

# Patulin

Inhibitor of protein farnesylation in a cell free assay. Inhibits incorporation of tritiated mevalonate into proteins in whole cells. Mycotoxin with anti-bacterial, potassium uptake inhibitory and possibly carcinogenic activities. Known contaminant of spoiled apples.

Citations: 8

[View Online »](#)

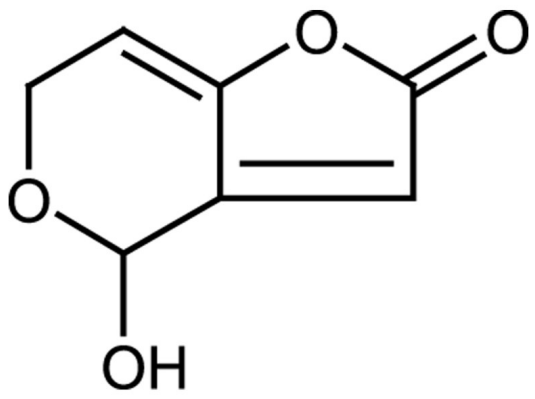
## Ordering Information

[Order Online »](#)

ALX-270-111-M005	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

Alternative Name	4-Hydroxy-4H-furo[3,2-c]pyran-2(6H)-one, Clairformin, Clavacin, Clavatin, Claviformin, Leucopin, Mycoin C3, Mycosin, Penicidin, Terinin
Appearance	White crystalline solid.
CAS	149-29-1
Couple Target	Farnesyltransferase
Couple Type	Inhibitor
Formula	$C_7H_6O_4$
MI	14: 7053
MW	154.1
Purity	≥98% (HPLC, TLC)
RTECS	LV2625000
Solubility	Soluble in DMSO, methanol, ethanol or water.
Source	Isolated from <i>Penicillium expansum</i> .



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

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

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

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

Last modified: May 29, 2024