

PRODUCT INFORMATION



4-Epianhydrochlortetracycline (hydrochloride)

Item No. 23146

CAS Registry No.: 158018-53-2
Formal Name: (4R,4aS,12aS)- 7-chloro-4-(dimethylamino)-1,4,4a,5,12a,12a-hexahydro-3,10,11,12a-tetrahydroxy-6-methyl-1,12-dioxo-2-naphthacenecarboxamide, monohydrochloride

MF: C₂₂H₂₁ClN₂O₇ • HCl

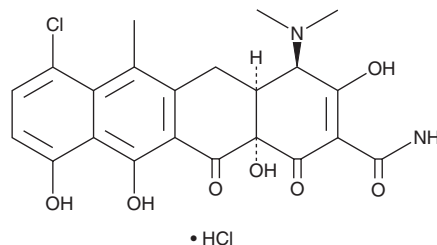
FW: 497.3

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

4-Epianhydrochlortetracycline (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the 4-epianhydrochlortetracycline (hydrochloride) in the solvent of choice, which should be purged with an inert gas. 4-Epianhydrochlortetracycline (hydrochloride) is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

4-Epianhydrochlortetracycline is a derivative of tetracycline (Item No. 14328).¹ It inhibits matrix metalloproteinase-9 (MMP-9) activity in a dose-dependent manner in an assay for collagen degradation. 4-Epianhydrochlortetracycline also completely inhibits neutrophil-derived MMP-9 degradation of denatured collagen at a concentration of 500 µg/ml. Unlike tetracycline, 4-epianhydrochlortetracycline lacks antibiotic activity.

Reference

1. Paemen, L., Martens, E., Norga, K., *et al.* The gelatinase inhibitory activity of tetracyclines and chemically modified tetracycline analogues as measured by a novel microtiter assay for inhibitors. *Biochem. Pharmacol.* **52(1)**, 105-111 (1996).

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, 11/03/2022

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