

Chaetocin

Methyltransferase inhibitor

Chaetocin, an antitumor antibiotic, is a thiodioxopiperazine natural product produced by *Chaetomium* species. Specific inhibitor of the lysine-specific methyltransferase SU (VAR)3-9 both *in vitro* and *in vivo*. It displays potent antimyeloma activity in IL-6-dependent myeloma cell lines. Its antimyeloma activity appears to be due to induction of oxidative stress and consequent apoptosis.

Citations: 6

[View Online »](#)

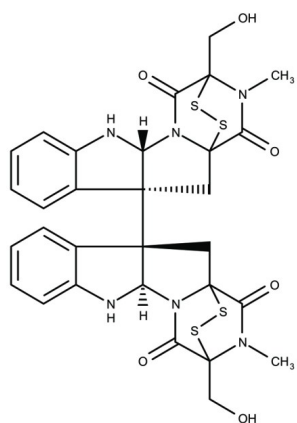
Ordering Information

[Order Online »](#)

BML-GR349-0200	200µg
----------------	-------

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.
Short Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	28097-03-2
Couple Target	Methyltransferase
Couple Type	Inhibitor
Formula	$C_{30}H_{28}N_6O_6S_4$
MW	696.9
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO.
Source	Isolated from <i>Chaetomium sp.</i>

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)