

Phallacidin

Actin ligand

Biological activity like phalloidin (Prod. No. ALX-350-265), but contains a carboxy group for coupling reactions.

Citations: 6

[View Online »](#)

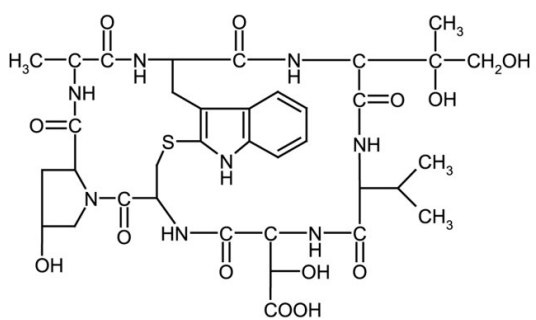
Ordering Information

[Order Online »](#)

ALX-350-269-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name EL-Phalloidin

Appearance White to off-white powder.

CAS 26645-35-2

Couple Target Actin

Couple Type Ligand

Formula $C_{37}H_{50}N_8O_{13}S$

MW 846.9

Purity ≥90%

Solubility Soluble in DMF, DMSO, 100% methanol, and water.

Source Isolated from *Amanita phalloides* 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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)