

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 μ 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 »](#)

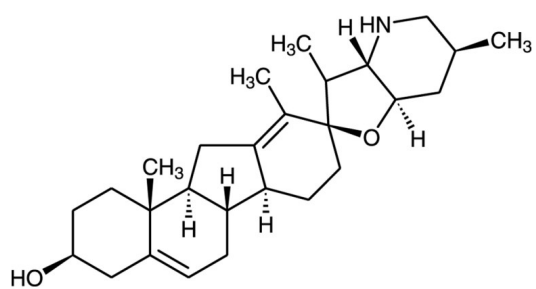
Ordering Information

[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.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com