

PRODUCT INFORMATION



Anhydrochlortetracycline (hydrochloride)

Item No. 19941

CAS Registry No.: 65490-24-6
Formal Name: (4S,4aS,12aS)-7-chloro-4-(dimethylamino)-1,4,4a,5,12,12a-hexahydro-3,10,11,12a-tetrahydroxy-6-methyl-1,12-dioxo-2-naphthacenecarboxamide, monohydrochloride

MF: C₂₂H₂₁ClN₂O₇ • HCl

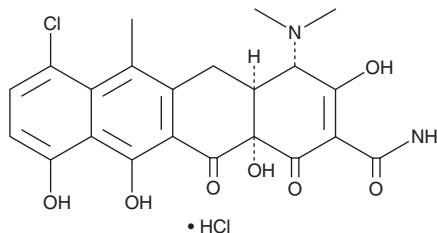
FW: 497.3

Purity: ≥90%

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

Anhydrochlortetracycline (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the anhydrochlortetracycline (hydrochloride) in the solvent of choice. Anhydrochlortetracycline (hydrochloride) is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Anhydrochlortetracycline is an impurity of the tetracycline antibiotic chlortetracycline formed by acid degradation.¹ It is a poor inhibitor of protein synthesis and has minor antimicrobial activity against various strains of actinomycetes.^{1,2}

References

1. Goodman, J.J., Matrishin, N.M., and Backus, E.J. The effect of anhydrochlortetracycline on the growth of actinomycetes. *J. Bacteriol.* **69(1)**, 70-72 (1955).
2. Rasmussen, B., Noller, H.F., Daubresse, G., *et al.* Molecular basis of tetracycline action: Identification of analogs whose primary target is not the bacterial ribosome. *Antimicrob. Agents Chemother.* **35(11)**, 2306-2311 (1991).

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, 12/22/2022

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