

Ansamitocin P-3

Microtubule inhibitor

Ansamitocin P-3 is a maytansine analog which displays potent cytotoxicity against the human solid tumor cell lines A-549 and HT-29. It shares the same binding site as rhizoxin and phomopsin A on porcine brain tubulin. Maytansine and its analogs cause extensive disassembly of the microtubule and totally prevent tubulin spiralization.

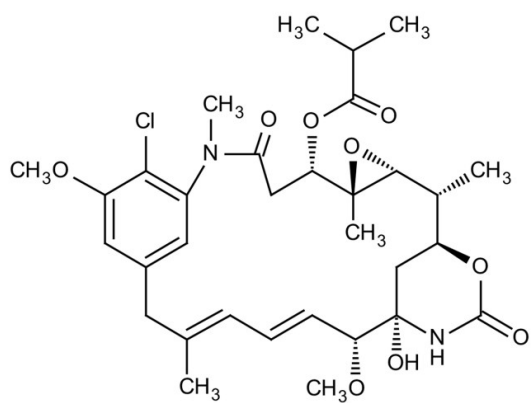
Ordering Information

[Order Online »](#)

BML-T121-0001	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 to off-white crystalline solid.

CAS 66547-09-9

Couple Target Microtubule

Couple Type Inhibitor

Formula $C_{32}H_{43}ClN_2O_9$

MW 635.1

Purity ≥93% (HPLC)

Solubility Soluble in DMSO.

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)