

PRODUCT INFORMATION



(-)-Triptonide

Item No. 29156

CAS Registry No.: 38647-11-9
Formal Name: (3bS,4aS,5aS,6aS,7aS,7bS,8aS,8bS)-3b,4,4a,7a,7b,8b,9,10-octahydro-8b-methyl-6a-(1-methylethyl)trioxireno[4b,5:6,7:8a,9]phenanthro[1,2-c]furan-1,6(3H,6aH)-dione

Synonyms: NSC 165677, Triptonide

MF: C₂₀H₂₂O₆

FW: 358.4

Purity: ≥98%

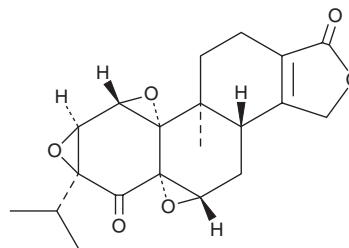
UV/Vis.: λ_{max}: 215 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Tripterygium wilfordii* Hook F



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

Laboratory Procedures

(-)-Triptonide is supplied as a crystalline solid. A stock solution may be made by dissolving the (-)-triptonide in the solvent of choice, which should be purged with an inert gas. (-)-Triptonide is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of (-)-triptonide in these solvents is approximately 10 mg/ml. (-)-Triptonide is slightly soluble in ethanol.

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

Description

(-)-Triptonide is a diterpenoid that has been found in *T. wilfordii* and has diverse biological activities, including immunosuppressant and anticancer properties.^{1,2} It inhibits IL-2 production and DNA synthesis in phytohemagglutinin-stimulated T cells (EC₅₀s = 2.46 and 3.2 ng/ml, respectively).¹ (-)-Triptonide inhibits the proliferation of PC3, DU145, and LNCaP cells (IC₅₀s = 11.96, 10.26, 12.01 nM, respectively).² It induces cell cycle arrest at the G₂/M phase and increases apoptosis in PC3 cells in a concentration-dependent manner. (-)-Triptonide (10 mg/kg per day) reduces tumor growth in a PC3 mouse xenograft model.

References

1. Tao, X., Cai, J.J., and Lipsky, P.E. The identity of immunosuppressive components of the ethyl acetate extract and chloroform methanol extract (T2) of tripterygium wilfordii Hook. F. *J. Pharmacol. Exp. Ther.* **272**(3), 1305-1312 (1995).
2. Dong, F., Yang, P., Wang, R., *et al.* Triptonide acts as a novel antiprostata cancer agent mainly through inhibition of mTOR signaling pathway. *Prostate* **79**(11), 1284-1293 (2019).

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, 11/02/2022

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