

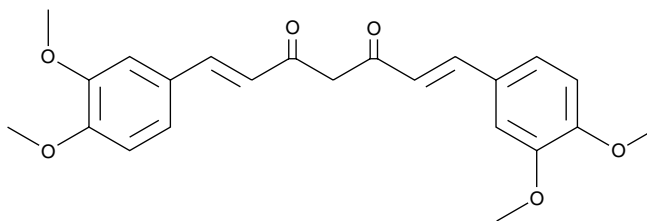
Product Information



Dimethoxycurcumin

Item No. 10009986

CAS Registry No.: 160096-59-3
Formal Name: 1,7-bis(3,4-dimethoxyphenyl)-1,6-heptadiene-3,5-dione
MF: C₂₃H₂₄O₆
FW: 396.4
Purity: ≥90%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that dimethoxycurcumin be stored as supplied at -20°C. It will be stable for at least two years.

Dimethoxycurcumin is supplied as a crystalline solid. A stock solution may be made by dissolving the dimethoxy curcumin in an organic solvent purged with an inert gas. Dimethoxycurcumin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of dimethoxy curcumin in these solvents is approximately 0.25, 5, and 10 mg/ml respectively.

Dimethoxycurcumin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, dimethoxycurcumin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Dimethoxycurcumin has a solubility of approximately 0.30 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.

Naturally occurring phytochemicals such as turmeric (curcumin) have been found to inhibit the growth of tumor cells. Dimethoxycurcumin is an analog of curcumin obtained by methylation of both free phenolic groups in the parent compound. A 30-fold increase in potency was observed in the growth suppression of HCT116 tumor cells following this modification.¹

Reference

1. Ohori, H., Yamakoshi, H., Tomizawa, M., *et al.* Synthesis and biological analysis of new curcumin analogues bearing an enhanced potential for the medicinal treatment of cancer. *Mol. Cancer Therapeutics* **5(10)**, 2563-2571 (2006).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10009986

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 10/14/2011

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com