

Mithramycin A

Transcription inhibitor

Mithramycin noncovalently interacts with G-C-rich duplex DNA in the presence of Mg^{2+} or Zn^{2+} . It induces differentiation of leukemic cells accompanied by an early decrease in c-myc expression. It selectively inhibits collagen-1 gene expression in human fibroblasts.

Citations: 13

[View Online »](#)

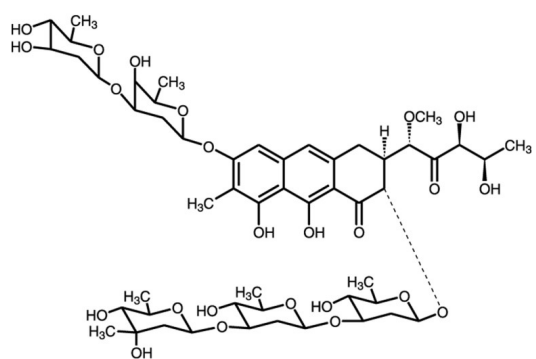
Ordering Information

[Order Online »](#)

BML-GR305-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Plicamycin
Appearance	Yellow solid.
CAS	18378-89-7
Formula	$C_{52}H_{76}O_{24}$
MI	14: 7536
MW	1085.1
Purity	≥95% (HPLC, TLC)
Solubility	Soluble in phosphate buffered saline (1mg/ml), 100% ethanol, DMSO or methanol
Source	Isolated from <i>Streptomyces argillaceus</i>

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)