

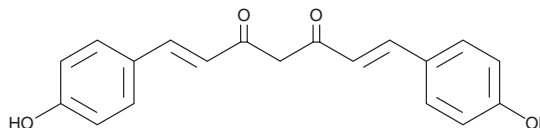
PRODUCT INFORMATION



Bisdemethoxycurcumin

Item No. 10960

CAS Registry No.: 33171-05-0
Formal Name: 1,7-bis(4-hydroxyphenyl)-1E,6E-heptadiene-3,5-dione
Synonym: BDMC
MF: C₁₉H₁₆O₄
FW: 308.3
Purity: ≥98%
UV/Vis.: λ_{max}: 419 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Bisdemethoxycurcumin (BDMC) is supplied as a crystalline solid. A stock solution may be made by dissolving the BDMC in the solvent of choice. BDMC is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of BDMC in these solvents is approximately 5, 10, and 14 mg/ml, respectively.

Description

BDMC is a natural demethoxy derivative of curcumin. It is a potent activator of macrophage phagocytosis, interacting with 1 α ,25-dihydroxy vitamin D₃ to stimulate amyloid- β clearance by macrophages (optimal stimulation at 100 nM BDMC).¹⁻³ More stable than curcumin in physiological media, BDMC suppresses proliferation in cancer cells.^{4,5} It down-regulates the transcriptional coactivator p300, suppressing the Wnt/ β -catenin pathway, and inhibits LPS induction of iNOS expression.^{6,7}

References

1. Fiala, M., Liu, P.T., Espinosa-Jeffrey, A., et al. *Proc. Natl. Acad. Sci. USA* **104**(31), 12849-12854 (2007).
2. Masoumi, A., Goldenson, B., Ghirmai, S., et al. *J. Alzheimers Dis.* **17**(3), 703-717 (2009).
3. Cashman, J.R., Ghirmai, S., Abel, K.J., et al. *BMC Neurosci.* **9**, (2008).
4. Sandur, S.K., Pandley, M.K., Ahn, K.S., et al. *Carcinogenesis* **28**(8), 1765-1773 (2007).
5. Basile, V., Ferrari, E., Lazzari, S., et al. *Biochem. Pharmacol.* **78**(10), 1305-1315 (2009).
6. Ryu, M.J., Cho, M., Song, J.Y., et al. *Biochem. Biophys. Res. Commun.* **377**(4), 1304-1308 (2008).
7. Mahattanadul, S., Nakamura, T., Panichayupakaranant, P., et al. *Phytomedicine* **16**(4), 342-351 (2009).

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, 03/28/2018

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