

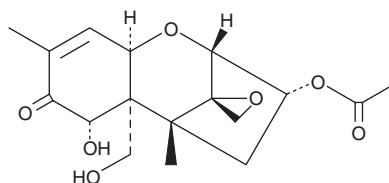
PRODUCT INFORMATION



3-Acetyldeoxy Nivalenol

Item No. 11429

CAS Registry No.: 50722-38-8
Formal Name: (3 α ,7 α)-3-(acetyloxy)-12,13-epoxy-7,15-dihydroxy-trichothec-9-en-8-one
Synonyms: 3-AcDON, DON 3-acetate, deoxy Nivalenol 3-acetate, NSC 267036
MF: C₁₇H₂₂O₇
FW: 338.4
Purity: $\geq 99\%$
UV/Vis.: λ_{max} : 218 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

3-Acetyldeoxy nivalenol is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-acetyldeoxy nivalenol in the solvent of choice, which should be purged with an inert gas. 3-Acetyldeoxy nivalenol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 3-acetyldeoxy nivalenol in ethanol and DMF is approximately 30 mg/ml and approximately 25 mg/ml in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 3-acetyldeoxy nivalenol can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 3-acetyldeoxy nivalenol in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

3-Acetyldeoxy nivalenol is a mycotoxin that has been found in *F. graminearum*.¹ *In vivo*, 3-acetyldeoxy nivalenol (40 mg/kg) induces duodenal and splenic cell necrosis, as well as lethality (LD₅₀ = 70 mg/kg) in mice.²

References

1. Jiao, F., Kawakami, A., and Nakajima, T. Effects of different carbon sources on trichothecene production and Tri gene expression by *Fusarium graminearum* in liquid culture. *FEMS Microbiol. Lett.* **285**(2), 212-219 (2008).
2. Schiefer, H.B., Nicholson, S., Kasali, O.B., *et al.* Pathology of acute 3-acetyldeoxynivalenol toxicity in mice. *Can. J. Comp. Med.* **49**(3), 315-318 (1985).

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, 10/20/2022

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