up to 3 months.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.