

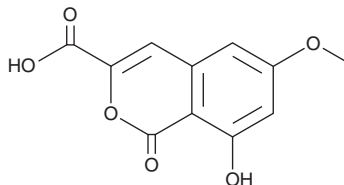
PRODUCT INFORMATION



6-Methoxy-8-hydroxyisocoumarin-3-carboxylic Acid

Item No. 43835

CAS Registry No.: 39051-96-2
Formal Name: 8-hydroxy-6-methoxy-1-oxo-1H-2-benzopyran-3-carboxylic acid
MF: $C_{11}H_8O_6$
FW: 236.2
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Synthetic



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

Description

6-Methoxy-8-hydroxyisocoumarin-3-carboxylic acid is an isocoumarin that has been found in *A. ochraceus* and has anti-angiogenic activity.^{1,2} It inhibits Sarcoma 180 cell-induced angiogenesis in an implanted dorsal air sac in mice when administered at a dose of 100 mg/kg per day.²

References

1. Yamazaki, M., Maebayashi, Y., and Miyaki, K. Isolation of a new metabolite, 6-methoxy-8-hydroxyisocoumarin-3-carboxylic acid from *Aspergillus ochraceus* Wilh. *Chem. Pharm. Bull.* **20**(10), 2276-2278 (1972).
2. Matsumoto, N., Nakashima, T., Isshiki, K., *et al.* Synthesis and biological evaluation of cytogenin derivatives. *J. Antibiot. (Tokyo)* **54**(3), 285-296 (2001).

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, 07/21/2026

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