

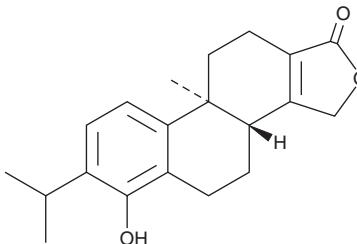
PRODUCT INFORMATION



Triptophenolide

Item No. 34390

CAS Registry No.: 74285-86-2
Formal Name: (3bR,9bS)-3b,4,5,9b,10,11-hexahydro-6-hydroxy-9b-methyl-7-(1-methylethyl)-phenanthro[1,2-c]furan-1(3H)-one
Synonyms: Hypolide, (+)-Triptophenolide
MF: C₂₀H₂₄O₃
FW: 312.4
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Salvia japonica*



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

Laboratory Procedures

Triptophenolide is supplied as a solid. A stock solution may be made by dissolving the triptophenolide in the solvent of choice, which should be purged with an inert gas. Triptophenolide is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of triptophenolide in these solvents is approximately 30 mg/ml.

Triptophenolide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, triptophenolide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Triptophenolide has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Triptophenolide is a diterpenoid that has been found in *T. wilfordii* and has antiandrogenic activity.¹ It is an antagonist of wild-type and mutant androgen receptors (ARs; IC₅₀s = 260, 388, 480, and 437 nM for wild-type, AR^{T877A}, AR^{F876L}, and AR^{W741C/T877A}, respectively, in transactivation assays). Triptophenolide (5 μM) inhibits the growth of, and reduces AR levels in, LNCaP prostate cancer cells.

Reference

1. He, Y., Wu, M., Liu, Y., *et al.* Identification of triptophenolide from *Tripterygium wilfordii* as a pan-antagonist of androgen receptor. *ACS Med. Chem. Lett.* **7**(12), 1024-1027 (2016).

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