

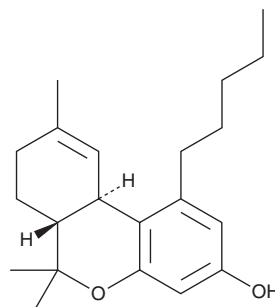
PRODUCT INFORMATION



Dihydrocurcumin

Item No. 39823

CAS Registry No.: 76474-56-1
Formal Name: (4Z,6E)-5-hydroxy-1,7-bis(4-hydroxy-3-methoxyphenyl)-4,6-heptadien-3-one
Synonym: Letestuianin B
MF: C₂₁H₂₂O₆
FW: 370.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/Zingiber officinale Rose



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

Laboratory Procedures

Dihydrocurcumin is supplied as a solid. A stock solution may be made by dissolving the dihydrocurcumin in the solvent of choice, which should be purged with an inert gas. Dihydrocurcumin is soluble in acetone, chloroform, dichloromethane, DMSO, and ethyl acetate.

Description

Dihydrocurcumin is an active metabolite of curcumin (Item Nos. 81025 | 81025.1).¹ It is formed from curcumin via reductases. It reduces triglyceride levels and increases glucose uptake in HepG2 and L-02 cellular models of steatosis induced by oleic acid (Item Nos. 90260 | 24659) when used at a concentration of 20 μM.²

References

1. Pandey, A., Chaturvedi, M., Mishra, S., *et al.* Reductive metabolites of curcumin and their therapeutic effects. *Heliyon* **6(11)**, (2020).
2. Yu, Q., Liu, Y., Wu, Y., *et al.* Dihydrocurcumin ameliorates the lipid accumulation, oxidative stress and insulin resistance in oleic acid-induced L02 and HepG2 cells. *Biomed. Pharmacother.* **103**, 1327-1336 (2018).

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, 11/03/2023

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