

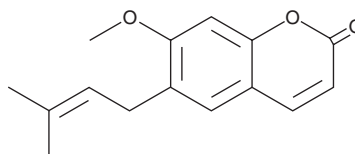
PRODUCT INFORMATION



Suberosin

Item No. 35035

CAS Registry No.: 581-31-7
Formal Name: 7-methoxy-6-(3-methyl-2-buten-1-yl)-2H-1-benzopyran-2-one
Synonym: NSC 31869
MF: $C_{15}H_{16}O_3$
FW: 244.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 223, 331 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Angelica pubescens*



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

Laboratory Procedures

Suberosin is supplied as a solid. A stock solution may be made by dissolving the suberosin in the solvent of choice, which should be purged with an inert gas. Suberosin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of suberosin in DMSO is approximately 3 mg/ml and approximately 10 mg/ml in ethanol and DMF.

Description

Suberosin is a prenylated coumarin that has been found in *C. articulata* and has diverse biological activities.¹⁻³ It inhibits platelet aggregation induced by arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607), collagen, platelet-activating factor (PAF), adenosine 5'-diphosphate (ADP; Item No. 21121), or U-46619 (Item No. 16450) in washed isolated rabbit platelets (IC_{50} s = 60, 132, 200, 234, and 200 μM , respectively).² Suberosin (50 or 100 μM) reduces arachidonic acid-induced formation of thromboxane B_2 (TXB₂; Item No. 19030) in the same model. It inhibits NF- κ B nuclear translocation induced by phytohemagglutinin (PHA) in, as well as PHA-induced proliferation of, human peripheral blood mononuclear cells (PBMCs) when used at a concentration of 100 μM .³

References

1. Paudel, S., Kim, Y., Choi, S.M., *et al.* Identification of suberosin metabolites in human liver microsomes by high-performance liquid chromatography combined with high-resolution quadrupole-orbitrap mass spectrometer. *J. Mass Spectrom.* **56**(4), e4623 (2021).
2. Teng, C.M., Li, H.L., Wu, T.S., *et al.* Antiplatelet actions of some coumarin compounds isolated from plant sources. *Thromb. Res.* **66**(5), 549-557 (1992).
3. Chen, Y.-C., Tsai, W.-J., Wu, M.-H., *et al.* Suberosin inhibits proliferation of human peripheral blood mononuclear cells through the modulation of the transcription factors NF-AT and NF- κ B. *Br. J. Pharmacol.* **150**(3), 298-312 (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, 11/21/2022

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