

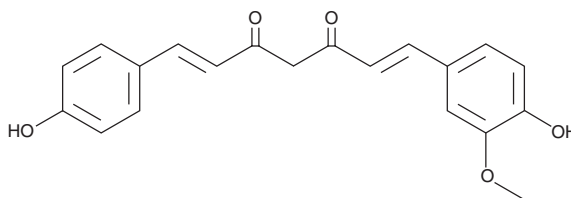
PRODUCT INFORMATION



Demethoxycurcumin

Item No. 10961

CAS Registry No.: 22608-11-3
Formal Name: 1-(4-hydroxy-3-methoxyphenyl)-
7-(4-hydroxyphenyl)-1E,6E-
heptadiene-3,5-dione
Synonym: DMC
MF: C₂₀H₁₈O₅
FW: 338.4
Purity: ≥98%
UV/Vis.: λ_{max}: 254, 421 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Demethoxycurcumin is supplied as a crystalline solid. A stock solution may be made by dissolving the demethoxycurcumin in the solvent of choice, which should be purged with an inert gas. Demethoxycurcumin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of demethoxycurcumin in these solvents is approximately 0.5, 10, and 14 mg/ml, respectively.

Description

Demethoxycurcumin (DMC) is a natural demethoxy derivative of curcumin. More stable than curcumin in physiological media, DMC suppresses proliferation in cancer cells at 50-100 μM.¹ It down-regulates the transcriptional coactivator p300, suppressing the Wnt/β-catenin pathway, and inhibits lipopolysaccharide induction of iNOS.²

References

1. Sandur, S.K., Pandley, M.K., Ahn, K.S., *et al.* Curcumin, demethoxycurcumin, bisdemethoxycurcumin, tetrahydrocurcumin and turmerones differentially regulate anti-inflammatory and anti-proliferative responses through a ROS-independent mechanism. *Carcinogenesis* **28(8)**, 1765-1773 (2007).
2. Ryu, M.J., Cho, M., Song, J.Y., *et al.* Natural derivatives of curcumin attenuate the Wnt/β-catenin pathway through down-regulation of the transcriptional coactivator p300. *Biochem. Biophys. Res. Commun.* **377(4)**, 1304-1308 (2008).

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, 02/21/2024

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