

PRODUCT INFORMATION



Teniposide

Item No. 14425

CAS Registry No.: 29767-20-2
Formal Name: (5R,5aR,8aR,9S)-5,8,8a,9-tetrahydro-5-(4-hydroxy-3,5-dimethoxyphenyl)-9-[[4,6-O-[(R)-2-thienylmethylene]-β-D-glucopyranosyl]oxy]-furo[3,4':6,7]naphtho[2,3-d]-1,3-dioxol-6(5aH)-one

Synonyms: NSC 122819, VM-26

MF: C₃₂H₃₂O₁₃S

FW: 656.7

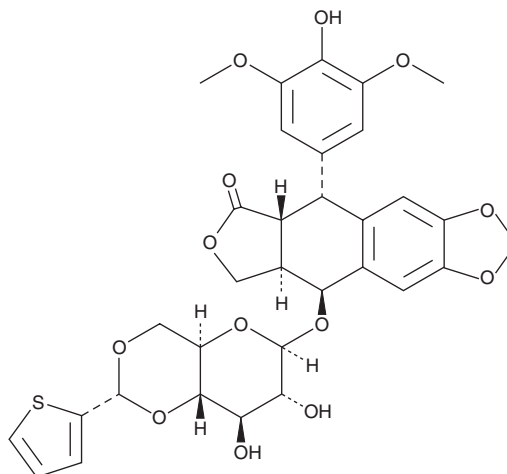
Purity: ≥95%

UV/Vis.: λ_{max}: 284 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

Teniposide is supplied as a crystalline solid. A stock solution may be made by dissolving the teniposide in the solvent of choice, which should be purged with an inert gas. Teniposide is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of teniposide in these solvents is approximately 5 mg/ml.

Teniposide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, teniposide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Teniposide has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Teniposide is a topoisomerase II poison.¹⁻⁴ It promotes calf thymus topoisomerase II-mediated DNA cleavage when used at concentrations ranging from 0.04 to 40 μg/ml.¹ Teniposide induces single- and double-stranded DNA breaks in a concentration-dependent manner.² It inhibits the proliferation of A549 lung cancer cells (IC₅₀ = 0.7 μM). Teniposide induces cell cycle arrest at the S phase and apoptosis in Tca8113 tongue squamous cell carcinoma cells when used at a concentration of 5 mg/L and induces cell cycle arrest at the G₂/M phase when used at 0.15 mg/L.³ It reduces tumor growth and inhibits metastasis in a 3LL murine Lewis lung carcinoma model when administered at a dose of 6.5 mg/kg.⁴ Formulations containing teniposide have been used in combination therapy in the treatment of refractory childhood acute lymphoblastic leukemia.

References

1. Ross, W., Rowe, T., Glisson, B., *et al.* Role of topoisomerase II in mediating epipodophyllotoxin-induced DNA cleavage. *Cancer Res.* **44(12 Pt. 1)**, 5857-5860 (1984).
2. Long, B.H., Musial, S.T., and Brattain, M.G. Single- and double-strand DNA breakage and repair in human lung adenocarcinoma cells exposed to etoposide and teniposide. *Cancer Res.* **45(7)**, 3106-3112 (1985).
3. Li, J., Chen, W., Zhang, P., *et al.* Topoisomerase II trapping agent teniposide induces apoptosis and G₂/M or S phase arrest of oral squamous cell carcinoma. *World J. Surg. Oncol.* **4(1)**, 41 (2006).
4. Colombo, T., Broggin, M., Vaghi, M., *et al.* Comparison between VP 16 and VM 26 in Lewis lung carcinoma of the mouse. *Eur. J. Cancer Clin. Oncol.* **22(2)**, 173-179 (1986).

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