

PRODUCT INFORMATION

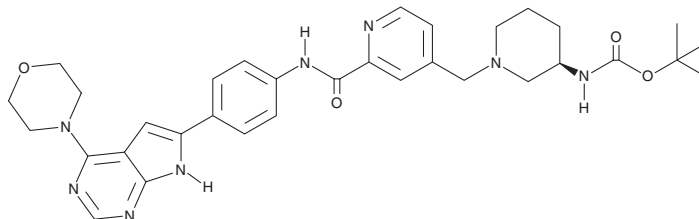


Menin-MLL Inhibitor 20

Item No. 40782

CAS Registry No.: 2448173-47-3
Formal Name: N-[(3R)-1-[[[2-[[[4-(4-morpholinyl)-7H-pyrrolo[2,3-d]pyrimidin-6-yl]phenyl]amino]carbonyl]-4-pyridinyl]methyl]-3-piperidinyl]-carbamic acid, 1,1-dimethylethyl ester

MF: C₃₃H₄₀N₈O₄
FW: 612.7
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

Menin-MLL inhibitor 20 is supplied as a solid. A stock solution may be made by dissolving the menin-MLL inhibitor 20 in the solvent of choice, which should be purged with an inert gas. Menin-MLL inhibitor 20 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Menin-MLL inhibitor 20 is a synthetic intermediate in the production of the irreversible inhibitor of the protein-protein interaction between MLL and menin, BMF-219, which induces cytotoxicity in cancer cells *in vitro*.¹

References

1. Butler, T., Palmer, J., Upasani, R., *et al.* Irreversible inhibitors of menin-MLL interaction. *Biomea Fusion, LLC*. WO 2020/142557 A1 (2020).

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, 05/06/2024

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