

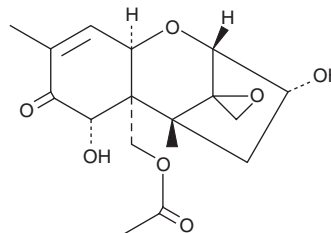
PRODUCT INFORMATION



15-Acetyldeoxy Nivalenol

Item No. 11430

CAS Registry No.: 88337-96-6
Formal Name: 15-(acetyloxy)-12,13-epoxy-3 α ,7 α -dihydroxy-trichothec-9-en-8-one
Synonyms: 15-Acetyl Vomitoxin, 15-AcDON, Deoxynivalenol 15-acetate, DON 15-acetate
MF: C₁₇H₂₂O₇
FW: 338.4
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

15-Acetyldeoxy nivalenol is supplied as a solid. A stock solution may be made by dissolving the 15-acetyldeoxy nivalenol in the solvent of choice, which should be purged with an inert gas. 15-Acetyldeoxy nivalenol is soluble in ethanol and chloroform.

Description

15-Acetyldeoxy nivalenol is a trichothecene mycotoxin produced by certain species of the fungus *Fusarium*, particularly those found on cereal crops.¹ It is an acetylated derivative of 4-deoxy nivalenol (DON; Item No. 11428) that can be converted to DON in the body. 15-Acetyldeoxy nivalenol induces vomiting and anorexia in livestock and experimental animals and its toxicity compared with DON is dependent upon the route of administration.^{2,3} It also changes brain bioamine neurotransmitter levels in pigs and chickens.⁴

References

1. Ajandouz, E.H., Berdah, S., Moutardier, V., *et al.* Hydrolytic fate of 3/15-acetyldeoxynivalenol in humans: Specific deacetylation by the small intestine and liver revealed using *in vitro* and *ex vivo* approaches. *Toxins (Basel)* **8(8)**, E232 (2016).
2. Pestka, J.J., Lin, W.-S., and Miller, E.R. Emetic activity of the trichothecene 15-acetyldeoxynivalenol in swine. *Food Chem. Toxicol.* **25(11)**, 855-888 (1987).
3. Payros, D., Alazzane-Kpembé, I., Pierron, A., *et al.* Toxicology of deoxynivalenol and its acetylated and modified forms. *Arch. Toxicol.* **90(12)**, 2931-2957 (2016).
4. Bonnet, M.S., Roux, J., Mounien, L., *et al.* Advances in deoxynivalenol toxicity mechanisms: The brain as a target. *Toxins (Basel)* **4(11)**, 1120-1138 (2012).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/07/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM