

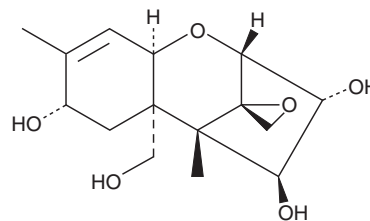
# PRODUCT INFORMATION



## T-2 Tetraol

Item No. 20432

**CAS Registry No.:** 34114-99-3  
**Formal Name:** 12,13-epoxy-trichothec-9-ene-3 $\alpha$ ,4 $\beta$ ,8 $\alpha$ ,15-tetrol  
**Synonym:** T-2 Toxin Tetraol  
**MF:** C<sub>15</sub>H<sub>22</sub>O<sub>6</sub>  
**FW:** 298.3  
**Purity:**  $\geq$ 98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years  
**Special Conditions:** Protect from light



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

### Laboratory Procedures

T-2 Tetraol is supplied as a solid. A stock solution may be made by dissolving the T-2 tetraol in the solvent of choice, which should be purged with an inert gas. A stock solution may be made by dissolving the T-2 tetraol in the solvent of choice. T-2 Tetraol is soluble in the organic solvent dichloromethane at a concentration of approximately 10 mg/ml.

### Description

T-2 tetraol is a trichothecene mycotoxin that has been found in *F. tricinctum* and an active metabolite of T-2 toxin (Item No. 11444).<sup>1</sup> It induces cell death (LC<sub>50</sub> = 9  $\mu$ g/ml), reduces proliferation (EC<sub>50</sub> = 0.25  $\mu$ g/ml), and decreases protein and DNA synthesis in primary human fibroblasts when used at a concentration of 10  $\mu$ g/ml. T-2 tetraol is toxic to newly hatched chicks (LD<sub>50</sub> = 33.79 mg/kg).<sup>3</sup> It reduces feed consumption and weight gain in newly hatched chicks when administered at a dose of 40 mg/kg. T-2 tetraol has been found in *Fusarium*-contaminated barley kernels.<sup>2</sup>

### References

1. Oldham, J.W., Allred, L.E., Milo, G.E., *et al.* The toxicological evaluation of the mycotoxins T-2 and T-2 tetraol using normal human fibroblasts *in vitro*. *Toxicol. Appl. Pharmacol.* **52(1)**, 159-168 (1979).
2. Perkowski, J., Kiecana, I., and Kaczmarek, Z. Natural occurrence and distribution of *Fusarium* toxins in contaminated barley cultivars. *Eur. J. Plant Pathol.* **109**, 331-339 (2003).
3. Chi, M.S., Robison, T.S., Mirocha, C.J., *et al.* Acute toxicity of 12,13-epoxytrichothecenes in one-day-old broiler chicks. *Appl. Environ. Microbiol.* **35(4)**, 636-640 (1978).

#### 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, 07/01/2025

#### 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