

PRODUCT INFORMATION



Epioxytetracycline

Item No. 27987

CAS Registry No.: 14206-58-7
Formal Name: (4R,4aR,5S,5aR,6S,12aS)-4-(dimethylamino)-1,4,4a,5,5a,6,11,12a-octahydro-3,5,6,10,12,12a-hexahydroxy-6-methyl-1,11-dioxo-2-naphthacene-carboxamide

Synonym: 4-*epi*-Oxytetracycline

MF: C₂₂H₂₄N₂O₉

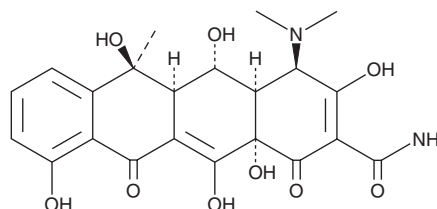
FW: 460.4

Purity: ≥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

Epioxytetracycline is supplied as a solid. A stock solution may be made by dissolving the epioxytetracycline in the solvent of choice, which should be purged with an inert gas. Epioxytetracycline is slightly soluble in methanol (warmed) and DMSO.

Description

Epioxytetracycline is a degradation product of the antibiotic oxytetracycline (Item No. 18076).^{1,2} It has been found in swine manure compost and wastewater and is considered a pollutant.

References

1. Wu, X., Wei, Y., Zheng, J., *et al.* The behavior of tetracyclines and their degradation products during swine manure composting. *Bioresour. Technol.* **102(10)**, 5924-5931 (2011).
2. Çelik, A., Casey, E., and Hasar, H. Degradation of oxytetracycline under autotrophic nitrifying conditions in a membrane aerated biofilm reactor and community fingerprinting. *J. Hazard. Mater.* (2018).

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, 12/02/2022

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