

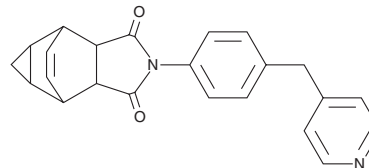
PRODUCT INFORMATION



DCZ0415

Item No. 42513

CAS Registry No.: 2242470-43-3
Formal Name: 4,4a,5,5a,6,6a-hexahydro-2-[4-(4-pyridinylmethyl)phenyl]-4,6-ethenocycloprop[f]isoindole-1,3(2H,3aH)-dione
MF: C₂₃H₂₀N₂O₂
FW: 356.4
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

DCZ0415 is supplied as a solid. A stock solution may be made by dissolving the DCZ0415 in the solvent of choice, which should be purged with an inert gas. DCZ0415 is sparingly soluble (1-10 mg/ml) in ethanol and soluble (≥10 mg/ml) in DMSO.

Description

DCZ0415 is an inhibitor of thyroid hormone receptor interactor 13 (TRIP13).¹ It binds to TRIP13 (K_d = 2.42 μM) and reduces non-homologous end joining (NHEJ) repair of DNA in a reporter assay. DCZ0415 (10 and 20 μM) impairs DNA repair in ARP-1 and NCI H929 multiple myeloma cells. It reduces the growth of a variety of multiple myeloma cells, including RPMI-8226, ARP-1, and ARK cells (IC₅₀s = 7.4, 5.12, and 3.25 μM, respectively) and induces apoptosis in ARP-1 cells. DCZ0415 (50 mg/kg per day) induces intratumoral apoptosis and impairs intratumoral DNA repair, reduces tumor volume, and increases survival in an NCI H929 mouse xenograft model.

Reference

1. Wang, Y., Huang, J., Li, B., *et al.* A small-molecule inhibitor targeting TRIP13 suppresses multiple myeloma progression. *Cancer Res.* **80**(3), 536-548 (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, 01/23/2025

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