

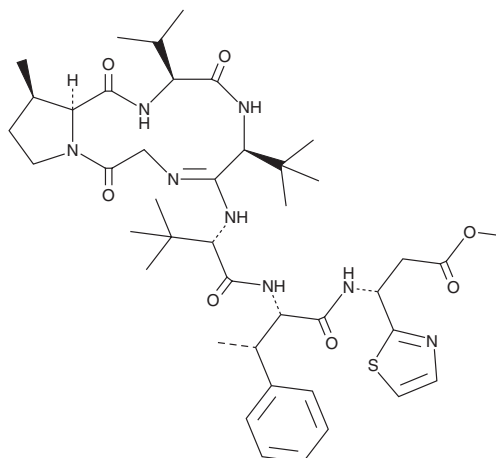
PRODUCT INFORMATION



Bottromycin A₂

Item No. 32921

CAS Registry No.: 15005-62-6
Formal Name: (3R)-N-[(2S)-2-[[glycyl-(3R)-3-methyl-L-prolyl-L-valyl]amino]-1-imino-3,3-dimethylbutyl]-3-methyl-L-valyl-(βS)-β-methyl-L-phenylalanyl-3-(2-thiazolyl)-β-alanine, cyclic (1→1')-imine, methyl ester
MF: C₄₂H₆₂N₈O₇S
FW: 823.1
Purity: ≥80%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces* No. 3668-L2



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

Laboratory Procedures

Bottromycin A₂ is supplied as a solid. A stock solution may be made by dissolving the bottromycin A₂ in the solvent of choice. Bottromycin A₂ is soluble in organic solvents such as methanol and ether, which should be purged with an inert gas. Bottromycin A₂ is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Bottromycin A₂ is an antibiotic originally isolated from *Streptomyces*.¹ It blocks the binding of aminoacyl-tRNA to the A site of the 50S ribosome, inhibiting protein synthesis.² Bottromycin A₂ is active against clinical isolates of methicillin-resistant *Staphylococcus aureus* (MRSA) and vancomycin-resistant *Enterococci* with MIC values of 1 and 0.5 μg/ml, respectively.

References

1. Nakamura, S., Yajima, T., Lin, Y., *et al.* Isolation and characterization of bottromycins A2, B2, C2. *J. Antibiot. (Tokyo)* **20(1)**, 1-5 (1967).
2. Gouda, H., Kobayashi, Y., Yamada, T., *et al.* Three-dimensional solution structure of bottromycin A₂: A potent antibiotic active against methicillin-resistant *Staphylococcus aureus* and vancomycin-resistant *Enterococci*. *Chem. Pharm. Bull. (Tokyo)* **60(2)**, 169-171 (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, 12/02/2022

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