

Enniatin A

Ionophore antibiotic

Ionophore antibiotic. Potential anticancer compound.

Enniatins are a family of depsipeptides produced by several species of the filamentous fungus *Fusarium*. Enniatins have initially been shown to act as ionophores. Later, their effects on acyl-CoA cholesterol transferase, ABC-transporters and the selectivity of their antitumor action came to the fore.

Citations: 11

[View Online »](#)

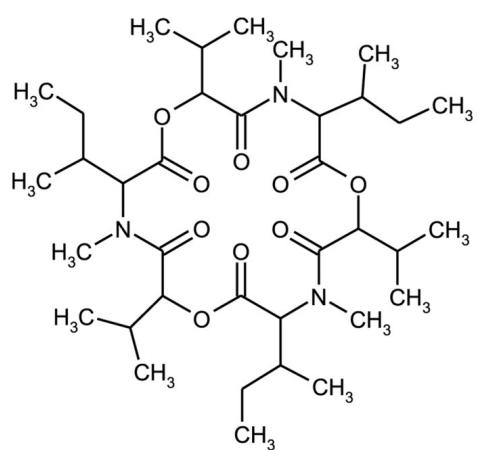
Ordering Information

[Order Online »](#)

ALX-380-310-MC05	0.5mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white crystalline solid.
CAS	2503-13-1
Formula	$C_{36}H_{63}N_3O_9$
Identity	Identity determined by 1H-NMR, 13C-NMR and MS.
MI	14: 3585
MW	681.9
Purity	≥97% (HPLC)
Solubility	Soluble in 100% ethanol, methanol, DMSO and dimethyl formamide; almost insoluble in water.
Source	Isolated from <i>Fusarium</i> sp.

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)