

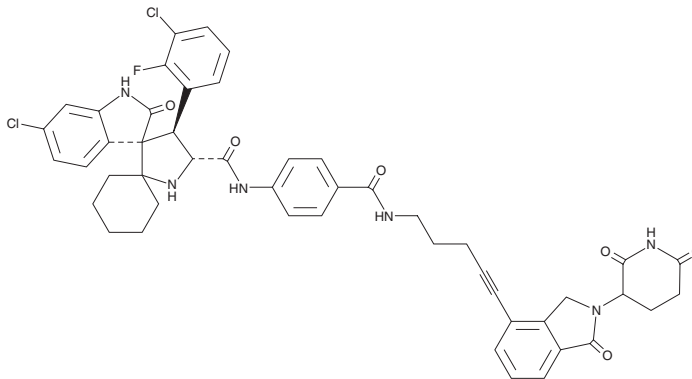
PRODUCT INFORMATION



MD-224

Item No. 40960

CAS Registry No.: 2136247-12-4
Formal Name: (3'R,4'S,5'R)-6''-chloro-4'-(3-chloro-2-fluorophenyl)-N-[4-[[[5-[2-(2,6-dioxo-3-piperidinyl)-2,3-dihydro-1-oxo-1H-indol-4-yl]-4-pentyn-1-yl]amino]carbonyl]phenyl]-1'',2''-dihydro-2''-oxo-dispiro[cyclohexane-1,2'-pyrrolidine-3',3''-[3H]indole]-5'-carboxamide



MF: C₄₈H₄₃Cl₂FN₆O₆
FW: 889.8
Purity: ≥95%
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

MD-224 is supplied as a solid. A stock solution may be made by dissolving the MD-224 in the solvent of choice, which should be purged with an inert gas. MD-224 is sparingly soluble (1-10 mg/ml) in acetonitrile and DMSO.

Description

MD-224 is a proteolysis-targeting chimera (PROTAC) that contains the murine double minute 2 (MDM2) inhibitor MI-1242 conjugated to lenalidomide (Item No. 14643).¹ It induces degradation of MDM2, which increases p53 levels, in RS4;11 acute lymphoblastic leukemia (ALL) cells when used at a concentration of 10 nM. MD-224 also inhibits the growth of RS4;11 cells (IC₅₀ = 1.5 nM). *In vivo*, MD-224 (50 mg/kg every other day for three weeks) induces complete tumor regression and does not affect body weight in an RS4;11 mouse xenograft model.

Reference

1. Li, Y., Yang, J., Aguilar, A., *et al.* Discovery of MD-224 as a first-in-class, highly potent, and efficacious proteolysis targeting chimera murine double minute 2 degrader capable of achieving complete and durable tumor regression. *J. Med. Chem.* **62**(2), 448-466 (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, 06/20/2025

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