

Chlortetracycline . hydrochloride

Protein synthesis inhibitor

Antibiotic. Inhibits protein synthesis in prokaryotes and eukaryotes. Inhibits binding of aminoacyl-tRNA to ribosomes. Used as a fluorescent Ca^{2+} probe.

Citations: 5

[View Online »](#)

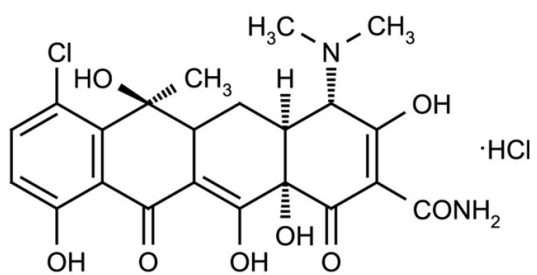
Ordering Information

[Order Online »](#)

ALX-350-238-G005	5g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Aureomycin . HCl

Appearance Yellow powder.

CAS 64-72-2

Formula $C_{22}H_{23}ClN_2O_8 \cdot HCl$

MI **14:** 2192

MW 515.4

Purity ≥85% (UHPLC)

Solubility Soluble in 1M NaOH (50mg/ml).

Source Isolated from *Streptomyces aureofaciens*.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)