

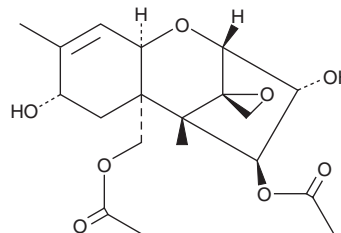
# PRODUCT INFORMATION



## Neosolaniol

Item No. 11436

**CAS Registry No.:** 36519-25-2  
**Formal Name:** (3 $\alpha$ ,4 $\beta$ ,8 $\alpha$ )-12,13-epoxy-trichothec-9-ene-3,4,8,15-tetrol 4,15-diacetate  
**Synonyms:** NSC 197212, 8-hydroxydiacetoxyl Scirpenol  
**MF:** C<sub>19</sub>H<sub>26</sub>O<sub>8</sub>  
**FW:** 382.4  
**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

Neosolaniol is supplied as a powder. A stock solution may be made by dissolving the neosolaniol in the solvent of choice, which should be purged with an inert gas. Neosolaniol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of neosolaniol in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Neosolaniol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, neosolaniol should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Neosolaniol 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

Neosolaniol is a type A trichothecene mycotoxin.<sup>1</sup> It is relatively less toxic than other type A trichothecenes, including T-2 toxin (Item No. 11444).<sup>2,3</sup>

### References

1. McCormick, S. P., Stanley, A. M., Stover, N. A., *et al.* Trichothecenes: From simple to complex mycotoxins. *Toxins (Basel)* **3**(7), 802-814 (2011).
2. Engler, K. H., Coker, R. D., and Evans, I. H. A colorimetric technique for detecting trichothecenes and assessing relative potencies. *Appl. Environ. Microbiol.* **65**(5), 1854-1857 (1999).
3. Islam, Z., Nagase, M., Ota, A., *et al.* Structure-function relationship of T-2 toxin and its metabolites in inducing thymic apoptosis *in vivo* in mice. *Biosci. Biotechnol. Biochem.* **62**(8), 1492-1497 (1998).

#### 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