

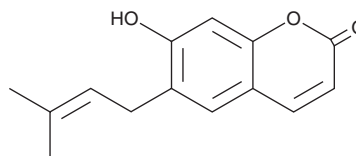
PRODUCT INFORMATION



7-Demethylsuberosin

Item No. 35282

CAS Registry No.: 21422-04-8
Formal Name: 7-hydroxy-6-(3-methyl-2-buten-1-yl)-2H-1-benzopyran-2-one
Synonym: Demethylsuberosin
MF: $C_{14}H_{14}O_3$
FW: 230.3
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 333 nm
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years
Item Origin: Plant/*Angelica dahurica*



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

Laboratory Procedures

7-Demethylsuberosin is supplied as a solid. A stock solution may be made by dissolving the 7-demethylsuberosin in the solvent of choice, which should be purged with an inert gas. 7-Demethylsuberosin is soluble in the organic solvent DMSO at a concentration of approximately 250 mg/ml.

Description

7-Demethylsuberosin is a prenylated coumarin that has been found in *Angelica tenuissima* and has anti-inflammatory and neuroprotective activities.^{1,2} It inhibits histamine release induced by phorbol 12-myristate 13-acetate (TPA; Item No. 10008014) and A23187 (Item No. 11016) from HMC-1 human mast cell leukemia cells when used at a concentration of 10 μM .¹ 7-Demethylsuberosin (0.1-10 μM) protects against glutamate-induced cytotoxicity in primary rat cortical cells.²

References

1. Choi, H.G., Je, I.-G., Kim, G.J., *et al.* Chemical constituents of the root of *Angelica tenuissima* and their anti-allergic inflammatory activity. *Nat. Prod. Commun.* **12**(5), 779-780 (2017).
2. Kang, S.Y. and Kim, Y.C. Neuroprotective coumarins from the root of *Angelica gigas*: Structure-activity relationships. *Arch. Pharm. Res.* **30**(11), 1368-1373 (2007).

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, 03/17/2025

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