

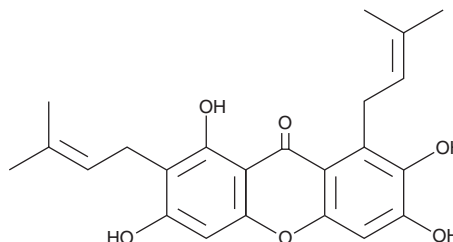
PRODUCT INFORMATION



γ -Mangostin

Item No. 26675

CAS Registry No.: 31271-07-5
Formal Name: 1,3,6,7-tetrahydroxy-2,8-bis(3-methyl-2-buten-1-yl)-9H-xanthen-9-one
MF: $C_{23}H_{24}O_6$
FW: 396.4
Purity: $\geq 98\%$
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/Mangosteen skin



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

Laboratory Procedures

γ -Mangostin is supplied as a crystalline solid. A stock solution may be made by dissolving the γ -mangostin in the solvent of choice. γ -Mangostin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of γ -mangostin in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

γ -Mangostin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, γ -mangostin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. γ -Mangostin has a solubility of approximately 0.02 mg/ml in a 1:40 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

γ -Mangostin is a xanthone that has been found in *G. mangostana* and has diverse biological activities.¹⁻⁴ It scavenges 2,2-diphenyl-1-picrylhydrazyl (DPPH; Item No. 14805) radicals *in vitro* with an IC_{50} value of 23.6 μM .¹ γ -Mangostin is cytotoxic to HL-60, SMMC-7721, A549, MCF-7, and SW480 cancer cells and BEAS-2B non-cancerous pulmonary epithelial cells *in vitro* (IC_{50}s = 7.39, 6.57, 10.07, 5.33, 8.4, and 7.43 μM , respectively).² It inhibits LPS-induced expression of IL-6, IL-10, and TNF- α in human macrophages in a concentration-dependent manner.³ Macrophage-conditioned media from macrophages pre-incubated with γ -mangostin (30 μM) prior to stimulation with LPS has a reduced ability to induce inflammation and insulin resistance in primary human adipocytes. γ -Mangostin (5 and 10 mg/kg) inhibits carbon tetrachloride-induced increases in serum levels of aspartate transaminase (AST) and alanine aminotransferase (ALT) and decreases in hepatic superoxide dismutase 2 (SOD2) activity and glutathione (GSH) content in a mouse model of liver fibrosis.⁴

References

1. Li, X. Comparative study of 1,1-diphenyl-2-picryl-hydrazyl radical (DPPH \bullet) scavenging capacity of the antioxidant xanthenes family. *ChemistrySelect* **3**(46), 13081-13086 (2018).
2. Chi, X.-Q., Zi, C.-T., Li, H.-M., *et al.* Design, synthesis and structure-activity relationships of mangostin analogs as cytotoxic agents. *RSC Adv.* **8**(72), 41377-41388 (2018).
3. Bumrungpert, A., Kalpravidh, R.W., Chuang, C.-C., *et al.* Xanthenes from mangosteen inhibit inflammation in human macrophages and in human adipocytes exposed to macrophage-conditioned media. *J. Nutr.* **140**(4), 842-847 (2010).
4. Wang, A., Zhou, F., Li, D., *et al.* γ -Mangostin alleviates liver fibrosis through Sirtuin 3-superoxide-high mobility group box 1 signaling axis. *Toxicol. Appl. Pharmacol.* **363**, 142-153 (2019).

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, 09/29/2022

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