

PRODUCT INFORMATION



PROTAC Mcl Degradar-1

Item No. 45150

CAS Registry No.: 2163793-38-0

Formal Name: N¹-[2-[6-[(4-bromophenyl)thio]-1,3-dioxo-1H-benz[de]isoquinolin-2(3H)-yl]ethyl]-N⁶-[6-[[2-(2,6-dioxo-3-piperidiny)]-2,3-dihydro-1,3-dioxo-1H-isoindol-4-yl]amino]hexyl]-hexanediamide

Synonym: Proteolysis-targeting Chimera Mcl Degradar-1

MF: C₄₅H₄₅BrN₆O₈S

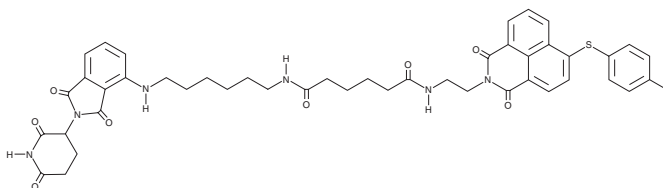
FW: 909.9

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

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

Description

PROTAC Mcl degrader-1 is a proteolysis-targeting chimera (PROTAC) composed of a myeloid cell leukemia-1 (Mcl-1) inhibitor conjugated to the cereblon (CRBN) inhibitor pomalidomide (Item No. 19877) via an alkyl linker.¹ It induces degradation of Mcl-1 in HeLa cells with a half-maximal degradation concentration (DC₅₀) value of 0.76 μM. PROTAC Mcl degrader-1 induces selective degradation of Mcl-1 over Bcl-2, Bcl-xL, murine double minute 2 (MDM2), and tubulin in HeLa cells when used at a concentration of 10 μM. It inhibits the growth of H23 non-small cell lung cancer (NSCLC) cells in a concentration-dependent manner.

Reference

1. Wang, Z., He, N., Guo, Z., *et al.* Proteolysis targeting chimeras for the selective degradation of Mcl-1/Bcl-2 derived from nonselective target binding ligands. *J. Med. Chem.* **62**(17), 8152-8163 (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, 03/05/2026

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