

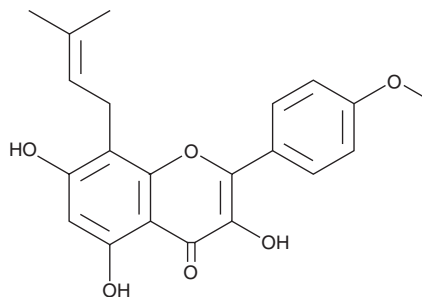
PRODUCT INFORMATION



Icaritin

Item No. 20236

CAS Registry No.: 118525-40-9
Formal Name: 3,5,7-trihydroxy-2-(4-methoxyphenyl)-8-(3-methyl-2-buten-1-yl)-4H-1-benzopyran-4-one
MF: C₂₁H₂₀O₆
FW: 368.4
Purity: ≥98%
UV/Vis.: λ_{max}: 273, 318, 374 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Icaritin is supplied as a crystalline solid. A stock solution may be made by dissolving the icaritin in the solvent of choice. Icaritin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of icaritin in these solvents is approximately 0.3, 5, and 3 mg/ml, respectively.

Icaritin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, icaritin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Icaritin has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Icaritin is a flavonoid first isolated from the Chinese herb *H. epimedii* that demonstrates anticancer activity against a variety of tumor cell lines.¹ It has been shown to inhibit fatty acid synthase, reducing IGF-1-induced activation of STAT3 in several melanoma cell lines.¹

Reference

1. Wu, J., Du, J., Fu, X., et al. Icaritin, a novel FASN inhibitor, exerts anti-melanoma activities through IGF-1R/STAT3 signaling. *Oncotarget* (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, 11/04/2022

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