

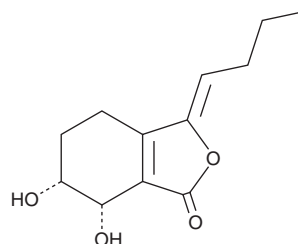
PRODUCT INFORMATION



Senkyunolide H

Item No. 35142

CAS Registry No.: 94596-27-7
Formal Name: (3Z,6R,7S)-rel-3-butylidene-4,5,6,7-tetrahydro-6,7-dihydroxy-1(3H)-isobenzofuranone
MF: C₁₂H₁₆O₄
FW: 224.3
Purity: ≥95%
UV/Vis.: λ_{max}: 274 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Angelica sinensis*



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

Laboratory Procedures

Senkyunolide H is supplied as a solid. A stock solution may be made by dissolving the senkyunolide H in the solvent of choice, which should be purged with an inert gas. Senkyunolide H is soluble in ethanol and DMSO.

Description

Senkyunolide H is a phthalide that has been found in *L. chuanxiong* and has diverse biological activities.¹⁻⁴ It increases levels of heme oxygenase-1 (HO-1) in HepG2 cells, as well as reduces hydrogen peroxide-induced increases in reactive oxygen species (ROS) and malondialdehyde (MDA) in the same cells when used at concentrations of 150 and 200 µg/ml.² Senkyunolide H inhibits the proliferation of primary mouse aorta smooth muscle cells (IC₅₀ = <0.1 µg/ml).³ It decreases bone resorption in ovariectomized mice when administered at a dose of 15 mg/kg.⁴

References

1. Yan, R., Li, S.-L., Chung, H.-S., *et al.* Simultaneous quantification of 12 bioactive components of *Ligusticum chuanxiong* Hort. by high-performance liquid chromatography. *J. Pharm. Biomed. Anal.* **37(1)**, 87-95 (2005).
2. Qi, H., Siu, S.O., Chen, Y., *et al.* Senkyunolides reduce hydrogen peroxide-induced oxidative damage in human liver HepG2 cells via induction of heme oxygenase-1. *Chem. Biol. Interact.* **183(3)**, 380-389 (2010).
3. Kobayashi, S., Mimura, Y., Notoya, K., *et al.* Antiproliferative effects of the traditional Chinese medicine *shimotsu-to*, its component cnidium rhizome and derived compounds on primary cultures of mouse aorta smooth muscle cells. *Jpn. J. Pharmacol.* **60(4)**, 397-401 (1992).
4. Yang, D., Liu, T., Jiang, G., *et al.* Senkyunolide H attenuates osteoclastogenesis and postmenopausal osteoporosis by regulating the NF-κB, JNK and ERK signaling pathways. *Biochem. Biophys. Res. Commun.* **533(3)**, 510-518 (2020).

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, 12/12/2022

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