

Herbimycin A

Tyrosine kinase inhibitor

Antibiotic. Cell permeable tyrosine kinase inhibitor. It is effective on Src, Yes, Fps, Ros, Abl and ErbB oncogene products and inhibits thrombin-induced tyrosine phosphorylation of phospholipase C. It also inhibits HSP90, NF- κ B activation and angiogenesis. Increases the sensitivity of certain cancer cells to chemotherapeutic agents. Blocks mitotic activity. Neurotoxic.

Citations: 21

[View Online »](#)

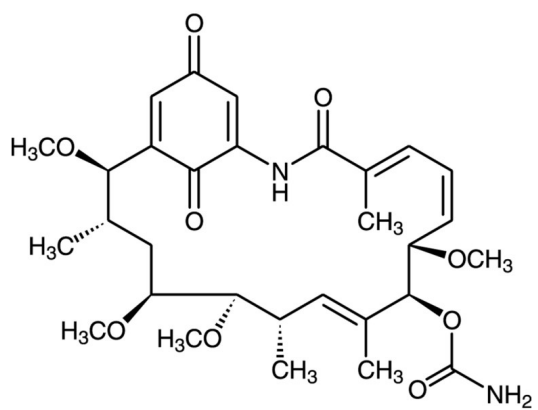
Ordering Information

[Order Online »](#)

BML-EI227-1000	1mg
BML-EI227-0100	100 μ g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	70563-58-5
Couple Target	Abl, c-Ros, EGFR, ErbB, Fps, HSP90, Src kinase, Yes
Couple Type	Inhibitor
Formula	$C_{30}H_{42}N_2O_9$
Formulation	Lyophilized.
MW	574.7
Purity	≥98% (HPLC)
RTECS	LX8930000
Solubility	Soluble in ethanol, methanol, DMF or 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)