

PRODUCT INFORMATION



Garcinol

Item No. 10566

CAS Registry No.: 78824-30-3
Formal Name: (1R,5R,7R)-3-(3,4-dihydroxybenzoyl)-4-hydroxy-8,8-dimethyl-1,7-bis(3-methyl-2-buten-1-yl)-5-[(2S)-5-methyl-2-(1-methylethenyl)-4-hexen-1-yl]-bicyclo[3.3.1]non-3-ene-2,9-dione

Synonyms: Camboginol, Guttiferone F

MF: C₃₈H₅₀O₆

FW: 602.8

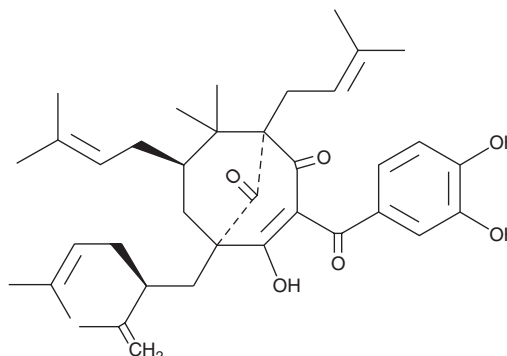
Purity: ≥95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years

Item Origin: Plant/*Garcinia indica*



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

Laboratory Procedures

Garcinol is supplied as a crystalline solid. A stock solution may be made by dissolving the garcinol in the solvent of choice, which should be purged with an inert gas. Garcinol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of garcinol in ethanol is approximately 20 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Garcinol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, garcinol should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Garcinol has a solubility of approximately 0.1 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Garcinol is a polyisoprenylated benzophenone that has been found in *G. indica* and has diverse biological activities.¹⁻⁵ It inhibits the histone acetyltransferases (HATs) p300 and p300/CREB-binding protein-associated factor (PCAF; IC₅₀s = ~7 and ~5 μM, respectively), as well as COX-1, microsomal prostaglandin E₂ (PGE₂) synthase-1 (mPGES-1) and 5-lipoxygenase (5-LO; IC₅₀s = 12, 0.1, and 0.3 μM, respectively).^{1,2} Garcinol (0.8 μM) is active against the fungus *C. neoformans in vitro*.³ It scavenges hydroxyl radicals and superoxide anions in cell-free assays.⁴ Garcinol (5 μM) induces neurite outgrowth of primary embryonic rat neural progenitor cells.⁵ It reduces ulcer area in rat models of gastric ulcers induced by water immersion stress or indomethacin (Item No. 70270) when administered at a dose of 200 mg/kg.⁴

References

1. Balasubramanyam, K., Altaf, M., Varier, R.A., et al. *J. Biol. Chem.* **279**(32), 33716-33726 (2004).
2. Koeberle, A., Northoff, H., and Werz, O. *Biochem. Pharmacol.* **77**(9), 1513-1521 (2009).
3. O'Meara, T.R., Hay, C., Price, M.S., et al. *Eukaryot. Cell* **9**(8), 1193-1202 (2010).
4. Yamaguchi, F., Saito, M., Ariga, T., et al. *J. Agric. Food Chem.* **48**(6), 2320-2325 (2000).
5. Weng, M.-S., Liao, C.-H., Yu, S.-Y., et al. *J. Agric. Food Chem.* **59**(3), 1031-1040 (2011).

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, 04/06/2023

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