

PRODUCT INFORMATION



Aurodox

Item No. 27584

CAS Registry No.: 12704-90-4
Formal Name: (α S,2R,3R,4R,6S)-N-[(2E,4E,6S,7R)-7-[(2S,3S,4R,5R)-5-[(1E,3E,5E)-7-(1,2-dihydro-4-hydroxy-1-methyl-2-oxo-3-pyridinyl)-6-methyl-7-oxo-1,3,5-heptatrien-1-yl]tetrahydro-3,4-dihydroxy-2-furanyl]-6-methoxy-5-methyl-2,4-octadien-1-yl]- α -ethyltetrahydro-2,3,4-trihydroxy-5,5-dimethyl-6-(1E,3Z)-1,3-pentadien-1-yl]-2H-pyran-2-acetamide

Synonyms: Antibiotic X 5108, Goldinodox, Goldinomycin, 1-methyl-Mocimycin

MF: C₄₄H₆₂N₂O₁₂

FW: 811.0

Purity: \geq 85% (mixture of tautomers)

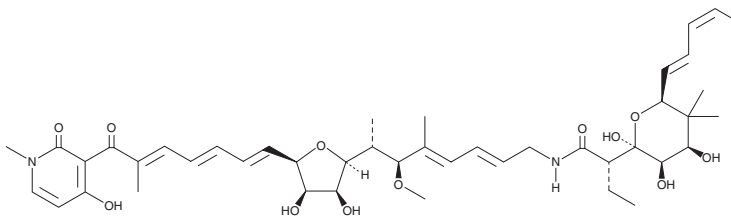
Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years

Item Origin: Bacterium/*Streptomyces* sp.

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



Laboratory Procedures

Aurodox is supplied as a solid. A stock solution may be made by dissolving the aurodox in the solvent of choice, which should be purged with an inert gas. Aurodox is soluble in organic solvents such as ethanol at a concentration of approximately 1 mg/ml. Aurodox is also soluble in methanol and DMSO.

Description

Aurodox is a polyketide antibiotic originally isolated from *S. goldiniensis*.¹ It is active against Gram-positive bacteria, including *B. megaterium*, *B. anthracis*, and *M. hominis* (MICs = 0.06, 0.6, and 3-10 μ g/ml, respectively), as well as *Pneumococcus* and *Streptomyces* species (MICs = 3-12 and 3-30 μ g/ml, respectively). It is effective against *S. pyogenes* infection in mice with a 50% curative dose (CD₅₀) value of 71 mg/kg. Aurodox inhibits bacterial protein synthesis by binding to bacterial elongation factor Tu (EF-Tu) and inhibiting its release from the ribosome.^{2,3}

References

- Berger, J., Lehr, H.A., Teitel, S., *et al.* A new antibiotic X-5108 OF *Streptomyces* origin. I. Production, isolation and properties. *J. Antibiot. (Tokyo)* **26(1)**, 15-22 (1973).
- Wolf, H., Chinali, G., and Parmeggiani, A. Mechanism of the inhibition of protein synthesis by kirromycin. Role of elongation factor Tu and ribosomes. *Eur. J. Biochem.* **75(1)**, 67-75 (1977).
- Bhuta, P. and Chládek, S. Stimulation of *Escherichia coli* elongation factor Tu-dependent GTP hydrolysis by aminoacyl oligonucleotides in the presence of aurodox. *FEBS Lett.* **122(1)**, 113-116 (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, 12/08/2022

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