

Doxorubicin . HCl

Topoisomerase inhibitor

A broad spectrum antitumor agent. Induces DNA damage by intercalating into DNA and inhibiting topoisomerase II. Binds covalently to DNA. Inhibits reverse transcriptase, RNA polymerase and the catalytic activity of Dnmt1. Immunosuppressive. Antineoplastic. Induces apoptosis. Down regulates c-myc and c-jun in doxorubicin-sensitive cells.

Citations: 27

[View Online »](#)

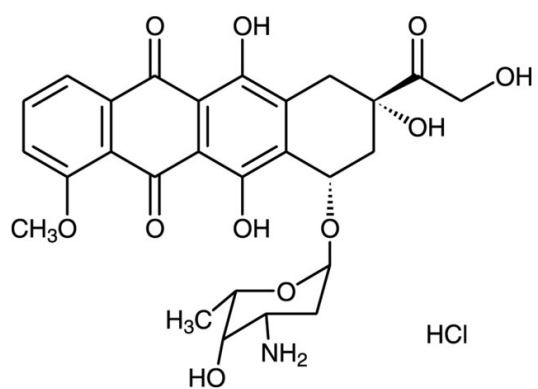
Ordering Information

[Order Online »](#)

BML-GR319-0005	5mg
BML-GR319-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at room temperature. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	DXR . HCl, 14-Hydroxydaunomycin . HCl, Adriamycin . HCl
Appearance	Red solid.
CAS	25316-40-9
Couple Target	c-Jun, c-Myc, DNA, Dnmt, Reverse transcriptase, Topoisomerase
Couple Type	Inhibitor, Ligand
Formula	$C_{27}H_{29}NO_{11} \cdot HCl$
MI	14: 3439
MW	580.0
Purity	≥95% (HPLC)
RTECS	QI9295900
Solubility	Soluble in water.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
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-fr@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