

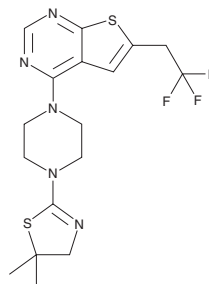
PRODUCT INFORMATION



MI-2-2

Item No. 41109

CAS Registry No.: 1454920-20-7
Formal Name: 4-[4-(4,5-dihydro-5,5-dimethyl-2-thiazolyl)-1-piperazinyl]-6-(2,2,2-trifluoroethyl)-thieno[2,3-d]pyrimidine
MF: C₁₇H₂₀F₃N₅S₂
FW: 415.5
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

MI-2-2 is supplied as a solid. A stock solution may be made by dissolving the MI-2-2 in the solvent of choice, which should be purged with an inert gas. MI-2-2 is soluble (≥10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

MI-2-2 is an inhibitor of the protein-protein interaction between menin and mixed-lineage leukemia 1 (MLL1).¹ It binds to menin (K_d = 22 nM) and inhibits the interaction between menin and MLL1 in a cell-free assay (IC₅₀ = 46 nM). MI-2-2 reduces cell growth (GI₅₀ = 3 μM) and induces cell cycle arrest at the G₁ phase in MV4-11 cells.

Reference

1. Shi, A., Murai, M.J., He, S., *et al.* Structural insights into inhibition of the bivalent menin-MLL interaction by small molecules in leukemia. *Blood* **120**(23), 4461-4469 (2012).

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, 06/11/2024

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