

PRODUCT INFORMATION

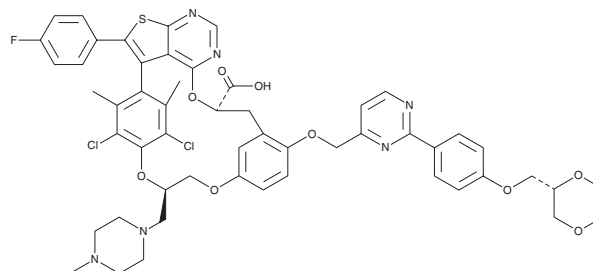


ABBV-467

Item No. 44289

CAS Registry No.: 2287186-66-5
Formal Name: (7R,16R)-19,23-dichloro-10-[[2-[4-[(2R)-1,4-dioxan-2-ylmethoxy]phenyl]-4-pyrimidinyl]methoxy]-1-(4-fluorophenyl)-7,8,15,16-tetrahydro-20,22-dimethyl-16-[(4-methyl-1-piperazinyl)methyl]-18,21-etheno-9,13-metheno-13H-6,14,17-trioxa-2-thia-3,5-diazacyclononadec[cd]indene-7-carboxylic acid

MF: C₅₃H₅₁Cl₂FN₆O₉S
FW: 1,038.0
Purity: ≥98%
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

ABBV-467 is supplied as a crystalline solid. A stock solution may be made by dissolving the ABBV-467 in the solvent of choice, which should be purged with an inert gas. ABBV-467 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in a 1:1 solution of acetonitrile:water.

Description

ABBV-467 is an inhibitor of myeloid cell leukemia-1 (Mcl-1; $K_i = <0.1$ nM).¹ It is selective for Mcl-1 over Bcl-2, Bcl-xL, Bcl-W, and Bcl-2-A1 (K_i s = >642, >376, >247, and >402 nM, respectively). ABBV-467 reduces the viability of AMO1, NCI H929, and MV4-11 myeloma cells (EC_{50} s = 0.16, 0.47, and 3.91 nM, respectively). *In vivo*, ABBV-467 (3.13, 6.25, and 12.5 mg/kg) reduces tumor growth in an AMO1 mouse xenograft model.

Reference

1. Yuda, J., Will, C., Phillips, D.C., *et al.* Selective MCL-1 inhibitor ABBV-467 is efficacious in tumor models but is associated with cardiac troponin increases in patients. *Commun. Med. (Lond)* **3(1)**, 154 (2023).

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, 08/25/2025

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