

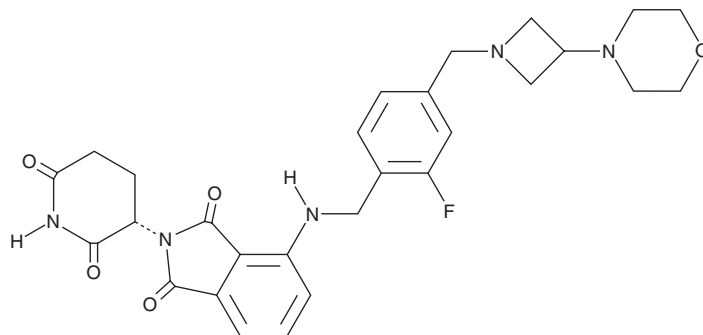
PRODUCT INFORMATION



Golcadomide

Item No. 44114

CAS Registry No.: 2379572-34-4
Formal Name: 2-[(3S)-2,6-dioxo-3-piperidiny]-4-[[[2-fluoro-4-[[3-(4-morpholinyl)-1-azetidiny]methyl]phenyl]methyl]amino]-1H-isoindole-1,3(2H)-dione
Synonym: CC-99282
MF: C₂₈H₃₀FN₅O₅
FW: 535.6
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

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

Description

Golcadomide is a cereblon (CRBN) E3 ligase modulator (CELMoD).¹ It decreases protein levels of Ikaros family zinc finger protein 1 (IKZF1) and IFN regulatory factor 4 (IRF4) in JSC-1, BCBL-1, Daudi, and BL-41 lymphoma cells when used at a concentration of 0.03 μM and decreases the viability of these cells in a concentration-dependent manner. Golcadomide (0.01 μM) increases the levels of intracellular adhesion molecule-1 (ICAM-1) and CD86 in Daudi, Raji, and Epstein-Barr virus-positive BL-41 cells.

Reference

1. Shrestha, P., Treco, E.N., Davis, D.A., *et al.* Novel cereblon-binding immunomodulators have increased potency against gammaherpesvirus-associated lymphomas in vitro. *J. Med. Virol.* **97(8)**, e70537 (2025).

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