

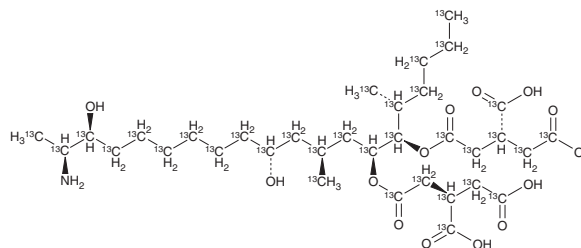
PRODUCT INFORMATION



Fumonisin B₃-¹³C₃₄ Item No. 31274

Formal Name: (2R,2'R)-2,2'-(((5R,6R,7S,9S,11R,18S,19S)-19-amino-11,18-dihydroxy-5,9-di(methyl-¹³C)icosane-6,7-diyl-1,2,3,4,5,6,7,8,9,10,11,12,13,14,15,16,17,18,19,20-¹³C₂₀)bis(oxy))bis(2-oxoethane-2,1-diyl-1,2-¹³C₂) disuccinic-¹³C₄ acid

Synonym: FB₃-¹³C₃₄
MF: [¹³C]₃₄H₅₉NO₁₄
FW: 739.6
Purity: ≥95%
Supplied as: A 10 µg/ml 1:1 solution in acetonitrile:water
Storage: -20°C
Stability: ≥4 years



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

Description

FB₃-¹³C₃₄ is intended for use as an internal standard for the quantification of fumonisin B₃ (Item No. 20434) by GC- or LC-MS. FB₃ is a mycotoxin that has been found in *Fusarium*.¹ It is cytotoxic to primary rat hepatocytes when used at concentrations of 125, 250, and 500 µM.² Dietary administration of FB₃ (0.05%) for 14 or 21 days induces hepatocyte nodules, a marker of cancer initiation, in rats.² FB₃ has been detected in corn and corn-based foods and livestock feeds.³

References

1. Rheeder, J.P., Marasas, W.F.O., and Vismer, H.F. Production of fumonisin analogs by *Fusarium* species. *Appl. Environ. Microbiol.* **68**(5), 2101-2105 (2002).
2. Gelderblom, W.C.A., Cawood, M.E., Snyman, S.D., et al. Structure-activity relationships of fumonisins in short-term carcinogenesis and cytotoxicity assays. *Food Chem. Toxicol.* **31**(6), 407-414 (1993).
3. Bullerman, L.B. Occurrence of *Fusarium* and fumonisins on food grains and in foods. *Fumonisin in food*. Jackson, L.S., DeVries, J.W., Bullerman, L.B., editors, 1st edition, Springer (1996).

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, 02/09/2026

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