

Jasplakinolide

F-actin probe

Cell permeable F-actin probe. Induces actin polymerization. Completely inhibits the binding of phalloidin to F-actin. Antifungal and anti-proliferative agent.

Citations: 26

[View Online »](#)

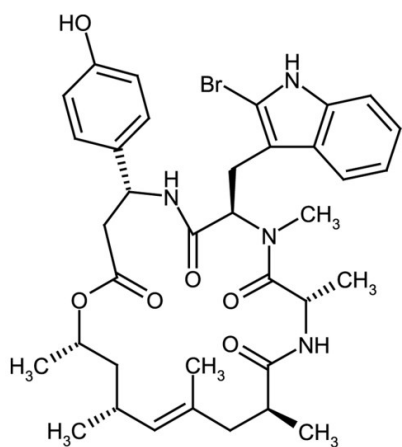
Ordering Information

[Order Online »](#)

ALX-350-275-C050	50µg
ALX-350-275-C100	100µg

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

Alternative Name	Jaspamide
Appearance	Off-white to yellow solid.
CAS	102396-24-7
Formula	$C_{36}H_{45}BrN_4O_6$
Identity	Determined by 1H -NMR and ESI-MS.
MW	709.7
Purity	≥97% (HPLC)
Solubility	Soluble in 100% ethanol, 83% methanol or DMSO.
Source	Isolated from <i>Jaspis splendens</i> .

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)