

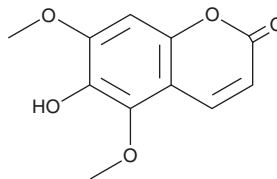
PRODUCT INFORMATION



Fraxinol

Item No. 28218

CAS Registry No.: 486-28-2
Formal Name: 6-hydroxy-5,7-dimethoxy-2H-1-benzopyran-2-one
MF: C₁₁H₁₀O₅
FW: 222.2
Purity: ≥98%
UV/Vis.: λ_{max}: 233, 310 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Fraxinus stylosa* Lingelsh.



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

Laboratory Procedures

Fraxinol is supplied as a crystalline solid. A stock solution may be made by dissolving the fraxinol in the solvent of choice, which should be purged with an inert gas. Fraxinol is soluble in the organic solvent DMSO.

Description

Fraxinol is a coumarin originally isolated from Ash tree bark.¹ It inhibits growth of GLC-4 small cell lung carcinoma and COLO 320 colorectal cancer cells (IC₅₀s = 193 and 165 μM, respectively).² Fraxinol potentiates barbiturate-induced sleep in rats and rabbits.³

References

1. Spath, E. and Jerzmanowska-Sienkiewiczowa, Z. Natural coumarins. XXV. Fraxinol, a new constituent of the bark of the ash. *Berichte der Deutschen Chemischen Gesellschaft [Abteilung] B: Abhandlungen* **70B**, 698-702 (1937).
2. Kolodziej, H., Kayser, O., Woerdenbag, H.J., et al. Structure-cytotoxicity relationships of a series of natural and semi-synthetic simple coumarins as assessed in two human tumour cell lines. *Z. Naturforsch. C* **52(3-4)**, 240-244 (1997).
3. Khadzhai, Y.I. Effect of coumarin, flavonoid, and khellin derivatives on the soporific effect of barbiturates. *Farmakologiya i Toksikologiya (Kiev)* **7**, 46-48 (1972).

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, 10/18/2022

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