

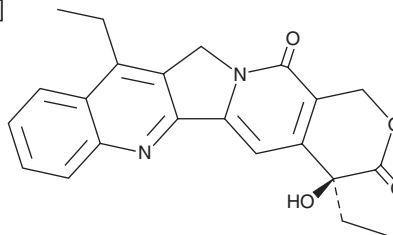
PRODUCT INFORMATION



7-ethyl Camptothecin

Item No. 34607

CAS Registry No.: 78287-27-1
Formal Name: (4S)-4,11-diethyl-4-hydroxy-1H-pyrano[3',4':6,7]indolizino[1,2-b]quinoline-3,14(4H,12H)-dione
Synonym: SN-22
MF: C₂₂H₂₀N₂O₄
FW: 376.4
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 254, 359 nm
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

7-ethyl Camptothecin is supplied as a solid. A stock solution may be made by dissolving the 7-ethyl camptothecin in the solvent of choice, which should be purged with an inert gas. 7-ethyl Camptothecin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 7-ethyl camptothecin in these solvents is approximately 1 mg/ml.

Description

7-ethyl Camptothecin is a derivative of camptothecin (Item No. 11694) and an anticancer agent.^{1,2} It inhibits wild-type but not Gly533 mutant topoisomerase I in RPMI-8402 and CPT-K5 cells, respectively (IC₅₀s = 5.48 and >2,500 nM, respectively).¹ 7-ethyl Camptothecin is cytotoxic to ADJ/PC6 SN-38-sensitive and PC6/SN2-5H2 SN-38-resistant cells (IC₅₀s = 0.85 and 2.98 nM, respectively). *In vivo*, 7-ethyl camptothecin (0.25-3 mg/kg per day) increases survival in Ehrlich, MM46, Meth A, L1210, P388, and B16/F10 mouse ascites tumor models.²

References

1. Yoshikawa, M., Ikegami, Y., Hayasaka, S., *et al.* Novel camptothecin analogues that circumvent ABCG2-associated drug resistance in human tumor cells. *Int. J. Cancer* **110**(6), 921-927 (2004).
2. Kunitomo, T., Nitta, K., Tanaka, T., *et al.* Antitumor activity of a new camptothecin derivative, SN-22, against various murine tumors. *J. Pharmacobiodyn.* **10**(3), 148-151 (1987).

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, 11/22/2022

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