

PRODUCT INFORMATION



diacetoxy Scirpenol-¹³C₁₉ Item No. 31287

Formal Name: ((2S,2'R,3'R,4'S,5'S,5a'R,9a'R)-4'-(acetoxymethanobenzo[b]oxepin)-5a'(6'H)-yl-2,2',3,3',4',5',5a',6',7',8',9',9a'-¹³C₁₂)methyl-¹³C₂ acetate-¹³C₂

Synonyms: Anguidin-¹³C₁₉, DAS-¹³C₁₉, Scirpenetriol 4,15-diacetate-¹³C₁₉

MF: [¹³C]₁₉H₂₆O₇

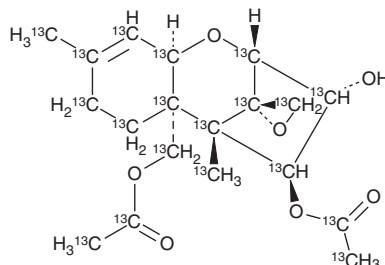
FW: 385.3

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

diacetoxy Scirpenol-¹³C₁₉ is intended for use as an internal standard for the quantification of diacetoxy scirpenol (Item No. 11427) by GC- or LC-MS. diacetoxy Scirpenol is a trichothecene mycotoxin originally isolated from *F. equiseti*.¹ It is toxic to pigs (LD₅₀ = 0.38 mg/kg, i.v.) and lethal to brine shrimp (LC₅₀ = 0.39 µg/ml).^{2,3} diacetoxy Scirpenol (1 µg/ml) suppresses phagocytosis by, and superoxide production in, isolated mouse macrophages in an *S. cerevisiae* model of phagocytosis.⁴ It is toxic to barley, wheat, and sorghum seedlings (LD₅₀s = 1.26, 3.98, and 10 mg/L, respectively).⁴ It has been found as a contaminant in grasses intended to be used as animal feed.⁵

References

1. Dawkins, A.W., Grove, J.F., and Tidd, B.K. Diacetoxyscirpenol and some related compounds. *Chem. Commun. (Lond.)* **2**, 27-28 (1965).
2. Weaver, G.A., Kurtz, H.J., Mirocha, C.J., et al. Acute toxicity of the mycotoxin diacetoxyscirpenol in swine. *Can. Vet. J.* **19(10)**, 267-271 (1978).
3. Perkowski, J., Foremska, E., and Latus-Ziętkiewicz, D. The yields of diacetoxyscirpenol produced by *Fusarium sambucinum* cultures isolated from potato tubers and their toxicity to brine shrimps (*Artemia salina*). *Mycotoxin Res.* **5(2)**, 61-67 (1989).
4. Ayral, A.M., Dubech, N., Le Bars, J., et al. In vitro effect of diacetoxyscirpenol and deoxynivalenol on microbicidal activity of murine peritoneal macrophages. *Mycopathologia* **120(2)**, 121-127 (1992).
4. Hasan, H.A. Phytotoxicity of pathogenic fungi and their mycotoxins to cereal seedling viability. *Mycopathologia* **148(3)**, 149-155 (1999).
5. Kononenko, G.P., Burkin, A.A., Gavrilova, O.P., et al. Fungal species and multiple mycotoxin contamination of cultivated forage crops. *Agric. Food Sci.* **24(4)**, 323-330 (2015).

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