

ML-324

Epigenetic Enzyme Inhibitor

Catalog # J48-904E

Lot # P3320-23

CAS # 1222800-79-4

Product Description

Molecular Formula: C₂₁H₂₃N₃O₂

Appearance: Light brown-colored powder

Molecular Weight: 349.43

Purity: >98% (TLC); NMR (Conforms)

Solubilization: Soluble in DMSO (20 mg/ml, warm); or ethanol (6 mg/ml, warm)

Alias

N-(3-(dimethylamino)propyl)-4-(8-hydroxyquinolin-6-yl)benzamide

Specific Activity

JMJD2 Inhibitor

Storage and Stability

Store desiccated as supplied at -20°C for up to 1 year.

Store solutions at -20°C for up to 1 month.

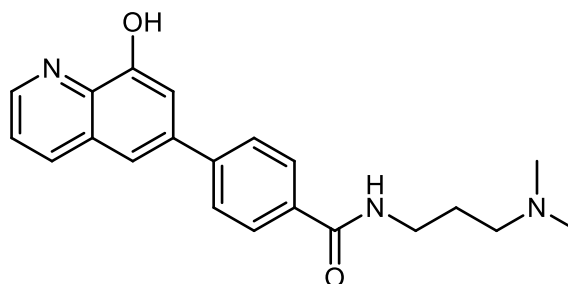
Scientific Background

Inhibits JMJD2 histone demethylases. (JMJD2 IC₅₀=920 nM) Represses the expression of viral (HCMV and HSV-1) immediate early genes (IC₅₀ ~ 10 μM) and abrogates infection. Represses the reactivation of HSV from the latent state in sensory neurons¹. Cell permeable.

References

1. Yu Liang et al. Sci. Transl. Med. 2013 5:167ra5

Molecular Structure



ML-324

Epigenetic Enzyme Inhibitor

Catalog #	J48-904E
CAS #	1222800-79-4
Lot #	P3320-23
Purity	>98% (TLC)
Format	Light brown-colored powder
Solubilization	Soluble in DMSO (20 mg/ml, warm); or ethanol (6 mg/ml, warm)
Storage & Shipping	Store desiccated as supplied at -20°C for up to 1 year. Store solutions at -20°C for up to 1 month.

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.