

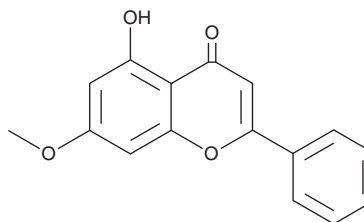
PRODUCT INFORMATION



Techtochrysin

Item No. 35791

CAS Registry No.: 520-28-5
Formal Name: 5-hydroxy-7-methoxy-2-phenyl-4H-1-benzopyran-4-one
Synonyms: 5-Hydroxy-7-methoxyflavone, NSC 80687, Techtochrysin
MF: C₁₆H₁₂O₄
FW: 268.3
Purity: ≥98%
UV/Vis.: λ_{max}: 269 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Pinus strobus*



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

Laboratory Procedures

Techtochrysin is supplied as a solid. A stock solution may be made by dissolving the techtochrysin in the solvent of choice, which should be purged with an inert gas. Techtochrysin is soluble in the organic solvent dimethyl formamide at a concentration of approximately 1 mg/ml. Techtochrysin is also slightly soluble in DMSO.

Description

Techtochrysin is a flavonoid that has been found in *T. indica* and has diverse biological activities.¹⁻³ It induces adipogenesis and glucose uptake in 3T3-L1 adipocytes when used at concentrations of 12.5, 25, and 50 µg/ml.¹ Techtochrysin is active against *S. aureus*, *L. monocytogenes*, *B. subtilis*, *S. enterica*, *E. coli*, *V. cholera*, *K. pneumoniae*, *S. dysenteriae*, and *P. aeruginosa* (MICs = 24-64 µg/ml).² It also scavenges peroxy radicals in a cell-free assay when used at concentrations of 1 and 10 µM.³

References

1. Hasan, M.M., Ahmed, Q.U., Soad, S.Z.M., *et al.* Flavonoids from *Tetracera indica* Merr. induce adipogenesis and exert glucose uptake activities in 3T3-L1 adipocyte cells. *BMC Complement. Altern. Med.* **17(1)**, 431 (2017).
2. Rajendran, N., Subramaniam, S., Christena, L.R., *et al.* Antimicrobial flavonoids isolated from Indian medicinal plant *Scutellaria oblonga* inhibit biofilms formed by common food pathogens. *Nat. Prod. Res.* **30(17)**, 2002-2006 (2016).
3. Thao, N.P., Luyen, B.T.T., Lee, S.H., *et al.* Anti-osteoporotic and antioxidant activities by rhizomes of *Kaempferia parviflora* Wall. ex Baker. *Nat. Prod. Sci.* **22(1)**, 13-19 (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, 12/02/2022

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