

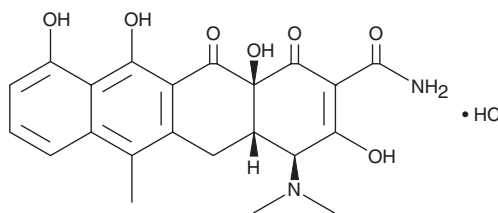
Product Information



Anhydrotetracycline (hydrochloride)

Catalog No. 10009542

CAS Registry No.: 13803-65-1
Formal Name: 4-(dimethylamino)-1,4S,4aS,5,12,12aS-hexahydro-3,10,11,12a-tetrahydroxy-6-methyl-1,12-dioxo-mono-hydrochloride-2-naphthacene-carboxamide
MF: C₂₂H₂₂N₂O₇ • HCl
FW: 462.9
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that anhydrotetracycline (hydrochloride) be stored as supplied at -20°C. It should be stable for at least two years.

Anhydrotetracycline (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the anhydrotetracycline (hydrochloride) in an organic solvent purged with an inert gas. Anhydrotetracycline (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of anhydrotetracycline (hydrochloride) in ethanol is approximately 2 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of anhydrotetracycline (hydrochloride) can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of anhydrotetracycline (hydrochloride) in PBS, pH 7.2, is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

The tetracycline repressor (TetR) is a transcriptional regulator which normally binds tightly to its palindromic tetO operator DNA, blocking gene expression.¹ Tet causes the repressor to dissociate from the DNA, allowing transcription to occur. A novel reverse TetR (revTetR) requires tetracycline as a co-repressor to bind tetO and block transcription.^{2,3} Anhydrotetracycline (hydrochloride) is a powerful effector in both the TetR and revTetR systems, binding the TetR 35-fold more strongly than Tet.^{1,4} Moreover, anhydrotetracycline poorly binds the 30S ribosomal subunit, compared to Tet,⁵ so it does not act as a general inhibitor of translation and is a poor antibiotic. Perhaps related to this, the concentration of anhydrotetracycline that inhibits eukaryotic cell growth is more than a 1,000-fold above the dose that alters transcription through TetR.¹

References

1. Gossen, M. and Bujard, H. *Nucleic Acids Res.* **21(18)**, 4411-4412 (1993).
2. Kamionka, A., Bogdanska-Urbaniak, J., Scholz, O., et al. *Nucleic Acids Res.* **32(2)**, 842-847 (2004).
3. Resch, M., Striegl, H., Henssler, E.M., et al. *Nucleic Acids Res.* **36(13)**, 4390-4401 (2008).
4. Degenkolb, J., Takahashi, M., Ellestad, G.A., et al. *Antimicrob. Agents Chemother.* **35(8)**, 1591-1595 (1991).
5. Rasmussen, B., Noller, H.F., Daubresse, G., et al. *Antimicrob. Agents Chemother.* **35(11)**, 2306-2311 (1991).

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/01/2009