

Mycalolide B

Depolymerizes actin

Mycalolide B is a novel marine natural product which binds to actin in a 1:1 molar ratio ($K_d=13\text{-}20\text{ nM}$). It selectively and completely depolymerizes F-actin to G-actin *via* a mechanism distinct from that of the cytochalasins. Unlike cytochalasins, it inhibits actin-activated Mg^{2+} -ATPase from rabbit skeletal muscle.

Citations: 16

[View Online »](#)

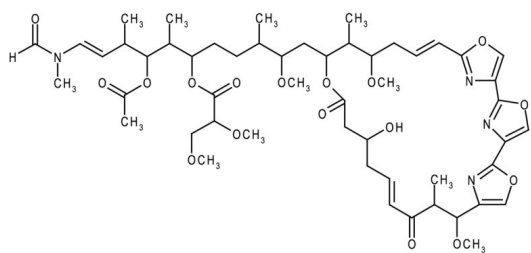
Ordering Information

[Order Online »](#)

BML-T123-0020	20µg
---------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Clear to white solid
CAS	122752-21-0
Couple Target	ATPase
Couple Type	Inhibitor
Formula	$C_{52}H_{74}N_4O_{17}$
MW	1027.2
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO, methanol or isopropanol
Source	Isolated from the marine sponge <i>Mycale</i> sp.

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)