

PRODUCT INFORMATION



Epimedin C

Item No. 34375

CAS Registry No.: 110642-44-9
Formal Name: 3-[[6-deoxy-2-O-(6-deoxy- α -L-mannopyranosyl)- α -L-mannopyranosyl]oxy]-7-(β -D-glucopyranosyloxy)-5-hydroxy-2-(4-methoxyphenyl)-8-(3-methyl-2-buten-1-yl)-4H-1-benzopyran-4-one

Synonym: Baohuoside VI

MF: C₃₉H₅₀O₁₉

FW: 822.8

Purity: \geq 98%

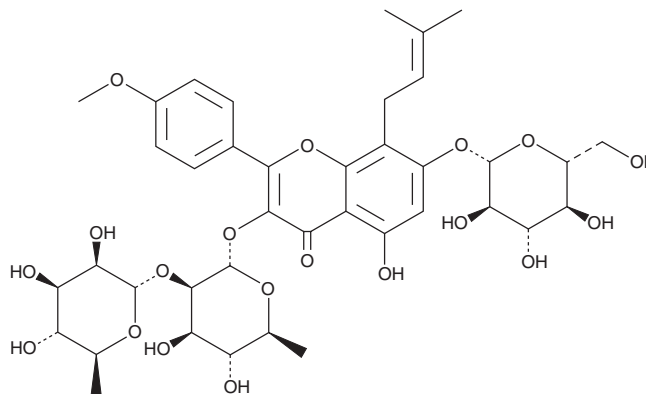
UV/Vis.: λ_{max} : 272 nm

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years

Item Origin: Plant/*Epimedii folium*



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

Laboratory Procedures

Epimedin C is supplied as a solid. A stock solution may be made by dissolving the epimedin C in the solvent of choice, which should be purged with an inert gas. Epimedin C is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of epimedin C in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of epimedin C can be prepared by directly dissolving the solid in aqueous buffers. The solubility of epimedin C in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Epimedin C is a flavonoid glycoside that has been found in *E. koreanum* and has diverse biological activities.¹⁻³ It scavenges DPPH (Item No. 14805) radicals in a cell-free assay (IC_{50} = 142 μ M).¹ Epimedin B increases proliferation of UMR-106 osteoblasts by 30.4 and 41.6% when used at concentrations of 0.1 and 10 μ M, respectively.² It inhibits ear edema induced by phorbol 12-myristate 13-acetate (TPA; Item No. 10008014) in mice (ED_{50} = 192 nmol/ear).³

References

1. Kim, E.S., Kim, M.K., Kang, H.K., *et al.* Flavonol glycosides with antioxidant activity from the aerial parts of *Epimedium koreanum* nakai. *Nat. Prod. Sci.* **14**(4), 233-238 (2008).
2. Meng, F.-H., M., Li, Y.-B., Xiong, A.-L., *et al.* Osteoblastic proliferative activity of *Epimedium brevicornum* maxim. *Phytomedicine* **12**(3), 189-93 (2005).
3. Yasukawa, K., Ko, S.-K., and Whang, W.-K. Inhibitory effects of the aerial parts of *Epimedium koreanum* on TPA-induced inflammation and tumour promotion in two-stage carcinogenesis in mouse skin. *J. Pharm. Nutr. Sci.* **6**(2), 38-42 (2016).

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, 09/30/2022

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