

PRODUCT INFORMATION

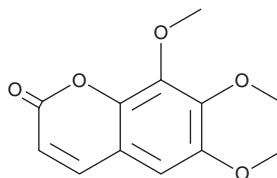


Dimethylfraxetin

Item No. 36485

CAS Registry No.: 6035-49-0
Formal Name: 6,7,8-trimethoxy-2H-1-benzopyran-2-one
Synonyms: Fraxetin dimethyl ether, 6,7,8-Trimethoxycoumarin

MF: $C_{12}H_{12}O_5$
FW: 236.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 227 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Fraxinus chinensis* Roxb. bark



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

Laboratory Procedures

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

Description

Dimethylfraxetin is a coumarin that has been found in *T. lucida* and has anti-inflammatory activity.¹ It reduces auricular edema induced by phorbol 12-myristate 13-acetate (TPA; Item No. 10008014) in mice when administered at a dose of 1 mg/ear.

Reference

1. Nayeli, M.-B., Maribel, H.-R., Enrique, J.-F., *et al.* Anti-inflammatory activity of coumarins isolated from *Tagetes lucida* Cav. *Nat. Prod. Res.* **34**(22), 3244-3248 (2020).

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, 12/19/2022

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