

# Blebbistatin(+/-)

## Inhibitor of non-muscle myosin II ATPase

Blebbistatin (so named because of its ability to block cell blebbing) is a rapid, selective, and high-affinity ( $IC_{50}$  approx. 4  $\mu$ M) inhibitor of non-muscle myosin II ATPase activity. During cell division, blebbistatin inhibits contraction of the cleavage furrow without disrupting mitosis. Blebbistatin impedes cell migration and cytokinesis. Blebbistatin also reduces infiltration of inflammatory cells, inhibits pancreatic adenocarcinoma cellular invasiveness, PDGF-stimulated fibroblast-collagen matrix contraction, and diminishes harm to vascular system in models of arterial thrombosis *in vivo*.

Citations: 25

[View Online »](#)

## Ordering Information

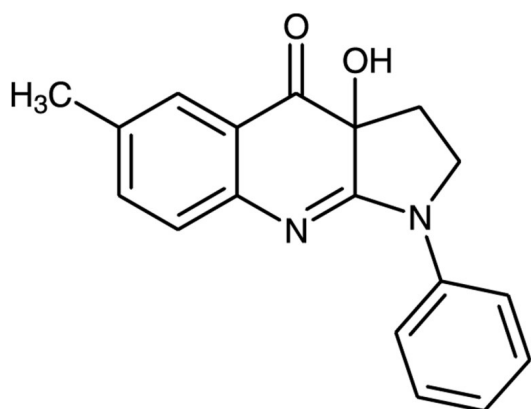
[Order Online »](#)

BML-EI315-0005	5mg
BML-EI315-0025	25mg

## Manuals, SDS & CofA

[View Online »](#)

- Rapid and selective inhibitor of Myosin II ATPase
- Used in cardiovascular research
- Highly cited



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Exposure to blue light (i.e. 450-490nm) may lead to the degradation and inactivation of blebbistatin via cytotoxic intermediates, which may be a source of concern when used in fluorescence applications.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	(±)-1,2,3,3a-Tetrahydro-3a-hydroxy-6-methyl-1-phenyl-4H-pyrrolo[2,3-b]quinolin-4-one, (±)-Blebbistatin
Appearance	Bright yellow solid.
CAS	674289-55-5
Couple Target	Myosin
Couple Type	Inhibitor
Formula	$C_{18}H_{16}N_2O_2$
Identity	Identity determined by NMR.
MW	292.3
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO (25mg/ml).

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)