

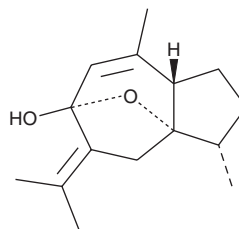
PRODUCT INFORMATION



Curcumenol

Item No. 26715

CAS Registry No.: 19431-84-6
Formal Name: (3S,3aS,6R,8aS)-1,2,3,4,5,8a-hexahydro-3,8-dimethyl-5-(1-methylethylidene)-6H-3a,6-epoxyazulen-6-ol
Synonym: (+)-Curcumenol
MF: C₁₅H₂₂O₂
FW: 234.3
Purity: ≥98%
UV/Vis.: λ_{max}: 262 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Curcuma* Rhizoma



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

Laboratory Procedures

Curcumenol is supplied as a crystalline solid. A stock solution may be made by dissolving the curcumenol in the solvent of choice. Curcumenol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of curcumenol in these solvents is approximately 3, 5, and 10 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of curcumenol can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of curcumenol in PBS, pH 7.2, is approximately 0.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Curcumenol is a sesquiterpene that has been found in *Curcuma* species.^{1,2} It is active against methicillin-resistant *S. aureus* (MRSA) and *P. aeruginosa* and cytotoxic to CEM-SS lymphoblastic leukemia cells (IC₅₀ = 11.9 μg/ml).¹ It also decreases cell viability of AGS human gastric carcinoma cells (IC₅₀ = 263.34 μM).²

References

- Aspollah Sukari, M., Wah, T.S., Saad, S.M., et al. Bioactive sesquiterpenes from *Curcuma ochrorhiza* and *Curcuma heyneana*. *Nat. Prod. Res.* **24**(9), 838-845 (2010).
- Lee, T.K., Lee, D., Lee, S.R., et al. Sesquiterpenes from *Curcuma zedoaria* rhizomes and their cytotoxicity against human gastric cancer AGS cells. *Bioorg. Chem.* **87**, 117-122 (2019).

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, 12/20/2022

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