

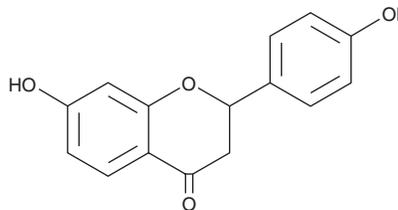
PRODUCT INFORMATION



(±)-Liquiritigenin

Item No. 28823

CAS Registry No.: 69097-97-8
Formal Name: 2,3-dihydro-7-hydroxy-2-(4-hydroxyphenyl)-4H-1-benzopyran-4-one
MF: C₁₅H₁₂O₄
FW: 256.3
Purity: ≥98%
UV/Vis.: λ_{max}: 215, 232, 276, 311 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Glycyrrhiza glabra*



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

Laboratory Procedures

(±)-Liquiritigenin is supplied as a solid. A stock solution may be made by dissolving the (±)-liquiritigenin in the solvent of choice, which should be purged with an inert gas. (±)-Liquiritigenin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (±)-liquiritigenin in ethanol is approximately 25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

(±)-Liquiritigenin is a flavonoid that has been found in *P. floribundum* and has anticancer activity.¹ It is cytotoxic to HL-60, HCT8, and B16 cancer cell lines (IC₅₀s = 22.5, 19.5, and 5 µg/ml, respectively).

Reference

1. Falcão, M.J., Pouliquem, Y.B., Lima, M.A., et al. Cytotoxic flavonoids from *Platymiscium floribundum*. *J. Nat. Prod.* **68**(3), 423-426 (2005).

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, 12/22/2022

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