

ML-324

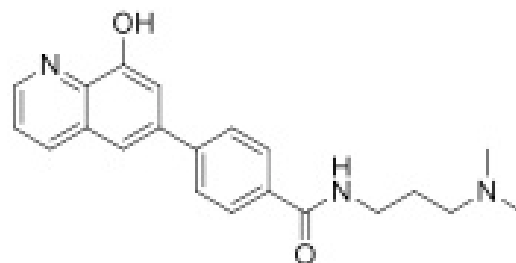
Catalog No: 14078, 14079**Format:** 25 mg, 5 mg**Chemical Properties:**

MW = 349.4

 $C_{21}H_{23}N_3O_2$

CAS Not found

Physical Properties: Light brown powder

Names: ML324, N-(3-(dimethylamino)propyl)-4-(8-hydroxyquinolin-6-yl)benzamide**Pharmacology:** Inhibits JMJD2 histone demethylases (JMJD2 IC_{50} = 920 nM). Represses the expression of viral (HCMV and HSV-1) immediate early genes (IC_{50} ~ 10 M) and abrogates infection. Represses the reactivation of HSV from the latent state in sensory neurons (ref 1). Cell permeable.**Solubilization:** May be dissolved in DMSO (20 mg/ml, warm); Ethanol (6 mg/ml, warm)**Fluorescent Properties:** N/A**Quality Control:**

>98% (TLC); NMR (Conforms)

References:

1. Yu Liang *et al. Sci. Transl. Med.* 2013, 5:167ra5**Storage and Guarantee:** Store desiccated as supplied at -20°C for up to 1 year. Store solutions at -20°C for up to 1 month. This product is guaranteed for 6 months from date of arrival.