

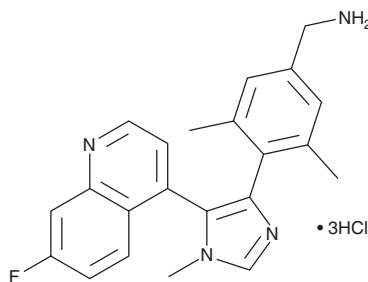
PRODUCT INFORMATION



BI-9321 (hydrochloride)

Item No. 27815

CAS Registry No.: 2387510-87-2
Formal Name: 4-[5-(7-fluoro-4-quinolinyl)-1-methyl-1H-imidazol-4-yl]-3,5-dimethylbenzenemethanamine, trihydrochloride
MF: C₂₂H₂₁FN₄ • 3HCl
FW: 469.8
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

BI-9321 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the BI-9321 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. BI-9321 (hydrochloride) is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

BI-9321 is an antagonist of the histone-lysine N-methyltransferase NSD3 (IC₅₀ = 0.203 μM in a time-resolved FRET (TR-FRET) assay).¹ It selectively binds to the NSD3 PWWP1 domain over 14 other PWWP domains, including NSD2-PWWP1 and NSD3-PWWP2, at 100 μM. BI-9321 is also selective for NSD3 over 20 methyllysine-binding proteins in differential static light scattering (DSL) assays at 500 μM, 35 DNA or RNA methyltransferases in enzyme activity assays at 10 μM, and 31 kinases in enzyme activity assays at 10 μM. It decreases expression of mRNA encoding Myc in MOLM-13 acute myeloid leukemia (AML) cells when used at a concentration of 10 μM for six hours and reduces the viability of MOLM-13 cells after six days (IC₅₀ = 26.8 μM).

Reference

1. Böttcher, J., Dilworth, D., Reiser, U., *et al.* Fragment-based discovery of a chemical probe for the PWWP1 domain of NSD3. *Nat. Chem. Biol.* **15**(8), 822-829 (2019).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/06/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM