

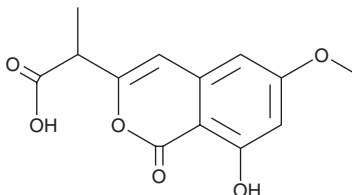
PRODUCT INFORMATION



NM-3

Item No. 40219

CAS Registry No.: 181427-78-1
Formal Name: 8-hydroxy-6-methoxy- α -methyl-1-oxo-1H-2-benzopyran-3-acetic acid
MF: $C_{13}H_{12}O_6$
FW: 264.2
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

NM-3 is supplied as a solid. A stock solution may be made by dissolving the NM-3 in the solvent of choice, which should be purged with an inert gas. NM-3 is soluble in methanol and DMSO.

Description

NM-3 is a synthetic isocoumarin and an inhibitor of angiogenesis.¹ It inhibits S180 cell-induced angiogenesis in an implanted dorsal air sac in mice when administered at doses of 1, 3, and 10 mg/kg. NM-3 (10, 20, and 40 mg/kg) induces intratumoral apoptosis and reduces tumor microvascular density in an SGC-7901 gastric cancer mouse xenograft model.² It decreases renal levels of VEGF, TGF- β 1, and angiopoietin-2 (ANG-2) and renal fibrosis in the *db/db* transgenic mouse model of diabetes when administered at a dose of 100 mg/kg.³

References

1. Nakashima, T., Hirano, S., Agata, N., *et al.* Inhibition of angiogenesis by a new isocoumarin, NM-3. *J. Antibiot. (Tokyo)* **52**(4), 426-428 (1999).
2. Ma, J., Zhu, Z., Zhu, J., *et al.* Study of angiogenesis inhibitor NM-3 on the growth of human gastric cancer. *Wannan Yixueyuan Xuebao* **25**(4), 265-267 (2006).
3. Ichinose, K., Maeshima, Y., Yamamoto, Y., *et al.* 2-(8-hydroxy-6-methoxy-1-oxo-1H-2-benzopyran-3-yl) Propionic acid, an inhibitor of angiogenesis, ameliorates renal alterations in obese type 2 diabetic mice. *Diabetes* **55**(5), 1232-1242 (2006).

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, 05/21/2025

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