

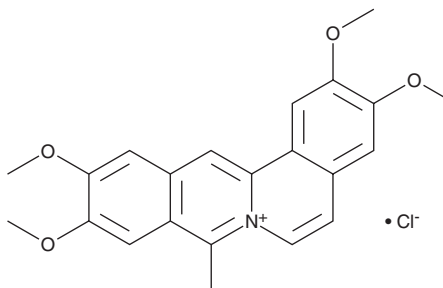
PRODUCT INFORMATION



Coralyne (chloride)

Item No. 45715

CAS Registry No.: 38989-38-7
Formal Name: 2,3,10,11-tetramethoxy-8-methyl-dibenzo[a,g]quinolizinium, monochloride
MF: $C_{22}H_{22}NO_4 \cdot Cl$
FW: 399.9
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Coralyne (chloride) is supplied as a solid. A stock solution may be made by dissolving the coralyne (chloride) in the solvent of choice, which should be purged with an inert gas. Coralyne (chloride) is slightly soluble (0.1-1 mg/ml) in a 1:1 solution of acetonitrile:water.

Description

Coralyne is a DNA topoisomerase I poison.¹ It inhibits DNA topoisomerase I via stabilization of the DNA-enzyme complex and inhibition of DNA relaxation. It is selective for DNA topoisomerase I over DNA topoisomerase II at 100 $\mu g/ml$.² Coralyne (100-400 mg/kg) increases survival in P388 and L1210 murine leukemia models.³

References

1. Wang, L.-K., Rogers, B.D., and Hecht, S.M. Inhibition of topoisomerase I function by coralyne and 5,6-dihydrocoralyne. *Chem. Res. Toxicol.* **9**(1), 75-83 (1996).
2. Gatto, B., Sanders, M.M., Yu, C., *et al.* Identification of topoisomerase I as the cytotoxic target of the protoberberine alkaloid coralyne. *Cancer Res.* **56**(12), 2795-2800 (1996).
3. Zee-Cheng, K.-Y., Paull, K.D., and Cheng, C.C. Experimental antileukemic agents. Coralyne, analogs, and related compounds. *J. Med. Chem.* **17**(3), 347-351 (1974).

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, 06/04/2026

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