

Leptomycin B

Nuclear export inhibitor

A potent and specific nuclear export inhibitor. Leptomycin B alkylates and inhibits CRM1/exportin 1, a protein required for nuclear export of proteins containing a nuclear export sequence (NES). In addition to antifungal and antibacterial activities, blocks the cell cycle and is a potent anti-tumor agent. At low nM concentrations, blocks the nuclear export of many proteins including HIV-1 Rev, MAPK/ERK, and NF- κ B/I κ B and it stabilizes the expression of p53. Other proteins that are influenced by leptomycin B are actin, c-Abl, cyclin B1, MDM2/p53, I κ B, MPF and PKA. Leptomycin B also inhibits the export and translation of many RNAs, including COX-2 and c-Fos mRNAs, by inhibiting export of ribonucleoproteins.

Citations: 12

[View Online »](#)

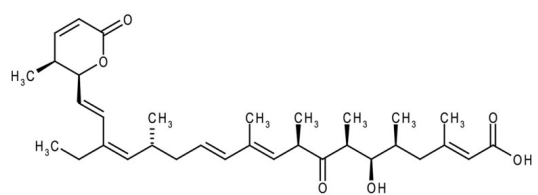
Ordering Information

[Order Online »](#)

ALX-380-100-C100	100 μ g
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Packaged under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	LMB
Appearance	Liquid.
CAS	87081-35-4
Couple Target	Exportin
Couple Type	Inhibitor
Formula	$C_{33}H_{48}O_6$
Formulation	Solution in 100% ethanol.
MW	540.7
Purity	≥96% (HPLC)
Solubility	Soluble in methanol, 100% ethanol, dichloromethane, or DMSO; insoluble in water.
Source	Isolated from <i>Streptomyces sp.</i>



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)