

# PRODUCT INFORMATION



## Triptolide

Item No. 11973

**CAS Registry No.:** 38748-32-2  
**Formal Name:** 3bS,4,4aS,6R,6aR,7aS,7bS,8bS,9,10-decahydro-6-hydroxy-8b-methyl-6a-(1-methylethyl)-trioxireno[4b,5aS:6,7:8aS,9]phenanthro[1,2-c]furan-1(3H)-one

**Synonyms:** NSC 163062, PG 490

**MF:** C<sub>20</sub>H<sub>24</sub>O<sub>6</sub>

**FW:** 360.4

**Purity:** ≥98%

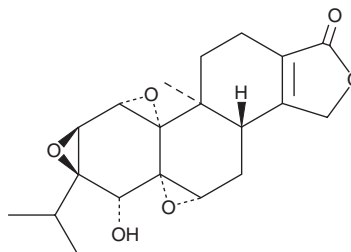
**UV/Vis.:** λ<sub>max</sub>: 217 nm

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Plant/*Tripterygium wilfordii*



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

### Laboratory Procedures

Triptolide is supplied as a crystalline solid. A stock solution may be made by dissolving the triptolide in the solvent of choice, which should be purged with an inert gas. Triptolide is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of triptolide in these solvents is approximately 11 and 12 mg/ml, respectively.

Triptolide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, triptolide should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Triptolide 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

Triptolide is a diterpenoid triepoxide that has been found in *T. wilfordii* and has diverse biological activities.<sup>1-4</sup> It inhibits dCTP pyrophosphatase 1 (K<sub>i</sub> = 168 μM), an enzyme that prevents halogenated nucleotides from entering DNA synthesis.<sup>1</sup> It reduces viability of A549, H1299, NCI H520, NCI H1650, and H1975 human non-small cell lung cancer (NSCLC) cells when used at a concentration of 50 nM.<sup>2</sup> Intranasal administration of liposomes containing triptolide (0.4 mg/kg) reduces tumor growth in a rat orthotopic model of lung cancer. Triptolide (0.2 mg/kg) reduces intestinal inflammation and prevents colon shortening in a mouse model of colitis induced by dextran sulfate (Item No. 23250).<sup>3</sup> It also inhibits skin allograft rejection and increases graft survival time in mice when administered at a dose of 0.1 mg/kg per day post-transplantation.<sup>4</sup>

### References

1. Corson, T.W., Cavga, H., Aberle, N., *et al.* Triptolide directly inhibits dCTP pyrophosphatase. *ChemBioChem* **12(11)**, 1767-1773 (2011).
2. Song, J.M., Molla, K., Anandharaj, A., *et al.* Triptolide suppresses the *in vitro* and *in vivo* growth of lung cancer cells by targeting hyaluronan-CD44/RHAMM signaling. *Oncotarget*. **8(16)**, 26927-26940 (2017).
3. Tang, B., Zhu, J., Zhang, B., *et al.* Therapeutic potential of triptolide as an anti-inflammatory agent in dextran sulfate sodium-induced murine experimental colitis. *Front. Immunol.* **11**, 592084 (2020).
4. Yang, S.-X., Gao, H.-L., Xie, S.-S., *et al.* Immunosuppression of triptolide and its effect on skin allograft survival. *Int. J. Immunopharmacol.* **14(6)**, 963-969 (1992).

#### 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, 12/01/2022

#### 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