## GSK-J4 (cell permeable)



Catalog No: 14070, 14071

Format: 25 mg, 5 mg

**Chemical Properties:** 

MW = 417.5

 $C_{24}H_{27}N_5O_2$  CAS 1373423-53-0

Physical Properties: Yellow wax

Names: N-[2-(2-Pyridinyl)-6-(1,2,4,5-tetrahydro-3H-3-benzazepin-3-yl)-4-

pyrimidinyl]-β-alanine ethyl ester

**Pharmacology:** Histone demethylase JMJD3/UTX inhibitor. Inhibits demethylation of histone H3K27. Reduces LPS-induced proinflammatory cytokine production in primary human macrophages (IC $_{50}$  = 9  $\mu$ M for the inhibition of TNF $\alpha$  release). Cell permeable, ethyl ester of GSK J1 (Catalog No. 14068)(ref1). Rescues newborn pups from embryonic lethality in BRAF knockin mice which recapitulate major features of RASopathies (ref2).

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

Fluorescent Properties: N/A

**Quality Control:** 

>99% (TLC); NMR (Conforms)

References:

1. L Kruidenier et al. Nature 2012, 488:404

2. SI Inoue et al. Hum. Mol. Genet. 2014 July 17, Epub.

**Storage and Guarantee:** Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 3 months. This product is guaranteed for 6 months from date of arrival.