

Thiolutin

RNA polymerase inhibitor

Sulfur containing antibiotic. Potent inhibitor of bacterial and yeast RNA polymerases. Inhibits mRNA chain elongation. Inhibits adhesion of endothelial cells (HUVECs) to vitronectin ($IC_{50}=0.83\text{ }\mu\text{M}$) thereby inhibiting tumor cell-induced angiogenesis *in vivo*. Inhibits transcription in *C. albicans*. Inhibits microbiological growth in beer. Active against soil borne pathogens.

Citations: 11

[View Online »](#)

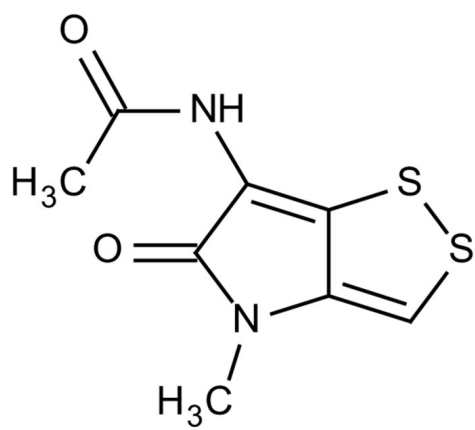
Ordering Information

[Order Online »](#)

BML-CT107-0001	1mg
BML-CT107-0005	5mg

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

Alternative Name 6-(Acetamido)-4-methyl-1,2-dithiolo[4,3-b]-pyrrol-5(4H)-one

Appearance Yellow crystalline solid.

CAS 87-11-6

Couple Target RNA polymerase

Couple Type Inhibitor

Formula $C_8H_8N_2O_2S_2$

MW 228.3

Purity ≥95% (HPLC, TLC)

Solubility Soluble in DMSO (5mg/ml), very slightly soluble in water

Source Isolated from *Streptomyces luteosporus*



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)