

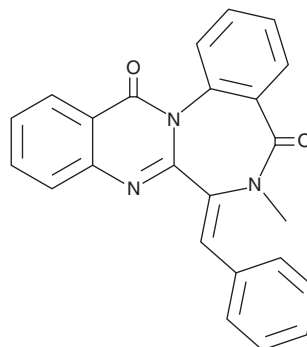
PRODUCT INFORMATION



Benzomalvin B

Item No. 25088

CAS Registry No.: 157047-97-7
Formal Name: (7E)-(+)-6,7-dihydro-6-methyl-7-(phenylmethylene)-quinazolino[3,2-a][1,4]benzodiazepine-5,13-dione
MF: C₂₄H₁₇N₃O₂
FW: 379.4
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 B is supplied as a solid. A stock solution may be made by dissolving the benzomalvin B in the solvent of choice, which should be purged with an inert gas. Benzomalvin B is soluble in organic solvents such as methanol, ethanol, DMSO, and dimethyl formamide.

Description

Benzomalvin B is a fungal metabolite originally isolated from *Penicillium*.¹ It acts as an antagonist of neurokinin-1 (NK₁) receptors, inhibiting binding of substance P (Item No. 24035) by 24% *in vitro* when used at a concentration of 100 µg/ml.

Reference

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).

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, 12/19/2022

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