

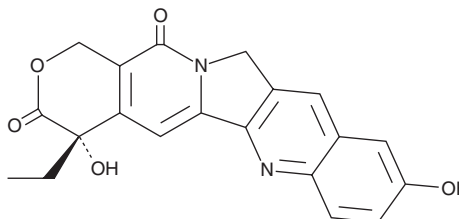
PRODUCT INFORMATION



(S)-10-hydroxy-Camptothecin

Item No. 14635

CAS Registry No.: 19685-09-7
Formal Name: (4S)-4-ethyl-4,9-dihydroxy-1H-pyrano[3',4':6,7]indolizino[1,2-b]quinoline-3,14(4H,12H)-dione
ChEMBL 273862, NSC 107124
Synonyms:
MF: C₂₀H₁₆N₂O₅
FW: 364.4
Purity: ≥98%
UV/Vis.: λ_{max}: 222, 267, 330, 383 nm
Supplied as: A crystalline 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

(S)-10-hydroxy-Camptothecin is supplied as a crystalline solid. A stock solution may be made by dissolving the (S)-10-hydroxy-camptothecin in the solvent of choice, which should be purged with an inert gas. (S)-10-hydroxy-Camptothecin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of (S)-10-hydroxy-camptothecin in these solvents is approximately 3 and 2 mg/ml, respectively.

(S)-10-hydroxy-Camptothecin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (S)-10-hydroxy-camptothecin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. (S)-10-hydroxy-Camptothecin has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(S)-10-hydroxy-Camptothecin is an inhibitor of topoisomerase I originally isolated from the Chinese tree *C. acuminata*. It is a member of the camptothecin family that demonstrates less toxicity than its parent compound.¹ (S)-10-hydroxy-Camptothecin has strong anti-tumor activity against a wide range of experimental tumors including L1210 leukemia cells (IC₅₀ = 1.15 μM).¹ *In vitro* treatment of human HepG2 cells with 5-20 μM (S)-10-hydroxy-camptothecin results in cell cycle arrest at the G₂/M phase.²

References

1. Yu, P., Xia, L., Zhao, J., *et al.* Synthesis and preliminary anticancer evaluation of 10-hydroxycamptothecin analogs. *Biol. Pharm. Bull.* **35**(8), 1295-1299 (2012).
2. Zhang, X.W., Jiang, J.F., and Xu, B. Differentiation-inducing action of 10-hydroxycamptothecin on human hepatoma Hep G2 cells. *Acta Pharmacol. Sin.* **21**(4), 364-368 (2000).

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, 11/23/2022

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