

BIX01294 (hydrochloride hydrate)

Catalog #: 27220

Lot #: 120626

Size: 10 mg

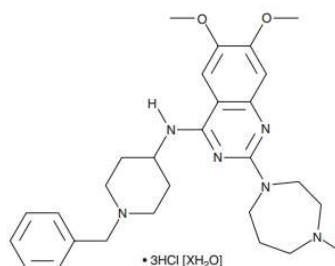
Structure:

CAS Registry #: -

Purity: ≥98%

Chemical Formula: C₂₈H₃₈N₆O₂ · 3HCl [XH₂O]

Molecular Weight: 600



Description: BIX01294 (hydrochloride hydrate) is a selective inhibitor of G9a, a histone methyltransferase (HMT). It can also inhibit G9a-like protein (GLP), but is less effective. There are no known inhibitory effects on other HMTs. Also, BIX01294 has been used in combination with BayK8644 to facilitate generation of induced pluripotent stem cells from somatic *in vitro*.

Appearance: A crystalline solid

Solubility: Dissolve in water or other aqueous buffers. Solubility in water is ~20 mg/ml. Do not store aqueous solutions for more than one day.

Biological Activity: BIX01294 inhibits G9a with an IC₅₀ = 1.7 μM, and GLP with an IC₅₀ = 38 μM.

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.
4. Kubicek, S., *et al.*, *Molecular Cell* 2007; **25**:473-481.
5. Shi, Y., *et al.*, *Cell Stem Cell* 2008; **3(5)**:568-574.