

HT-2 Toxin

Mycotoxin

Trichothecene group mycotoxin. Induces apoptosis

Not for sale in Turkey and South Africa.

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

Citations: 4

[View Online »](#)

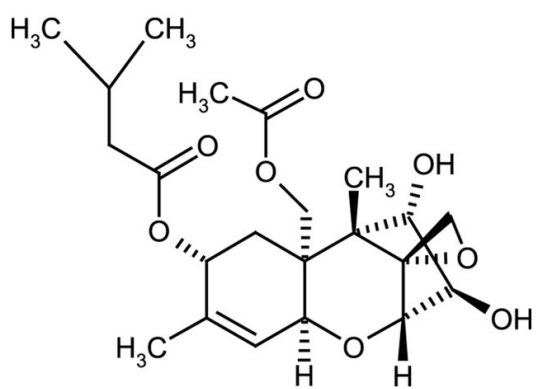
Ordering Information

[Order Online »](#)

ALX-630-113-M001	1mg
ALX-630-113-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 Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 15-Acetoxy-3 α ,4 β -dihydroxy-8 α -(3-methylbutyryloxy)-12,13-epoxytrichothec-9-ene

Appearance White to slight yellow solid.

CAS 26934-87-2

Formula C₂₂H₃₂O₈

MW 424.5

Purity ≥98% (TLC)

RTECS YD0050000

Solubility Soluble in 100% ethanol, DMSO or ethyl acetate. Slightly soluble in petroleum ether; poorly soluble in water.

Source Semisynthetic, derived from T2 toxin from *Fusarium tricinctum*.

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-
be@enzolifesciences.com](mailto:info-be@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)