

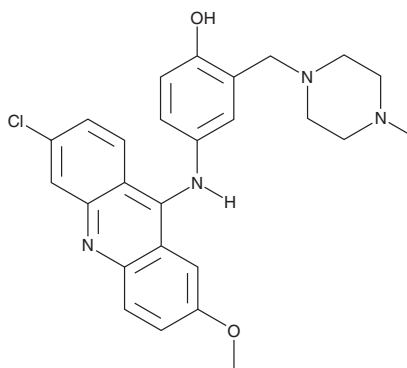
PRODUCT INFORMATION



HM03

Item No. 40908

CAS Registry No.: 500565-15-1
Formal Name: 4-[[6-chloro-2-methoxy-9-acridinyl)amino]-2-[[4-methyl-1-piperazinyl)methyl]-phenol
MF: C₂₆H₂₇ClN₄O₂
FW: 463.0
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

HM03 is supplied as a solid. A stock solution may be made by dissolving the HM03 in the solvent of choice, which should be purged with an inert gas. HM03 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

HM03 is an inhibitor of the protein-protein interaction between the DNA excision repair protein ERCC-1 and the DNA repair endonuclease XPF (IC₅₀ = 1.86 μM).¹ It is also an inhibitor of D-alanine-D-alanine ligase B (DdlB; K_i = 218 μM for the *E. coli* enzyme).² HM03 is active against *E. coli* and *S. aureus* (MICs = 64 and 32 μg/ml, respectively). It potentiates cytotoxicity induced by cisplatin (Item No. 13119) in A549 lung and HCT116 colon cancer cells when used at a concentration of 100 μM.³

References

1. Elmenoufy, A.H., Gentile, F., Jay, D., *et al.* Targeting DNA repair in tumor cells via inhibition of ERCC1-XPF. *J. Med. Chem.* **62**(17), 7684-7696 (2019).
2. Kovac, A., Konc, J., Vehar, B., *et al.* Discovery of new inhibitors of D-alanine:D-alanine ligase by structure-based virtual screening. *J. Med. Chem.* **51**(23), 7442-7448 (2008).
3. Jordheim, L.P., Barakat, K.H., Heinrich-Balard, L., *et al.* Small molecule inhibitors of ERCC1-XPF protein-protein interaction synergize alkylating agents in cancer cells. *Mol. Pharmacol.* **84**(1), 12-24 (2013).

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, 07/10/2024

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