

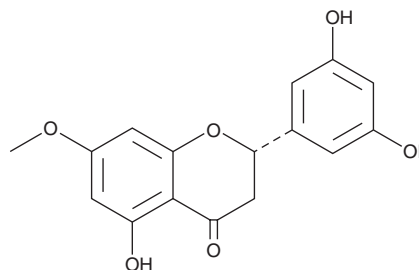
PRODUCT INFORMATION



Blumeatin

Item No. 34809

CAS Registry No.: 118024-26-3
Formal Name: 2S-(3,5-dihydroxyphenyl)-2,3-dihydro-5-hydroxy-7-methoxy-4H-1-benzopyran-4-one
MF: C₁₆H₁₄O₆
FW: 302.3
Purity: ≥98%
UV/Vis.: λ_{max}: 288 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Blumea balsamifera*



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

Laboratory Procedures

Blumeatin is supplied as a solid. A stock solution may be made by dissolving the blumeatin in the solvent of choice, which should be purged with an inert gas. Blumeatin is soluble in DMSO.

Description

Blumeatin is a flavonoid that has been found in *Blumea* and has antioxidant and anticancer activities.¹⁻³ It scavenges DPPH (Item No. 14805) radicals in a cell-free assay with a 50% scavenging capacity (SC₅₀) value of 90.8 µg/ml.² Blumeatin (10 and 100 µM) decreases carbon tetrachloride-induced lipid peroxidation in isolated rhesus monkey hepatocytes.¹ It is also cytotoxic to KB oral carcinoma cells (IC₅₀ = 47.72 µg/ml).³

References

1. Zhao, J.H., Xu, S.-B., Wang, Z.-L., *et al.* Protective actions of Blumea flavanones on primary cultured hepatocytes and liver subcellular organelle against lipid peroxidation. *J. Chin. Pharma. Sci.* **7(3)**, 152-156 (1998).
2. Nessa, F., Ismail, Z., Mohamed, N., *et al.* Free radical-scavenging activity of organic extracts and of pure flavonoids of *Blumea balsamifera* DC leaves. *Food Chem.* **88(2)**, 243-252 (2004).
3. Saewan, N., Koysoomboon, S., and Chantrapromma, K. Anti-tyrosinase and anti-cancer activities of flavonoids from *Blumea balsamifera* DC. *J. Med. Plant Res.* **5(6)**, 1018-1025 (2011).

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, 12/12/2022

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