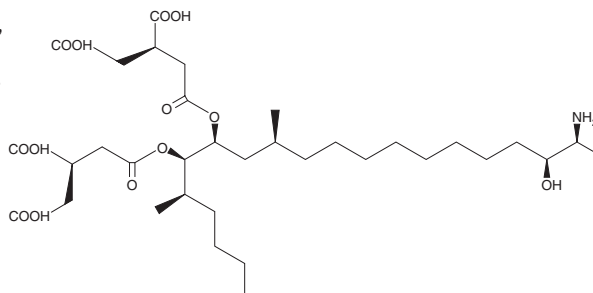


PRODUCT INFORMATION



Fumonisin B₄ Item No. 35683

CAS Registry No.: 136379-60-7
Formal Name: (2R,2'R)-1,2,3-propanetricarboxylic acid, 1,1'-[(1S,2R)-1-[(2S,11S,12S)-12-amino-11-hydroxy-2-methyltridecyl]-2-[(1R)-1-methylpentyl]-1,2-ethanediyl] ester
Synonym: FB₄
MF: C₃₄H₅₉NO₁₃
FW: 689.8
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Fumonisin B₄ (FB₄) is supplied as a solid. A stock solution may be made by dissolving the FB₄ in the solvent of choice, which should be purged with an inert gas. FB₄ is soluble in methanol and DMSO.

Description

FB₄ is a mycotoxin that has been found in *F. verticillioides*.¹ It has also been found in corn and poultry feed.^{2,3}

References

1. Gelderblom, W.C.A., Sewram, V., Shephard, G.S., *et al.* Structure and natural occurrence of stereoisomers of the fumonisin B series mycotoxins. *J. Agric. Food Chem.* **55(11)**, 4388-4394 (2007).
2. Abbas, H.K., Williams, W.P., Windham, G.L., *et al.* Aflatoxin and fumonisin contamination of commercial corn (*Zea mays*) hybrids in Mississippi. *J. Agric. Food Chem.* **50(18)**, 5246-5254 (2002).
3. Ezekiel, C.N., Bandyopadhyay, R., Sulyok, M., *et al.* Fungal and bacterial metabolites in commercial poultry feed from Nigeria. *Food Addit. Contam. Part A Chem. Anal. Control Expo. Risk Assess.* **29(8)**, 1288-1299 (2012).

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, 01/02/2024

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