

PRODUCT INFORMATION



Azathramycin

Item No. 46202

CAS Registry No.: 76801-85-9

Formal Name: (2R,3S,4R,5R,8R,10R,11R,12S,13S,14R)-13-[[[2,6-dideoxy-3-C-methyl-3-O-methyl- α -L-ribohexopyranosyl]oxy]-2-ethyl-3,4,10-trihydroxy-3,5,8,10,12,14-hexamethyl-11-[[[3,4,6-trideoxy-3-(dimethylamino)- β -D-xylohexopyranosyl]oxy]-1-oxa-6-azacyclopentadecan-15-one

Synonyms: Azaerythromycin A, N-desmethyl Azithromycin

MF: $C_{37}H_{70}N_2O_{12}$

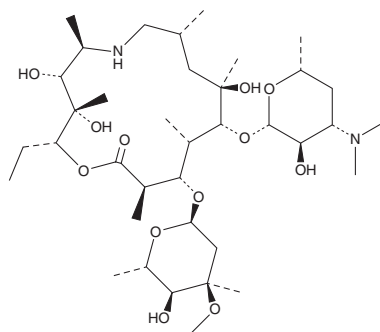
FW: 735.0

Purity: $\geq 95\%$

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

Azathramycin is supplied as a solid. A stock solution may be made by dissolving the azathramycin in the solvent of choice, which should be purged with an inert gas. Azathramycin is sparingly soluble (1-10 mg/ml) in ethanol.

Description

Azathramycin is a macrolide antibiotic.¹ It is active against *S. pyogenes*, *S. aureus*, *E. coli*, and several strains of *H. influenzae* (MICs = ≤ 0.025 , 0.78, 0.78, and 0.78-1.56 $\mu\text{g/ml}$, respectively). Azathramycin is also active against *T. gondii* (IC₅₀ = 65 μM).² Subcutaneous administration of azathramycin protects mice against an acutely fatal *S. aureus* infection with a half-maximal protective dose (PD₅₀) value of 9 mg/kg.¹

References

1. Bright, G.M., Nagel, A.A., Bordner, J., et al. Synthesis, *in vitro* and *in vivo* activity of novel 9-deoxo-9a-AZA-9a-homoerythromycin A derivatives; a new class of macrolide antibiotics, the azalides. *J. Antibiot.* **41**(8), 1029-1047 (1988).
2. Lee, Y., Choi, J.Y., Fu, H., et al. Chemistry and biology of macrolide antiparasitic agents. *J. Med. Chem.* **54**(8), 2792-2804 (2011).

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, 08/19/2026

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