

Fumagillin

Methionine aminopeptidase 2 inhibitor

Inhibitor of angiogenesis and endothelial cell proliferation. Specific inhibitor of methionine aminopeptidase type II (MetAP-II) by covalently modifying a conserved active-site histidine. Precursor to TNP-470. Exhibits antitumor activity in estrogen-induced pituitary adenoma. Antineoplastic. Anti-infective.

Citations: 13

[View Online »](#)

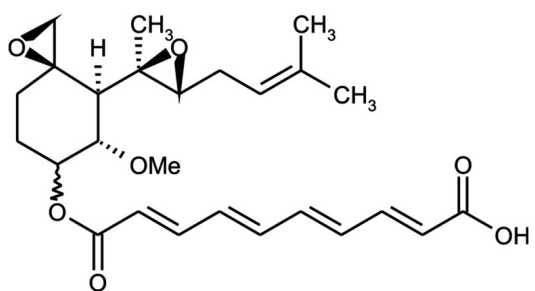
Ordering Information

[Order Online »](#)

BML-CT100-0005	5mg
----------------	-----

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 at -20°C. Stock solutions are stable for up to 3 months at -20°C. Unstable in aqueous solutions with pH >9.0 or pH <5.0.
Handling	Protect from light and oxygen. Keep under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Fumidil B, Fumadil B, TNP-470
Appearance	Off-white to light yellow solid.
CAS	23110-15-8
Couple Target	Aminopeptidase
Couple Type	Inhibitor
Formula	C ₂₆ H ₃₄ O ₇
MI	14: 4286
MW	458.5
Purity	≥95% (TLC)
Solubility	Soluble in DMSO (25mg/ml) or diluted aqueous bases; slightly soluble in methanol or other alcoholic solvents; insoluble in water.
Source	Isolated from <i>Aspergillus fumigatus</i> .
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)