

PRODUCT INFORMATION



Taccalonolide A

Item No. 40798

CAS Registry No.: 108885-68-3

Formal Name: (1 α ,2 α ,3 α ,5 α ,7 β ,11 α ,12 α ,15 α ,16 β ,24 β ,25S)-1,11,12,15-tetrakis(acetyloxy)-2,3-epoxy-7,23,25-trihydroxy-6-oxo- γ -lactone, 16,24-cycloergost-22-en-26-oic acid

MF: C₃₆H₄₆O₁₄

FW: 702.7

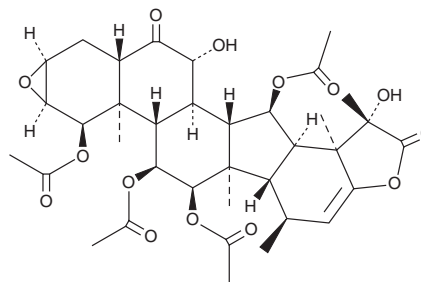
Purity: $\geq 98\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Plant/*Marsdenia officinalis*



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

Laboratory Procedures

Taccalonolide A is supplied as a solid. A stock solution may be made by dissolving the taccalonolide A in the solvent of choice, which should be purged with an inert gas. Taccalonolide A is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

Taccalonolide A is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Taccalonolide A is a taccalonolide that has been found in *T. chantrieri* and has anticancer activity.¹ It induces microtubule bundling, formation of abnormal mitotic spindles, and cell cycle arrest at the G₂/M phase in HeLa cervical cancer cells when used at a concentration of 3.5 μ M. Taccalonolide A decreases the proliferation of HeLa cells (IC₅₀ = 594 nM). *In vivo*, taccalonolide A increases the number of tumor free mice but reduces body weight in a 16/C murine breast cancer model.

Reference

1. Peng, J., Risinger, A.L., Fest, G.A., *et al.* Identification and biological activities of new taccalonolide microtubule stabilizers. *J. Med. Chem.* **54**(17), 6117-6124 (2011).

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, 07/15/2024

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