

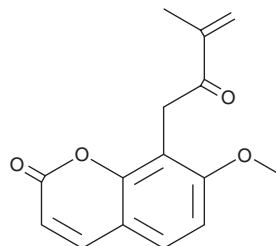
PRODUCT INFORMATION



Murrayone

Item No. 34767

CAS Registry No.: 19668-69-0
Formal Name: 7-methoxy-8-(3-methyl-2-oxo-3-buten-1-yl)-2H-1-benzopyran-2-one
MF: C₁₅H₁₄O₄
FW: 258.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Murraya paniculata*



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

Laboratory Procedures

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

Description

Murrayone is a coumarin that has been found in *L. japonicus* and has antiplatelet activity.^{1,2} It inhibits ADP-induced platelet aggregation *in vitro*.¹

References

1. Yang, H., Zhou, Q.-m., Peng, C., *et al.* Coumarins from *Leonurus japonicus* and their anti-platelet aggregative activity. *Zhongguo Zhong Yao Za Zhi* **39(22)**, 4356-4359 (2014).
2. Song, P.-P., Zhao, J., Liu, Z.-L., *et al.* Evaluation of antifungal activities and structure-activity relationships of coumarin derivatives. *Pest Manag. Sci.* **73(1)**, 94-101 (2017).

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, 08/29/2023

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