

# Zearalenone

## Mycotoxin

Zearalenone (ZEN) is an estrogenic mycotoxin in animals and a phytohormone in plants. Inducer of sister chromatid exchange and chromosomal aberration. Acts as a protonophoric uncoupler in plant mitochondria.

Citations: 5

[View Online »](#)

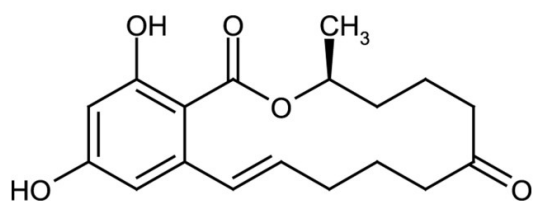
## Ordering Information

[Order Online »](#)

|                  |      |
|------------------|------|
| ALX-630-105-M010 | 10mg |
|------------------|------|

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Benzoxacyclotetradecin-1,7(8H)-dione, ZEN

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

**CAS** 17924-92-4

**Formula**  $C_{18}H_{22}O_5$

**MI** **14:** 10116

**MW** 318.4

**Purity** ≥99% (HPLC, TLC)

**RTECS** DM2550000

**Solubility** Soluble in methanol (5mg/ml).

**Source** Isolated from *Giberella zeae*

**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)