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 μM) 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 »

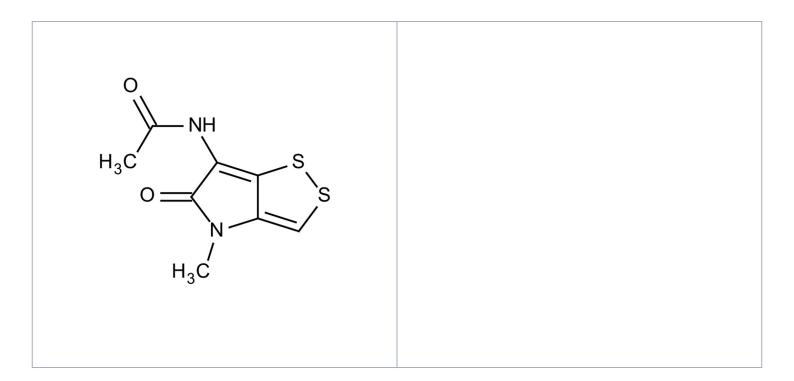
Ordering Information

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 luteosporeus

Last modified: May 29, 2024

