

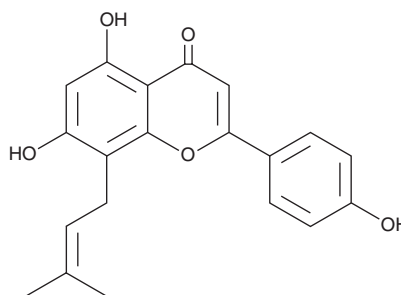
PRODUCT INFORMATION



Licoflavone C

Item No. 35031

CAS Registry No.: 72357-31-4
Formal Name: 5,7-dihydroxy-2-(4-hydroxyphenyl)-8-(3-methyl-2-buten-1-yl)-4H-1-benzopyran-4-one
Synonym: 8-Prenylapigenin
MF: C₂₀H₁₈O₅
FW: 338.4
Purity: ≥98%
UV/Vis.: λ_{max}: 274, 328 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Glycyrrhiza inflata*



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

Laboratory Procedures

Licoflavone C is supplied as a solid. A stock solution may be made by dissolving the licoflavone C in the solvent of choice, which should be purged with an inert gas. Licoflavone C is soluble in the organic solvent ethanol.

Description

Licoflavone C is a prenylated flavone that has been found in *G. ephedroides* and has diverse biological activities.¹⁻³ It induces estrogen receptor-dependent gene expression in *S. cerevisiae* expressing human estrogen receptor α (ER α) in a reporter assay when used at a concentration of 0.1 μ M.¹ Licoflavone C inhibits LPS-induced production of nitric oxide (NO) in RAW 264.7 macrophages (IC₅₀ = 20.4 μ M).² It is active against *E. coli*, *P. aeruginosa*, and various strains of *Candida* (MICs = 7.81-15.62 μ g/ml) and is cytotoxic to HepG2 cells (IC₅₀ = 9 μ g/ml).³

References

1. Garritano, S., Pinto, B., Giachi, I., *et al.* Assessment of estrogenic activity of flavonoids from Mediterranean plants using an *in vitro* short-term test. *Phytomedicine* **12(1-2)**, 143-147 (2005).
2. Han, A.-R., Kang, Y.-J., Windono, T., *et al.* Prenylated flavonoids from the heartwood of *Artocarpus communis* with inhibitory activity on lipopolysaccharide-induced nitric oxide production. *J. Nat. Prod.* **69(4)**, 719-721 (2006).
3. Edziri, H., Mastouri, M., Mahjoub, M.A., *et al.* Antibacterial, antifungal and cytotoxic activities of two flavonoids from *Retama raetam* flowers. *Molecules* **17(6)**, 7284-7293 (2012).

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, 10/19/2022

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