

# Ophiobolin A

Calmodulin antagonist. PDE inhibitor.

Cell permeable antagonist of calmodulin. Inhibits  $\text{Ca}^{2+}$ /calmodulin-dependent phosphodiesterase. Inhibits P-glycoprotein-mediated transport. Exhibits antibacterial, antitumor and nematocidal activities. Phytotoxic.

Citations: 12

[View Online »](#)

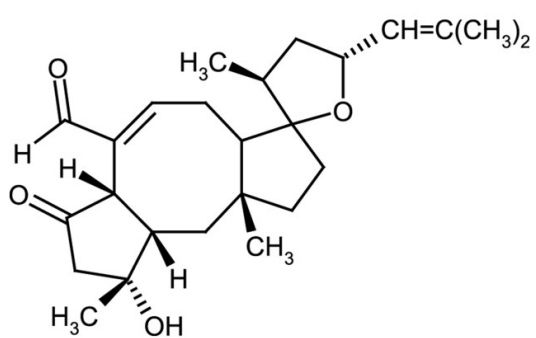
## Ordering Information

[Order Online »](#)

|                  |     |
|------------------|-----|
| ALX-270-109-M001 | 1mg |
|------------------|-----|

## 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

**Alternative Name** Cochliobolin A, Ophiobalin

**Appearance** White to off-white solid.

**CAS** 4611-05-6

**Couple Target** PDE

**Couple Type** Inhibitor

**Formula**  $C_{25}H_{36}O_4$

**MW** 400.6

**Purity** ≥95% (HPLC)

**RTECS** RL1576000

**Solubility** Soluble in dimethyl formamide, DMSO, 100% ethanol or methanol.

**Source** Isolated from *Bipolaris leersia* MST-FP107.

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)