

Latrunculin A

Ultra-pure; inhibits actin polymerization

Inhibits actin polymerization, microfilament organization and microfilament-mediated processes. Acts via a different mechanism and 10- to 20-fold more potently than cytochalasins. Inhibits phagocytosis by macrophages. Exposure to 0.2 mg/ml for one hour produces maximal change in cell morphology.

Citations: 63

[View Online »](#)

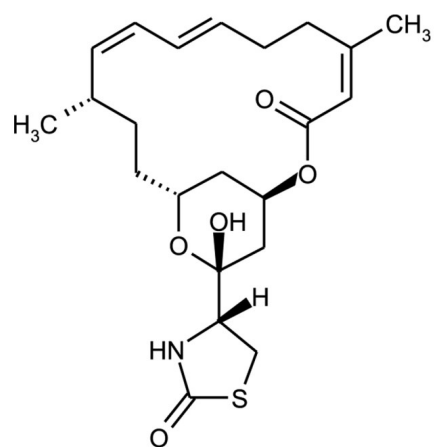
Ordering Information

[Order Online »](#)

BML-T119-0100	100µg
BML-T119-0500	500µg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for 1 year after receipt when stored at -20°C. Stock solutions are stable for up to 3 months when stored at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to yellow waxy solid.
CAS	76343-93-6
Formula	$C_{22}H_{31}NO_5S$
MI	14: 5378
MW	421.6
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/ml).

Technical Info / Product Notes Note: Product is not sterile.



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)