

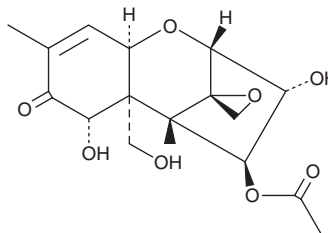
# PRODUCT INFORMATION



## Fusarenon X

Item No. 11432

**CAS Registry No.:** 23255-69-8  
**Formal Name:** 4 $\beta$ -(acetyloxy)-12,13-epoxy-3 $\alpha$ ,7 $\alpha$ ,15-trihydroxy-trichothec-9-en-8-one  
**Synonyms:** 4-acetyl Nivalenol, NSC 197211  
**MF:** C<sub>17</sub>H<sub>22</sub>O<sub>8</sub>  
**FW:** 354.4  
**Purity:**  $\geq$ 99%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 2 years



**Special Conditions:** Protect from light; Hygroscopic

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

### Laboratory Procedures

Fusarenon X is supplied as a crystalline solid. A stock solution may be made by dissolving the fusarenon X in the solvent of choice. Fusarenon X is soluble in the organic solvent dichloromethane, which should be purged with an inert gas.

### Description

Fusarenon X is a type B trichothecene mycotoxin typically derived from *Fusarium* species. It is primarily found in contaminated cereals.<sup>1</sup> Fusarenon X inhibits protein synthesis, which leads to disruption of DNA synthesis.<sup>1</sup> As this occurs in actively proliferating cells, fusarenon X causes immunosuppression, intestinal malabsorption, developmental toxicity, and genotoxicity.<sup>1</sup> Fusarenon X can be metabolized to fusarenon X-glucoside by infected plants.<sup>2</sup>

### References

1. Aupanun, S., Poapolathep, S., Girogi, M. *et al.* An overview of the toxicology and toxicokinetics of fusarenon-X, a type B trichothecene mycotoxin. *J. Vet. Med. Sci.* (2016).
2. Berthiller, F., Crews, C., Dall'Asta, C., *et al.* Masked mycotoxins: A review. *Mol. Nutr. Food Res.* **57**, 165-186 (2013).

#### 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, 10/26/2017

#### 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