

γ -Amanitin

RNA polymerase II inhibitor

Biological activity like α -amanitin.

Citations: 10

[View Online »](#)

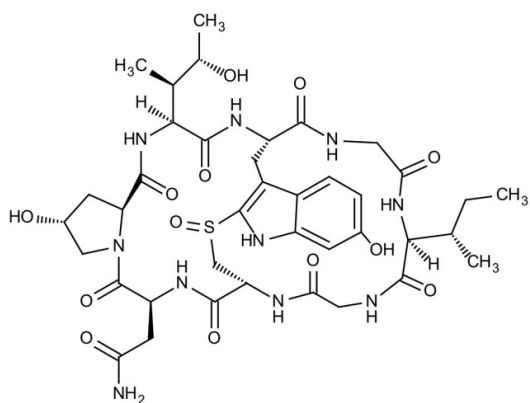
Ordering Information

[Order Online »](#)

ALX-350-272-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Dilute with buffer immediately prior to use. Product is destroyed by concentrated acid or base.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	EL-Amanitin
Appearance	White to off-white powder.
CAS	21150-23-2
Couple Target	RNA polymerase
Couple Type	Inhibitor
Formula	$C_{39}H_{54}N_{10}O_{13}S$
MI	14: 373
MW	903.0
Purity	≥90%
Solubility	Soluble in DMF, DMSO, and 100% methanol.
Source	Isolated from <i>Amanita phalloides</i> mushrooms.

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)