

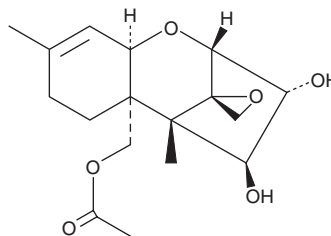
PRODUCT INFORMATION



15-acetoxy Scirpenol

Item No. 11431

CAS Registry No.: 2623-22-5
Formal Name: 12,13-epoxy-trichothec-9-ene-3 α ,4 β ,15-triol, 15-acetate
Synonyms: 4-Deacetylanguidin, NSC 267030
MF: C₁₇H₂₄O₆
FW: 324.4
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-acetoxy Scirpenol is supplied as a powder. A stock solution may be made by dissolving the 15-acetoxy scirpenol in the solvent of choice, which should be purged with an inert gas. 15-acetoxy Scirpenol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 15-acetoxy scirpenol in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

15-acetoxy Scirpenol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 15-acetoxy scirpenol should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 15-acetoxy Scirpenol 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

15-acetoxy Scirpenol is a trichothecene mycotoxin that is produced predominantly by members of the *Fusarium* genus.¹ It can induce apoptotic cell death in human Jurkat T cells.²

References

1. Peraica, M., Radić, B., Lucić, A., *et al.* Toxic effects of mycotoxins in humans. *Bull. World Health Organ.* **77(9)**, 754-766 (1999).
2. Lee, D.H., Park, T. and Kim, H.W. Induction of apoptosis by disturbing mitochondrial-membrane potential and cleaving PARP in Jurkat T cells through treatment with acetoxyscirpenol mycotoxins. *Biol. Pharm. Bull.* **29(4)**, 648-654 (2006).

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, 07/21/2025

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