

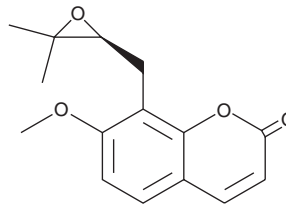
PRODUCT INFORMATION



Meranzin

Item No. 34743

CAS Registry No.: 23971-42-8
Formal Name: 8-[[[(2S)-3,3-dimethyl-2-oxiranyl]methyl]-7-methoxy-2H-1-benzopyran-2-one
MF: C₁₅H₁₆O₄
FW: 260.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/Citrus maxima peel



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

Laboratory Procedures

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

Description

Meranzin is a coumarin that has been found in *Murraya exotica* and has diverse biological activities.^{1,2} It is cytotoxic to NSCLC-N6 and KB cancer cells (IC₅₀s = 14.6 and 12.3 μg/ml, respectively).² Meranzin (10 mg/kg) increases the latency to paw withdrawal in the hot plate test and reduces xylene-induced ear edema in mice.¹ It decreases paw edema induced by carrageenan in rats when administered at the same dose.

References

1. Wu, L., Li, P., Wang, X., *et al.* Evaluation of anti-inflammatory and antinociceptive activities of *Murraya exotica*. *Pharm. Biol.* **48(12)**, 1344-1353 (2010).
2. Kofinas, C., Chinou, L., Loukis, A., *et al.* Cytotoxic coumarins from the aerial parts of *Tordylium apulum* and their effects on a non-small-cell bronchial carcinoma cell line. *Planta Med.* **64(2)**, 174-176 (1998).

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, 12/08/2022

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