

PRODUCT INFORMATION



3-Acetyldeoxy Nivalenol-¹³C₁₇ Item No. 31258

CAS Registry No.: 1217476-81-7

Formal Name: (2S,2'R,3'R,5'R,5a'R,6'S,9a'R)-6'-hydroxy-5a'-(hydroxymethyl-¹³C)-5',8'-di(methyl-¹³C)-7'-oxo-2',3',4',5',5a',6',7',9a'-octahydrospiro[oxirane-2,10'-[2,5]methanobenzo[b]oxepin]-3'-yl-2,2',3,3',4',5',5a',6',7',8',9',9a'-¹³C₁₂ acetate-¹³C₂
Synonyms: 3-AcDON-¹³C₁₇, DON 3-acetate-¹³C₁₇, Deoxy Nivalenol-3-acetate

MF: [¹³C]₁₇H₂₂O₇

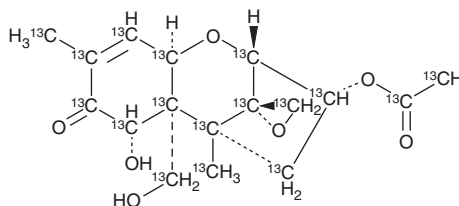
FW: 355.2

Purity: ≥98%

Supplied as: A 25 µg/ml solution in acetonitrile

Storage: -20°C

Stability: ≥4 years



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

Description

3-Acetyldeoxy nivalenol-¹³C₁₇ is intended for use as an internal standard for the quantification of 3-acetyldeoxy nivalenol (Item No. 11429) by GC- or LC-MS. 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, 02/09/2026

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