

# Cyclopamine

## Blocks Hedgehog signaling

Cyclopamine is a naturally occurring steroidal alkaloid that causes cyclopia by blocking sonic hedgehog signaling. It specifically inhibits the hedgehog (Hh) pathway by interaction with the Hh signaling protein smoothened. Cyclopamine blocks the oncogenic effects of mutations of Ptch in fibroblasts and inhibits the growth of cells lacking Ptch function. It arrested cell cycle at G0/G1 and induced apoptosis in several small-cell lung cancer cell lines at 5  $\mu$ M. Cyclopamine also induces expression of the transcription factor pdx-1, which stimulates insulin secretion in differentiation protocols for mouse embryonic stem cells

Citations: 24

[View Online »](#)

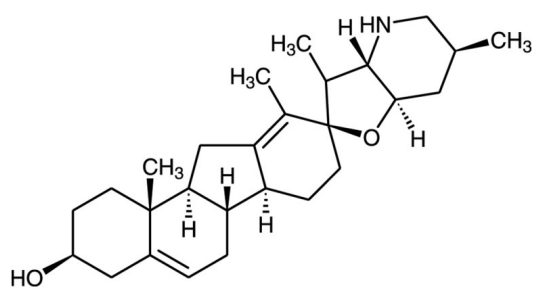
## Ordering Information

[Order Online »](#)

BML-GR334-0001	1mg
BML-GR334-0005	5mg

## 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.
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 powder
CAS	4449-51-8
Formula	$C_{27}H_{41}NO_2$
MW	411.6
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol (20mg/ml), DMSO (4mg/ml), dimethyl formamide (10mg/ml), methanol (7mg/ml). Insoluble in water.



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)