

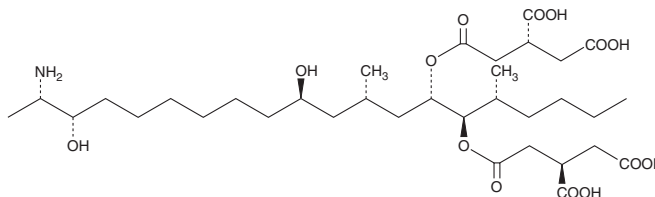
# PRODUCT INFORMATION



## Fumonisin B<sub>3</sub> Item No. 20434

**CAS Registry No.:** 1422359-85-0  
**Formal Name:** (2R,2'R)-1,1'-[(1S,2R)-1-[(2S,4R,11S,12S)-12-amino-4,11-dihydroxy-2-methyltridecyl]-2-[(1R)-1-methylpentyl]-1,2-ethanediyl]ester, 1,2,3-propanetricarboxylic acid

**Synonym:** FB<sub>3</sub>  
**MF:** C<sub>34</sub>H<sub>59</sub>NO<sub>14</sub>  
**FW:** 705.8  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥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

Fumonisin B<sub>3</sub> is supplied as a solid. A stock solution may be made by dissolving the fumonisin B<sub>3</sub> in the solvent of choice. Fumonisin B<sub>3</sub> is soluble in methanol at a concentration of approximately 5 mg/ml.

### Description

Fumonisin B<sub>3</sub> is a mycotoxin produced from *F. moniliforme*, a prevalent fungus of corn and other grains.<sup>1</sup> Outbreaks of food poisoning in livestock and humans following the consumption of *Fusarium* infested corn are caused by fumonisins.<sup>2</sup> Fumonisin function as inhibitors of ceramide synthase (sphingosine N-acyltransferase).<sup>3</sup>

### References

1. Newkirk, D.K., Benson, R.W., Howard, P.C., *et al.* On-line immunoaffinity capture, coupled with HPLC and electrospray ionization mass spectrometry, for automated determination of fumonisins. *J. Agric. Food Chem.* **46**, 1677-1688 (1998).
2. Gelderblom, W.C.A., Jaskiewicz, K., Marasas, W.F.O., *et al.* Fumonisin - novel mycotoxins with cancer-promoting activity produced by *Fusarium moniliforme*. *Appl. Environ. Microbiol.* **54**, 1806-1811 (1988).
3. Wang, E., Norred, W.P., Bacon, C.W., *et al.* Inhibition of sphingolipid biosynthesis by fumonisins. Implications for diseases associated with *Fusarium moniliforme*. *J. Biol. Chem.* **266**, 14486-14490 (1991).

**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