

# PRODUCT INFORMATION

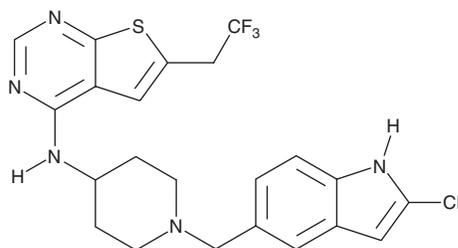


## MI-136

Item No. 19245

**CAS Registry No.:** 1628316-74-4  
**Formal Name:** 5-[[4-[[6-(2,2,2-trifluoroethyl)thieno[2,3-d]pyrimidin-4-yl]amino]-1-piperidinyl]methyl]-1H-indole-2-carbonitrile

**MF:** C<sub>23</sub>H<sub>21</sub>F<sub>3</sub>N<sub>6</sub>S  
**FW:** 470.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 220, 280 nm  
**Supplied as:** A crystalline 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-136 is supplied as a crystalline solid. A stock solution may be made by dissolving the MI-136 in the solvent of choice. MI-136 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of MI-136 in ethanol is approximately 10 mg/ml and approximately 1 mg/ml in DMSO and DMF.

### Description

MI-136 is an inhibitor of the menin-MLL interaction (IC<sub>50</sub> = 31 nM; K<sub>d</sub> = 23.6 nM).<sup>1,2</sup> By blocking MLL recruitment predominantly at the level of the menin-MLL interaction, it can inhibit androgen receptor (AR)-mediated transcription.<sup>3</sup> At 40 mg/kg, MI-136 has been shown to block AR signaling, inhibiting the growth of castration-resistant tumors in mice.<sup>3</sup>

### References

1. Borkin, D., Pollock, J., Kempinska, K., *et al.* Property focused structure-based optimization of small molecule inhibitors of the protein-protein interaction between menin and mixed lineage leukemia (MLL). *J. Med. Chem.* **59**(3), 892-913 (2016).
2. Borkin, D., He, S., Miao, H., *et al.* Pharmacologic inhibition of the menin-MLL interaction blocks progression of MLL leukemia *in vivo*. *Cancer Cell* **27**, 589-602 (2015).
3. Malik, R., Khan, A.P., Asangani, I.A., *et al.* Targeting the MLL complex in castration resistant prostate cancer. *Nat. Med.* **21**(4), 344-352 (2015).

#### 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, 12/19/2022

#### 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