

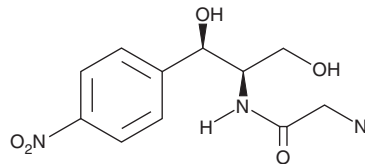
PRODUCT INFORMATION



Azidamfenicol

Item No. 25449

CAS Registry No.: 13838-08-9
Formal Name: 2-azido-N-[(1R,2R)-2-hydroxy-1-(hydroxymethyl)-2-(4-nitrophenyl)ethyl]-acetamide
MF: C₁₁H₁₃N₅O₅
FW: 295.3
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

Azidamfenicol is supplied as a solid. A stock solution may be made by dissolving the azidamfenicol in the solvent of choice, which should be purged with an inert gas. Azidamfenicol is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Azidamfenicol is a broad-spectrum chloramphenicol-like antibiotic that is active against Gram-negative and Gram-positive bacteria.¹ It inhibits ribosomal peptidyltransferase ($K_i = 22 \mu\text{M}$).²

References

- Poitschek, C., Thurner, J., and Haller, I. Determination of antimicrobial activity of a combination of azidoamphenicol, clotrimazole and dexamethasone in vitro. *Arzneimittel-Forschung* **28(2)**, 232-234 (1978).
- Bhuta, P., Chung, H.L., Hwang, J.-S., et al. Analogues of chloramphenicol: Circular dichroism spectra, inhibition of ribosomal peptidyltransferase, and possible mechanism of action. *J. Med. Chem.* **23(12)**, 1299-1305 (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, 10/04/2022

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