

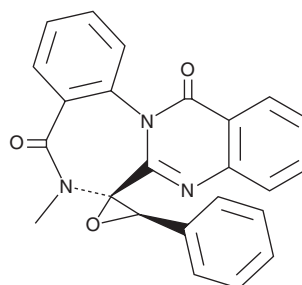
PRODUCT INFORMATION



Benzomalvin C

Item No. 22064

CAS Registry No.: 157047-98-8
Formal Name: (2R,3S)-rel-(+)-6'-methyl-3-phenyl-spiro[oxirane-2,7'(13'H)-quinazolino[3,2-a][1,4]benzodiazepine]-5',13'(6'H)-dione
MF: C₂₄H₁₇N₃O₃
FW: 395.4
Purity: ≥95%
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

Benzomalvin C is supplied as a solid. A stock solution may be made by dissolving the benzomalvin C in the solvent of choice, which should be purged with an inert gas. Benzomalvin C is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Benzomalvin C is a weak antagonist of the neurokinin-1 (NK₁) receptor inhibiting binding of substance P by 46% when used at 100 µg/ml *in vitro*.¹ It is also a weak inhibitor of indoleamine 2,3-dioxygenase (IDO) with an IC₅₀ value of 130 µM for recombinant IDO.² It was isolated from *Penicillium* and contains an epoxide group at C-19 and C-20, which is not present in benzomalvins A, B, or E.

References

1. Sun, H.H., Barrow, C.J., Sedlock, D.M., *et al.* Benzomalvins, new substance P inhibitors from a *Penicillium* sp. *J. Antibiot. (Tokyo)* **47(5)**, 515-522 (1994).
2. Jang, J.-P., Jang, J.-H., Soung, N.-K., *et al.* Benzomalvin E, an indoleamine 2,3-dioxygenase inhibitor isolated from *Penicillium* sp. FN070315. *J. Antibiot. (Tokyo)* **65(4)**, 215-217 (2012).

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, 11/14/2022

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