

Deoxynivalenol

Mycotoxin

Mycotoxin found in cereals. Binds to the ribosome and inhibits protein synthesis.

Emetic and refusal activity in pig.

Citations: 16

[View Online »](#)

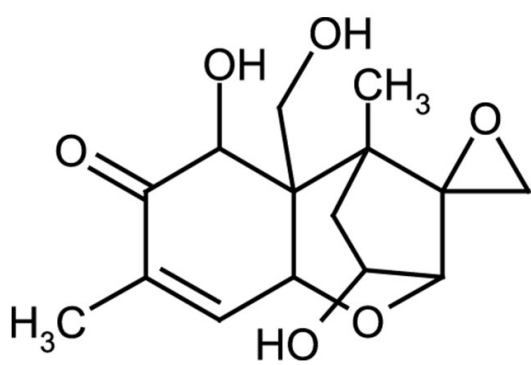
Ordering Information

[Order Online »](#)

ALX-630-115-M001	1mg
ALX-630-115-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 +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name DON, 3 α ,7 α ,15-Trihydroxy-12,13-epoxytrichothec-9-en-8-one, Vomitoxin

Appearance White to off-white powder.

CAS 51481-10-8

Formula C₁₅H₂₀O₆

MI **14:** 10032

MW 296.3

Purity ≥98% (TLC, HPLC)

RTECS YD0167000

Solubility Soluble in methanol or ethyl acetate; slightly soluble in water.

Source Isolated from *Fusarium* sp.

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)