

# T-2 Toxin

## Monoamine oxidase inhibitor

Mycotoxin. Induces DNA damage and apoptosis. Increases blood-brain barrier permeability and inhibits monoamine oxidase activity in the brain.

*Not for sale in Turkey and South Africa.*

*May require a license for import, please [contact us](#) for more information.*

Citations: 12

[View Online »](#)

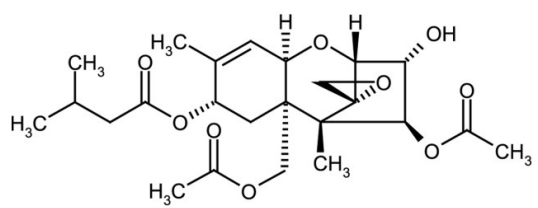
## Ordering Information

[Order Online »](#)

|                  |     |
|------------------|-----|
| ALX-630-101-M001 | 1mg |
| ALX-630-101-M005 | 5mg |

## 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** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 4β,15-Diacetoxy-3α-hydroxy-8α-(3-methylbutyryloxy)-12,13-epoxytrichothec-9-ene

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

**CAS** 21259-20-1

**Couple Target** Monoamine oxidase

**Couple Type** Inhibitor

**Formula** C<sub>24</sub>H<sub>34</sub>O<sub>9</sub>

**MW** 466.5

**Purity** ≥98% (TLC)

**RTECS** YD0100000

**Solubility** Soluble in methylene chloride (5mg/ml) or chloroform.

**Source** Isolated from *Fusarium Sporotrichoides*.

**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-  
de@enzolifesciences.com](mailto:info-de@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)