

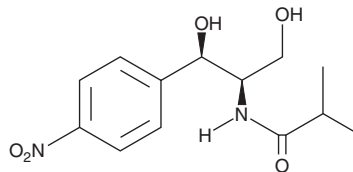
# PRODUCT INFORMATION



## Corynecin III

Item No. 25450

**CAS Registry No.:** 18048-95-8  
**Formal Name:** N-[(1R,2R)-2-hydroxy-1-(hydroxymethyl)-2-(4-nitrophenyl)ethyl]-2-methyl-propanamide  
**MF:** C<sub>13</sub>H<sub>18</sub>N<sub>2</sub>O<sub>5</sub>  
**FW:** 282.3  
**Purity:** ≥99%  
**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 III is supplied as a solid. A stock solution may be made by dissolving the corynecin III in the solvent of choice, which should be purged with an inert gas. Corynecin III is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Corynecin III is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

Corynecin III is a chloramphenicol-like antibiotic originally isolated from *Corynebacterium*.<sup>1</sup> It inhibits the growth of Gram-positive and Gram-negative bacteria, including *S. faecalis*, *S. aureus*, *B. subtilis*, *E. coli*, *P. aeruginosa*, *P. vulgaris*, *S. sonnei*, *S. typhosa*, and *K. pneumoniae* (MICs = 2.6-83 µg/ml).

### Reference

1. 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, 12/20/2022

#### 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