

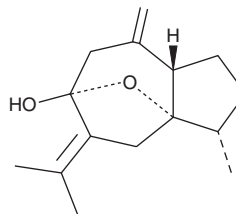
PRODUCT INFORMATION



Isocurcumenol

Item No. 30710

CAS Registry No.: 24063-71-6
Formal Name: octahydro-3S-methyl-8aS-methylene-5-(1-methylethylidene)-6H-3aS,6-epoxyazulen-6R-ol
Synonym: (+)-Isocurcumenol
MF: C₁₅H₂₂O₂
FW: 234.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Curcuma phaeocaulis*



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

Laboratory Procedures

Isocurcumenol is supplied as a solid. A stock solution may be made by dissolving the isocurcumenol in the solvent of choice, which should be purged with an inert gas. Isocurcumenol is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of isocurcumenol in these solvents is approximately 1 mg/ml. Isocurcumenol is also slightly soluble in ethanol.

Description

Isocurcumenol is a sesquiterpene that has been found in *C. zedoaria* rhizomes and has diverse biological activities.¹⁻³ Isocurcumenol inhibits LPS-induced production of nitric oxide (NO) in isolated mouse peritoneal macrophages.³ It reduces LPS/D-galactosamine-induced increases in serum levels of glutamic oxaloacetic transaminase (GOT) and glutamic pyruvic transaminase (GPT) in a mouse model of liver injury when administered at a dose of 50 mg/kg

References

1. Chen, C., Long, L., Zhang, F., *et al.* Antifungal activity, main active components and mechanism of *Curcuma longa* extract against *Fusarium graminearum*. *PLoS One* **13(3)**, e0194284 (2018).
2. Lakshmi, S., Padmaja, G., and Remani, P. Antitumour effects of isocurcumenol isolated from *Curcuma zedoaria* rhizomes on human and murine cancer cells. *Int. J. Med. Chem.* **253962** (2011).
3. Matsuda, H., Ninomiya, K., Morikawa, T., *et al.* Inhibitory effect and action mechanism of sesquiterpenes from zedoariae rhizoma on D-galactosamine/lipopolysaccharide-induced liver injury. *Bioorg. Med. Chem. Lett.* **8(4)**, 339-344 (1998).

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, 08/29/2023

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