

PRODUCT INFORMATION



10-hydroxy Camptothecin

Item No. 46297

CAS Registry No.: 64439-81-2

Formal Name: 4-ethyl-4,9-dihydroxy-1H-pyrano[3',4':6,7]indolizino[1,2-b]quinoline-3,14(4H,12H)-dione

Synonyms: DL-10-hydroxy Camptothecin, 10-HCPT

MF: $C_{20}H_{16}N_2O_5$

FW: 364.4

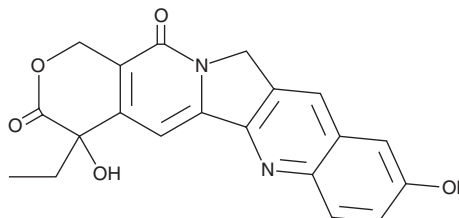
Purity: $\geq 98\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Plant/*Camptotheca acuminata*



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

Laboratory Procedures

10-hydroxy Camptothecin is supplied as a solid. A stock solution may be made by dissolving the 10-hydroxy camptothecin in the solvent of choice, which should be purged with an inert gas. 10-hydroxy Camptothecin is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

10-hydroxy Camptothecin is an alkaloid originally isolated from *C. acuminata* that has DNA topoisomerase I inhibitory and anticancer activities.^{1,2} It inhibits DNA topoisomerase I ($IC_{50} = 1 \mu\text{M}$), reduces the proliferation of L1210 leukemia cells ($IC_{50} = 18 \text{ nM}$), and increases survival by 78% in an L1210 murine leukemia model at $21 \mu\text{mol/kg}$.²

References

1. Wani, M.C. and Wall, M.E. Plant antitumor agents. II. Structure of two new alkaloids from *Camptotheca acuminata*. *J. Org. Chem.* **34**(5), 1364-1367 (1968).
2. Kingsbury, W.D., Boehm, J.C., Jakas, D.R., et al. Synthesis of water-soluble (aminoalkyl)camptothecin analogues: Inhibition of topoisomerase I and antitumor activity. *J. Med. Chem.* **34**(1), 98-107 (1991).

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, 08/04/2026

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