

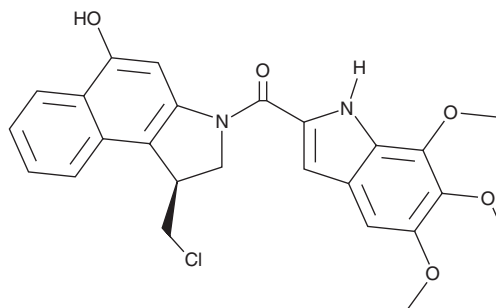
PRODUCT INFORMATION



Duocarmycin TM

Item No. 43594

CAS Registry No.: 157922-77-5
Formal Name: 1S-(chloromethyl)-2,3-dihydro-3-[[5,6,7-trimethoxy-1H-indol-2-yl]carbonyl]-1H-benz[e]indol-5-ol
MF: C₂₅H₂₃ClN₂O₅
FW: 466.9
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

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

Description

Duocarmycin TM is a DNA alkylating agent.¹ It is cytotoxic to L1210 murine skin lymphocytic leukemia cells (IC₅₀ = 30 pM). Duocarmycin TM is also active against *P. falciparum* (IC₅₀ = 2-8 pM).²

References

1. Ham, Y.-W. and Boger, D.L. A powerful selection assay for mixture libraries of DNA alkylating agents. *J. Am. Chem. Soc.* **126**(30), 9194-9195 (2004).
2. Alder, A., Struck, N.S., Xu, M., et al. A non-reactive natural product precursor of the duocarmycin family has potent and selective antimalarial activity. *Cell Chem. Biol.* **29**(5), 840-853 (2022).

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, 08/20/2025

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