

# Cytochalasin B

Ultra-pure; inhibits actin polymerization

Cytochalasin B is a cell-permeable fungal toxin which binds to the barbed end of actin filaments inhibiting both the association and dissociation of subunits. It causes the disruption of actin filaments and inhibition of actin polymerization. It also inhibits monosaccharide transport across cell membranes. inhibits glucose transport and platelet aggregation. Blocks adenosine-induced apoptotic body formation without affecting activation of endogenous ADP-ribosylation in leukemia HL-60 cells.

Citations: 10

[View Online »](#)

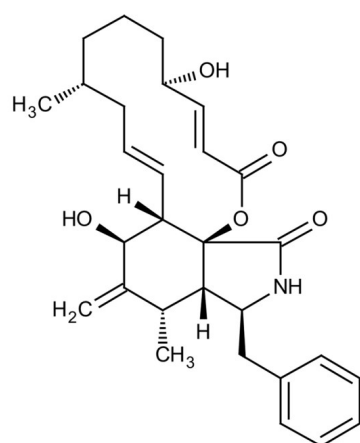
## Ordering Information

[Order Online »](#)

BML-T108-0005	5mg
---------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	White to off-white solid.
CAS	14930-96-2
Formula	$C_{29}H_{37}NO_5$
MW	479.6
Purity	≥98% (HPLC, TLC)
Solubility	Soluble in methanol or DMSO (20mg/ml).
Source	Isolated from <i>Drechslera dematoidea</i>



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)