

Idarubicin . hydrochloride

Topoisomerase II α inhibitor

More potent, lipophilic and antineoplastic analog of daunorubicin. DNA-damaging effect includes DNA oxidation and methylation, DNA intercalation, inhibition of DNA synthesis, induction of DNA strand breaks and delay of cell cycle progression. Inhibitor of topoisomerase II α . Produces endonucleolytic cleavage and is a marker of apoptosis.

Citations: 5

[View Online »](#)

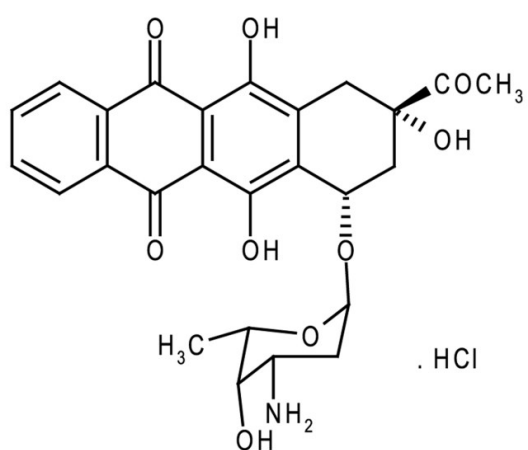
Ordering Information

[Order Online »](#)

ALX-380-260-M005	5mg
------------------	-----

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 Idamycin, DMDR, 4-Demethoxydaunorubicin

Appearance Orange crystalline solid.

CAS 57852-57-0

Couple Target Topoisomerase

Couple Type Inhibitor

Formula $C_{26}H_{27}NO_9 \cdot HCl$

Identity Determined by IR.

MI 14: 4886

MW 497.5 . 36.5

Purity ≥98% (HPLC)

RTECS HB7877000

Solubility Soluble in water or methanol.

Source Synthetic.

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)