

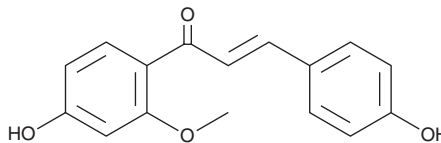
PRODUCT INFORMATION



2'-O-Methylisoliquiritigenin

Item No. 36490

CAS Registry No.: 51828-10-5
Formal Name: 1-(4-hydroxy-2-methoxyphenyl)-3-(4-hydroxyphenyl)-2-propen-1-one
MF: $C_{16}H_{14}O_4$
FW: 270.3
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Caesalpinia sinensis*



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

Laboratory Procedures

2'-O-Methylisoliquiritigenin is supplied as a solid. A stock solution may be made by dissolving the 2'-O-methylisoliquiritigenin in the solvent of choice, which should be purged with an inert gas. 2'-O-Methylisoliquiritigenin is sparingly soluble (1-10 mg/ml) in methanol and DMSO.

Description

2'-O-Methylisoliquiritigenin is a chalcone that has been found in *C. sappan* and has diverse biological activities.^{1,2} It reduces colony formation and induces apoptosis and cell cycle arrest at the G₂/M phase in HCC827 non-small cell lung cancer (NSCLC) cells when used at a concentration of 12 μM .¹ 2'-O-Methylisoliquiritigenin induces reactive oxygen species (ROS) production and mitochondrial membrane depolarization in HCC827 cells. It inhibits UV radiation-induced increases in dorsal skin thickness in mice when applied topically at doses of 0.5 or 1 $\mu\text{mol}/\text{mouse}$.² 2'-O-Methylisoliquiritigenin (10 or 20 mg/kg) decreases tumor volume and weight in a SK-MEL-2 melanoma mouse xenograft model.

References

1. Lee, J.-Y., Lee, S.-O., Kwak, A.-W., *et al.* 3-Deoxysappanchalcone inhibits cell growth of gefitinib-resistant lung cancer cells by simultaneous targeting of EGFR and MET kinases. *Biomol. Ther. (Seoul)* **31(4)**, 446-455 (2023).
2. Fu, X., Zhao, R., Yoon, G., *et al.* 3-Deoxysappanchalcone inhibits skin cancer proliferation by regulating T-lymphokine-activated killer cell-originated protein kinase *in vitro* and *in vivo*. *Front. Cell Dev. Biol.* **9:638174**, (2021).

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, 02/05/2025

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