

Daunorubicin . hydrochloride

Antitumor antibiotic

Anthracycline anticancer antibiotic. Induces DNA damage by intercalating into DNA and inhibiting topoisomerase I and II. Inhibits RNA and DNA synthesis. Induces DNA single strand breaks and apoptosis in tumor cells.

Citations: 15

[View Online »](#)

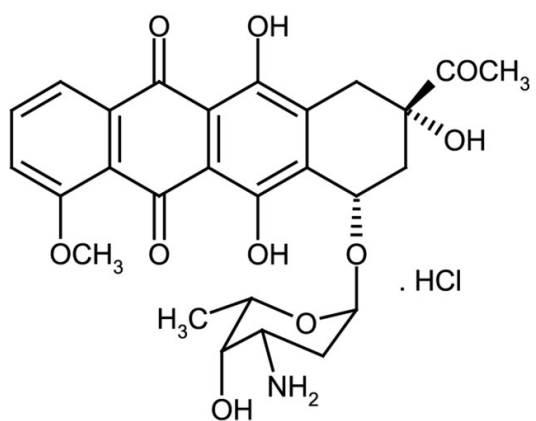
Ordering Information

[Order Online »](#)

ALX-380-043-M010	10mg
------------------	------

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 Daunomycin . HCl

Appearance Orange to red crystalline powder.

CAS 23541-50-6

Couple Target Topoisomerase

Couple Type Inhibitor

Formula $C_{27}H_{29}NO_{10} \cdot HCl$

MI 14: 2832

MW 527.5 . 36.5

Purity ≥90% (Assay)

Solubility Soluble in water or methanol.

Source Isolated from *Streptomyces spec.*

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-
fr@enzolifesciences.com](mailto:info-fr@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)