

## BIX-01294 3HCl

Epigenetic Enzyme Inhibitor

**Catalog # S344-903**

Lot # P3320-5

CAS # 935693-62-2

### Product Description

Molecular Formula: C<sub>28</sub>H<sub>38</sub>N<sub>6</sub>O<sub>2</sub>·3HCl

Appearance: Off-white crystalline powder

Molecular Weight: 600.02

Purity: >98%; NMR (Conforms)

Solubilization: May be dissolved in water (50 mg/ml) or DMSO (50 mg/ml)

### Alias

2-(Hexahydro-4-methyl-1H-1,4-diazepin-1-yl)-6,7-dimethoxy-N-[1-(phenylmethyl)-4-piperidiny]-4-quinazolinamine trihydrochloride

### Specific Activity

Histone Methyl Transferase Inhibitor

### Storage and Stability

Store desiccated as supplied at ambient for up to 2 years. Store solutions at -20°C for up to 3 months.

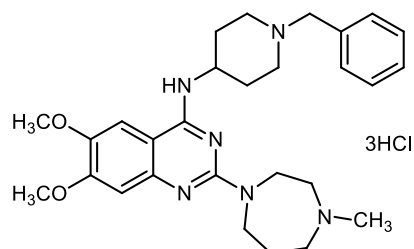
### Scientific Background

BIX-01294 is a selective inhibitor of G9a histone methyltransferase (G9aHMTase; IC<sub>50</sub> = 1.7 μM) as well as GLP HMTase (IC<sub>50</sub> = 38 μM) leading to a decrease in H3K9Me2 (histone H3 lysine 9 methylation) in vitro<sup>1</sup>. It facilitates the reactivation of pluripotency genes and induces passive demethylation, thus promoting reprogramming<sup>2</sup>. BIX-01294, in combination with BAY K8644 (a calcium channel agonist), was found to improve reprogramming efficiencies of Oct4-Klf4-(OK)-infected neural progenitor cells<sup>3</sup>. It induced apoptosis and decreased proliferation, mobility and invasion in human neuroblastoma cells<sup>4</sup>.

### References

1. S Kubicek et al. Mol. Cell 2007 25:473
2. D Huangfu et al. Nat. Biotechnol. 2008 26:795
3. Y Shi et al. Cell Stem Cell 2008 2:525
4. Z Lu et al. Anticancer Drugs 2013 24:484

### Molecular Structure



## BIX-01294 3HCl

Epigenetic Enzyme Inhibitor

Catalog #	S344-903
CAS #	935693-62-2
Lot #	P3320-5
Purity	>98%
Format	Off-white crystalline powder
Solubilization	May be dissolved in water (50 mg/ml) or DMSO (50 mg/ml)
Storage & Shipping	Store desiccated as supplied at ambient for up to 2 years. Store solutions at -20°C for up to 3 months.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**