

PRODUCT INFORMATION



Mazethramycin B

Item No. 43834

CAS Registry No.: 68373-95-5

Formal Name: N-methyl-3-(5,10,11,11a-tetrahydro-9-hydroxy-11-methoxy-8-methyl-5-oxo-1H-pyrrolo[2,1-c][1,4]benzodiazepin-2-yl)-2-propenamide

Synonym: NSC 326670

MF: $C_{18}H_{21}N_3O_4$

FW: 343.4

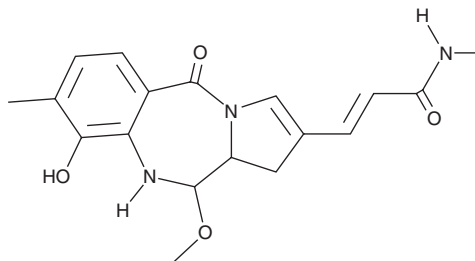
Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Bacterium/*Streptomyces thioluteus* ME561-L4



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

Description

Mazethramycin B is an antitumor antibiotic that has been isolated from *S. thioluteus*.¹ It is active against nine bacteria, including *S. aureus*, *B. subtilis*, and *K. pneumoniae* (MICs = 3.12, 1.56, and 3.12 mg/ml, respectively), and the fungus *C. pseudotropicalis* (MIC = 6.25 mg/ml). Mazethramycin B (31.3 mg/kg per day) increases survival in a L1210 murine leukemia model.

Reference

1. Kunimoto, S., Masuda, T., Kanbayashi, N., *et al.* Mazethramycin, a new member of anthramycin group antibiotics. *J. Antibiot. (Tokyo)* **33**(6), 665-667 (1980).

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, 07/09/2025

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