## GSK-J1 (cell impermeable)



Catalog No: 14068, 14069

Format: 25 mg, 5 mg

**Chemical Properties:** 

MW = 389.5 $C_{22}H_{23}N_5O_2$ 

CAS 1373422-53-7

Physical Properties: Pale yellow powder

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

pyrimidinyl]-β-alanine

**Pharmacology:** A potent and selective inhibitor of jumonji H3K27 histone demethylases JMJD3 and UTX (IC $_{50}$  = 60 nM, human JMJD3). This is the first known inhibitor selective for the H3K27me3-specific JMJ subfamily which binds to the active catalytic site of the enzyme. The COOH group confers cell impermeability and as such is useful as a standard *in vitro* assay. A cell permeable ethyl ester analog is also available (Catalog No. 14070).

Solubilization: May be dissolved in DMSO (20 mg/ml, warm)

Fluorescent Properties: N/A

**Quality Control:** 

>99% (TLC); NMR (Conforms)

References:

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

**Storage and Guarantee:** Store desiccated as supplied at room temperature 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.

