

Doxycycline . hyclate

MMP inhibitor

Broad spectrum antibiotic and bacteriostatic. Shows antiprotozoal properties. Potent inhibitor of MMPs (matrix metalloproteinases) in vivo. Inhibits collagen synthesis. Increases and promotes smooth muscle cell adhesion. Inhibits proliferation and migration. Protects microvasculature by inhibiting plasminogen systems.

Citations: 16

[View Online »](#)

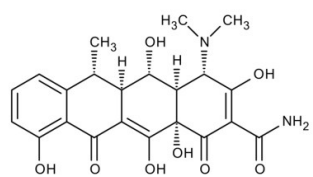
Ordering Information

[Order Online »](#)

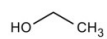
ALX-380-273-G001	1g
ALX-380-273-G005	5g

Manuals, SDS & CofA

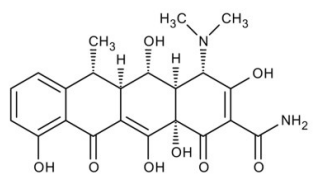
[View Online »](#)



HCl



H_2O



HCl

Handling & Storage

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

Handling Hygroscopic.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Doxycycline . hydrochloride . hemiethanolate . hemihydrate

Appearance Yellow crystalline solid.

CAS 24390-14-5

Couple Target MMP

Couple Type Inhibitor

Formula $C_{46}H_{58}Cl_2N_4O_{18}$

MI **14:** 3440

MW 1025.87 g/mol

Purity ≥98% (HPLC)

RTECS QI8925000

Solubility Soluble in water (50mg/ml).

Source Semisynthetic from oxytetracycline.



ENZO LIFE SCIENCES,
INC.
May 29, 2024
info-
usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-
de@enzolifesciences.com

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