

Phomopsin A

A potent mycotoxin and anti-mitotic compound. Causes lupinosis in livestock fed infected lupins. Binds selectively to dimeric tubulin, inhibiting the formation of the microtubule spindle to block cell division. Binds to tubulin at a site overlapping that of vinblastine (Prod. No. ALX-350-257) and maytansine. Protects tubulin from decay.

Citations: 3

[View Online »](#)

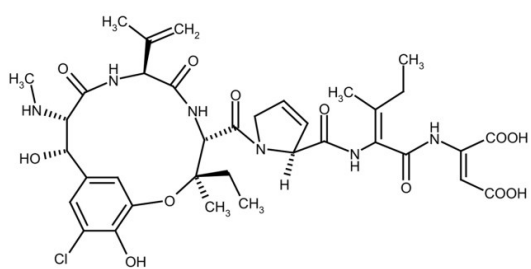
Ordering Information

[Order Online »](#)

ALX-350-417-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 -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White solid.

CAS 64925-80-0

Formula $C_{36}H_{45}ClN_6O_{12}$

MW 789.2

Purity ≥95% (HPLC)

RTECS SY2593000

Solubility Soluble in 100% ethanol, methanol, dimethylformamide or DMSO; slightly soluble in water.

Source Isolated from *Phomopsis leptostromiformis* MST-FP2068.

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)