

Monensin . sodium salt

Na⁺ ionophore

Antibiotic that functions as Na⁺ ionophore. Blocks glycoprotein secretion. Inhibits proliferation of several lymphoma cell lines through cell cycle arrest and apoptosis. Blocks ceramide transport through the Golgi apparatus.

Citations: 16

[View Online »](#)

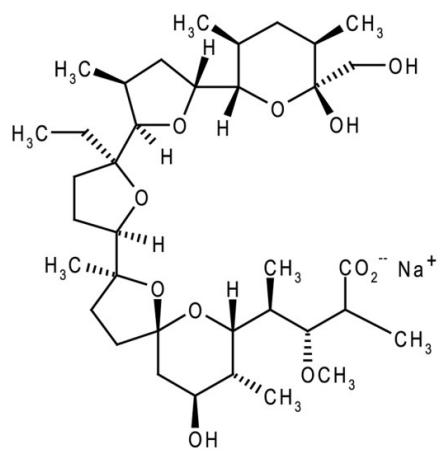
Ordering Information

[Order Online »](#)

ALX-380-026-G001	1g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Store tightly sealed in the dark.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to light yellow powder.
CAS	22373-78-0
Formula	$C_{36}H_{61}O_{11} \cdot Na$
MI	14: 6246
MW	669.9 . 23.0
Purity	≥90%
Solubility	Soluble in 100% ethanol (25mg/ml), methanol (50mg/ml), chloroform or DMSO (25mg/ml). Slightly soluble in water.

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)