

UNC0224

Catalog #: 27303

Size: 5 mg

CAS Registry #: 1197196-48-7

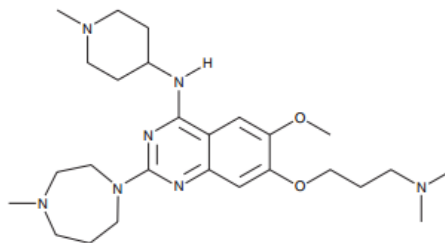
Purity: ≥98%

Chemical Formula: C₂₆H₄₃N₇O₂

Molecular Weight: 485.67

Lot #: 120627

Structure:



Description: UNC0224 is a potent and selective inhibitor of G9a, a histone methyltransferase (HMT). It also inhibits GLP (G9a-Like Protein), a closely-related H3K9 HMT, but is more than 1,000-fold selective against SET7/9 and SET8.

Appearance: A crystalline solid

Solubility: Soluble in DMSO, DMF and ethanol, purged with an inert gas. Solubility in DMSO and DMF is approximately 2 mg/ml, and solubility in ethanol is ~15 mg/ml. Do not store aqueous solutions for more than one day.

Biological Activity: UNC0224 inhibits G9a with an IC₅₀ value of 15 nM. It has been shown to inhibit GLP with assay-dependent IC₅₀ values of 20-58 nM.

Storage/Stability: Store at or below -20°C for up to two years.

Quality Control: The purity was determined by HPLC analysis.

References:

1. Tachibana, M., *et al.*, *Genes Dev.* 2002; **16**:1779-1791.
2. Wagschal, A., *et al.*, *Mol. Cell. Biol.* 2008; **28(3)**:1104-1113.
3. Thomas, L.R., *et al.*, *J. Immunol.* 2008; **181**:485-493.