

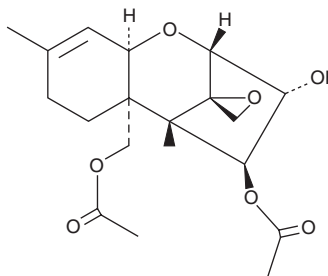
PRODUCT INFORMATION



diacetoxy Scirpenol

Item No. 11427

CAS Registry No.: 2270-40-8
Formal Name: (3 α ,4 β)-12,13-epoxy-4,15-diacetate-trichothec-9-ene-3,4,15-triol
Synonyms: Anguidin, DAS, NSC 141537, NSC 177378, Scirpenetriol 4,15-diacetate
MF: C₁₉H₂₆O₇
FW: 366.4
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

diacetoxy Scirpenol (DAS) is supplied as a crystalline solid. A stock solution may be made by dissolving the DAS in the solvent of choice. DAS is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of DAS in these solvents is approximately 20 mg/ml.

DAS is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, DAS should first be dissolved in DMF and then diluted with the aqueous buffer of choice. DAS has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

DAS is a potent, trichothecene mycotoxin produced by certain *Fusarium* strains, which play an important role as plant pathogens, causing a wide range of diseases.¹ It is toxic to fungi, plants, animals, and various mammalian cell cultures, inhibiting *de novo* protein synthesis.² In a biological activity assay, DAS was shown to inhibit expression of the *Arabidopsis* MAMP-responsive reporter gene *WRKY29p::GUS* with an IC₅₀ value of 50 nM, thereby suppressing a rapid cell death immune response important for plant defense.³

References

1. Bauer, J., Bollwahn, W., Gareis, M., *et al.* Kinetic profiles of diacetoxyscirpenol and two of its metabolites in blood serum of pigs. *Appl. Environ. Microbiol.* **49(4)**, 842-845 (1985).
2. Shams, M., Mitterbauer, R., Corradini, R., *et al.* Isolation and characterization of a new less-toxic derivative of the *Fusarium* mycotoxin diacetoxyscirpenol after thermal treatment. *J. Agric. Food Chem.* **59(17)**, 9709-9714 (2011).
3. Serrano, M., Hubert, D.A., Dangl, J.L., *et al.* A chemical screen for suppressors of the *avrRpm1*-RPM1-dependent hypersensitive cell death response in *Arabidopsis thaliana*. *Planta.* **231(5)**, 1013-1023 (2010).

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, 02/28/2019

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