

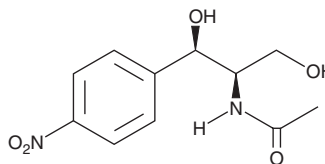
PRODUCT INFORMATION



Corynecin I

Item No. 23237

CAS Registry No.: 4423-58-9
Formal Name: N-[(1R,2R)-2-hydroxy-1-(hydroxymethyl)-2-(4-nitrophenyl)ethyl]-acetamide
MF: C₁₁H₁₄N₂O₅
FW: 254.2
Purity: ≥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

Corynecin I is supplied as a solid. A stock solution may be made by dissolving the corynecin I in the solvent of choice, which should be purged with an inert gas. Corynecin I is soluble in organic solvents such as methanol and acetonitrile.

Description

Corynecin I is a chloramphenicol-like antibiotic originally isolated from *Corynebacterium*.¹ It is active against Gram-positive and Gram-negative bacteria with MIC values of 5.2-83 µg/ml for *S. sonnei*, *P. vulgaris*, *K. pneumoniae*, and *S. aureus* among others.²

References

1. Shirhata, K., Hayashi, T., Deguchi, T., *et al.* The structures of corynecins; chloramphenicol analogues produced by a *n*-paraffin-grown bacterium. *Agric. Biol. Chem.* **36(12)**, 2229-2232 (1972).
2. Suzuki, T., Honda, H., and Katsumata, R. Production of antibacterial compounds analogous to chloramphenicol by a *n*-paraffin-grown bacterium. *Agr. Biol. Chem.* **36(12)**, 2223-2228 (1972).

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, 11/09/2023

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