

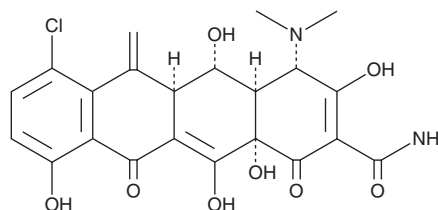
PRODUCT INFORMATION



Meclocycline

Item No. 27940

CAS Registry No.: 2013-58-3
Formal Name: (4S,4aR,5S,5aR,12aS)-7-chloro-4-(dimethylamino)-1,4,4a,5,5a,6,11,12a-octahydro-3,5,10,12,12a-pentahydroxy-6-methylene-1,11-dioxo-2-naphthacencarboxamide
Synonyms: GS-2989, NSC 78502
MF: C₂₂H₂₁ClN₂O₈
FW: 476.9
Purity: ≥95%
UV/Vis.: λ_{max}: 244, 354 nm
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

Meclocycline is supplied as a solid. A stock solution may be made by dissolving the meclocycline in the solvent of choice, which should be purged with an inert gas. Meclocycline is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Description

Meclocycline is a tetracycline antibiotic.¹ It has broad antimicrobial activity *in vitro*, including growth inhibition of the bacterium *S. epidermidis* when used at a concentration of 0.8 µg/ml.^{1,2} Meclocycline inhibits lactate dehydrogenase (LDH) release and cell toxicity induced by expression of *htt-N63-148Q*, a mutant form of the huntingtin gene, without altering *htt-N63-148Q* protein levels in PC12 cells.³

References

1. Blackwood, R.K., Rennhard, H.H., Beereboom, J.J., *et al.* 6-Methylenetetracyclines. I. A new class of tetracycline antibiotics. *J. Am. Chem. Soc.* **83(12)**, 2773-2775 (1961).
2. Pulverer, G. and Jeljaszewicz, J. Susceptibility of *Staphylococcus epidermidis* to lysostaphin and major antimicrobial agents. *Chemotherapy* **19(2)**, 109-114 (1973).
3. Wang, W., Duan, W., Igarashi, S., *et al.* Compounds blocking mutant huntingtin toxicity identified using a Huntington's disease neuronal cell model. *Neurobiol. Dis.* **20(2)**, 500-508 (2005).

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, 10/28/2022

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