

Triptolide

NF- κ B activation inhibitor

Catalog # ant-tpl

For research use only

Version # 12B24-MM

PRODUCT INFORMATION

Content:

Triptolide is supplied as a white solid.

- **ant-tpl:** 1 mg

Storage and stability:

- Triptolide is shipped at room temperature. Store as supplied at -20°C in a tightly sealed vial. Protect from light. Triptolide as a solid is stable for 6 months when properly stored.
- Once solubilized, prepare aliquots of triptolide and store at -20°C . Avoid repeated freeze-thaw cycles. Protect from light. Solubilized triptolide is stable for 3 months when properly stored.

Quality control

Purity : >98%

DESCRIPTION

Triptolide, a diterpenoid isolated from *Tripterygium wilfordii* hook F, has been used for centuries in traditional Chinese medicine to treat immune-related disorders. In addition to its anti-inflammatory and immunosuppressive activities, triptolide possesses potent antitumor properties. In a broad range of human tumor cells, Triptolide suppresses cell proliferation and induces apoptosis through caspase activation¹. At a molecular level, Triptolide inhibits global gene transcription by inducing degradation of RNA polymerase II (Pol II)², and by inhibiting the ATPase activity of XPB³, a subunit of the general transcription factor TFIIH. Triptolide interferes with a number of transcription factors including p53⁴, NF- κ B⁵, nuclear factor of activated T-cells (NFAT)⁵ and heat shock factor protein 1 (HSF-1)⁶.

1. **Carter BZ. et al., 2006.** Triptolide induces caspase-dependent cell death mediated via the mitochondrial pathway in leukemic cells. *Blood* 108: 630 - 637. 2. **Wang Y. et al., 2011.** Triptolide (TPL) inhibits global transcription by inducing proteasome-dependent degradation of RNA polymerase II (Pol II). *PLoS One*. 6(9):e23993. 3. **Titov D. et al., 2011.** XPB, a subunit of TFIIH, is a target of the natural product triptolide. *Nat Chem Biol*. 7(3):182-8. 4. **Chang, W.T. et al., 2001.** Triptolide and chemotherapy cooperate in tumor cell apoptosis. A role for the p53 pathway. *J. Biol. Chem.* 276, 2221–2227. 5. **Qiu D. et al., 1999.** Immunosuppressant PG490 (triptolide) inhibits T-cell interleukin-2 expression at the level of purine-box/nuclear factor of activated T-cells and NF-kappaB transcriptional activation. *J. Biol. Chem.* 274, 13443–13450. 6. **Westerheide SD. et al., 2006.** Triptolide, an Inhibitor of the Human Heat Shock Response That Enhances Stress-induced Cell Death. *J. Biol. Chem.*, 281: 9616 - 9622.

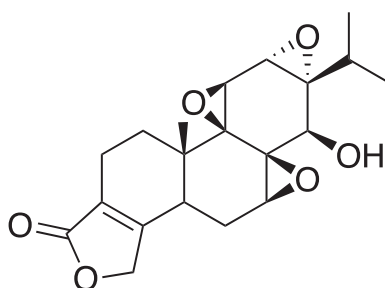
CHEMICAL PROPERTIES

CAS number: 38748-32-2

Formula: C₂₀H₂₄O₆

Molecular weight: 360.4

Solubility: DMSO, ethanol (10 mg/ml)



METHOD

Preparation of sterile stock solution (10 mM)

To obtain a 10 mM stock solution:

1. Add 280 μl DMSO to 1 mg Triptolide.
2. Vortex until complete solubilization.
3. Prepare aliquots of Triptolide and store at -20°C .

Working concentration: 10-100 nM

RELATED PRODUCT

Product	Catalog Code
Bay11-7082 (I κ B- α inhibitor)	tlrl-b82
Celastrol (NF- κ B inhibitor)	ant-cls

TECHNICAL SUPPORT

Toll free (US): 888-457-5873

Outside US: (+1) 858-457-5873

Europe: +33 562-71-69-39

E-mail: info@invivogen.com

Website: www.invivogen.com



3950 Sorrento Valley Blvd. Suite 100
San Diego, CA 92121 - USA