

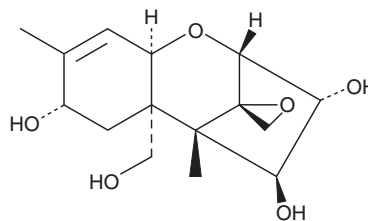
PRODUCT INFORMATION



T-2 Tetraol

Item No. 20432

CAS Registry No.: 34114-99-3
Formal Name: 12,13-epoxy-trichothec-9-ene-3 α ,4 β ,8 α ,15-tetrol
Synonym: T-2 Toxin Tetraol
MF: C₁₅H₂₂O₆
FW: 298.3
Purity: \geq 98%
Supplied as: A powder
Storage: -20°C
Stability: \geq 2 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 powder. A stock solution may be made by dissolving the T-2 tetraol in the solvent of choice. 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, which should be purged with an inert gas, at a concentration of approximately 10 mg/ml.

Description

T-2 Tetraol is a group A trichothecene mycotoxin derived by the metabolism of T-2 toxin (Item No. 11444). It is much less toxic than the parent compound and is unable to induce apoptosis in thymic cells in mice.^{1,2}

References

- 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).
- 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, 06/23/2022

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