

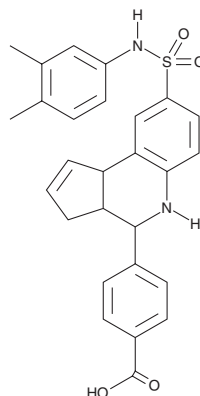
PRODUCT INFORMATION



MX69

Item No. 27224

CAS Registry No.: 1005264-47-0
Formal Name: 4-[8-[[[(3,4-dimethylphenyl)amino]sulfonyl]-3a,4,5,9b-tetrahydro-3H-cyclopenta[c]quinolin-4-yl]-benzoic acid
MF: C₂₇H₂₆N₂O₄S
FW: 474.6
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

MX69 is supplied as a solid. A stock solution may be made by dissolving the MX69 in the solvent of choice, which should be purged with an inert gas. MX69 is soluble in the organic solvent DMSO.

Description

MX69 is an inhibitor of the interaction between the ubiquitin ligase MDM2 protein and XIAP mRNA.¹ It binds to the RING domain of MDM2 ($K_d = 2.34 \mu\text{M}$) and decreases the levels of MDM2 and XIAP in a concentration-dependent manner without affecting the levels of Bcl-2, cIAP-1, and cIAP-2. It induces autoubiquitination and degradation of MDM2 and increases the half-life of p53 in EU-1 cells. MX69 is cytotoxic to a variety of acute lymphoblastic leukemia (ALL) and neuroblastoma cancer cell lines but not to human normal bone marrow mononuclear (NBMM) cells. It increases survival in an EU-1 mouse model of leukemia when administered at a dose of 100 mg/kg.

Reference

1. Gu, L., Zhang, H., Liu, T., *et al.* Discovery of dual inhibitors of MDM2 and XIAP for cancer treatment. *Cancer Cell* **30(4)**, 623-636 (2016).

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, 11/08/2022

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