

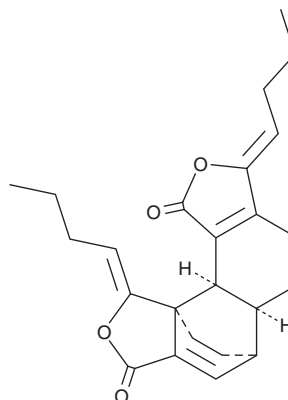
PRODUCT INFORMATION



Levistolide A

Item No. 34386

CAS Registry No.: 88182-33-6
Formal Name: (1Z,5S,5aS,8Z,10bS,10cS)-1,8-dibutylidene-5,5a,6,7,8,10b-hexahydro-1H-5,10c-ethanonaphtho[1,2-c:7,8-c']difuran-3,10-dione
Synonyms: (Z,Z')-Diligustilide, Z-Ligustilide Dimer
MF: C₂₄H₂₈O₄
FW: 380.5
Purity: ≥98%
UV/Vis.: λ_{max}: 231, 277 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Ligusticum chuanxiong*



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

Laboratory Procedures

Levistolide A is supplied as a solid. A stock solution may be made by dissolving the levistolide A in the solvent of choice, which should be purged with an inert gas. Levistolide A is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of levistolide A in these solvents is approximately 10, 20, and 30, respectively.

Levistolide A is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, levistolide A should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Levistolide A has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Levistolide A is a phthalide that has been found in *Angelica tenuissima* and has diverse biological activities.¹⁻⁴ It inhibits histamine release induced by phorbol 12-myristate 13-acetate (TPA; Item No. 10008014) and A23187 from HMC-1 human mast cell leukemia cells when used at a concentration of 10 μM.¹ Levistolide A (100 μM) induces the production of reactive oxygen species (ROS), endoplasmic reticulum (ER) stress, and apoptosis in HCT116 colon cancer cells.² It prevents ethanol-induced gastric ulcers in rats when administered at a dose of 10 mg/kg.³ Levistolide A (2 mg/kg) decreases aggregation of amyloid-β in brain and improves spatial and learning memory in the APP/PS1 transgenic mouse model of Alzheimer's disease.⁴

References

1. Choi, H.G., Je, I.-G., Kim, G.J., *et al.* Chemical constituents of the root of *Angelica tenuissima* and their anti-allergic inflammatory activity. *Nat. Prod. Commun.* **12(5)**, 779-780 (2017).
2. Yang, Y., Zhang, Y., Wang, L., *et al.* Levistolide A induces apoptosis via ROS-mediated ER stress pathway in colon cancer cells. *Cell. Physiol. Biochem.* **42(3)**, 929-938 (2017).
3. Velázquez-Moyado, J.A., Martínez-González, A., Linares, E., *et al.* Gastroprotective effect of diligustilide isolated from roots of *Ligusticum porteri* coulter & rose (Apiaceae) on ethanol-induced lesions in rats. *J. Ethnopharmacol.* **174**, 403-409 (2015).
4. Qu, X., Guan, P., Han, L., *et al.* Levistolide A attenuates Alzheimer's pathology through activation of the PPARγ pathway. *Neurotherapeutics* **18(1)**, 326-339 (2021).

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, 11/04/2022

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