

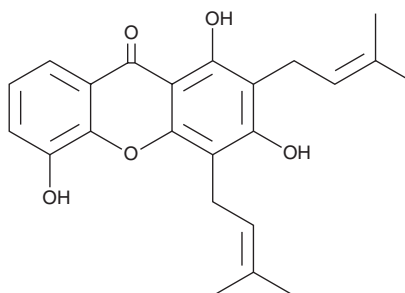
PRODUCT INFORMATION



8-Desoxygartanin

Item No. 40275

CAS Registry No.: 33390-41-9
Formal Name: 1,3,5-trihydroxy-2,4-bis(3-methyl-2-buten-1-yl)-9H-xanthen-9-one
Synonyms: 8-Deoxygartanin, NSC 692944
MF: C₂₃H₂₄O₅
FW: 380.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Garcinia mangostana* L.



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

Laboratory Procedures

8-Desoxygartanin is supplied as a solid. A stock solution may be made by dissolving the 8-desoxygartanin in the solvent of choice, which should be purged with an inert gas. 8-Desoxygartanin is soluble in the organic solvent acetonitrile.

Description

8-Desoxygartanin is a xanthone that has been found in *G. mangostana* and has anti-inflammatory and enzyme inhibitory activities.^{1,2} It inhibits NF-κB p65 binding to DNA in a cell-free assay (IC₅₀ = 11.3 μM).¹ 8-Desoxygartanin inhibits butyrylcholinesterase (BChE) and acetylcholinesterase (AChE; IC₅₀s = 2.46 and 7.75 μg/ml).²

References

1. Han, A.-R., Kim, J.-A., Lantvit, D.D., *et al.* Cytotoxic xanthone constituents of the stem bark of *Garcinia mangostana* (mangosteen). *J. Nat. Prod.* **72**(11), 2028-2031 (2009).
2. Khaw, K.Y., Choi, S.B., Tan, S.C., *et al.* Prenylated xanthenes from mangosteen as promising cholinesterase inhibitors and their molecular docking studies. *Phytomedicine* **21**(11), 1303-1309 (2014).

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, 02/19/2024

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