
BIX 01294 trihydrochloride hydrate

(Catalog No. M64000)

Alternative Names

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

Description

Selectively impairs the G9a histone lysine methyltransferase (HMTase) and the generation of H3K9me2 in vitro. BIX-01294 enables neural progenitor cells (NPCs) to be reprogrammed in the absence of exogenous Oct4.

CAS Number

935693-62-2

Chemical Formula

$C_{28}H_{38}N_6O_2$

Molecular Weight

490.6

Size

1 mg

Purity

≥95%

Solubility

Soluble in DMSO (5mg/ml).

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
BIX 01294 trihydrochloride hydrate	1 mg	M64000 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.