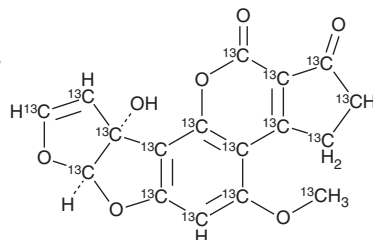


PRODUCT INFORMATION



Aflatoxin M₁-¹³C₁₇ Item No. 31264

CAS Registry No.: 2707441-97-0
Formal Name: (6aR,9aR)-9a-hydroxy-4-(methoxy-¹³C)-2,3,6a,9a-tetrahydrocyclopenta[c]furo[3',2':4,5]furo[2,3-h]chromene-1,11-dione-1,2,3,3a,3b,4,5,5a,6a,8,9,9a,9b,9c,11,11a-¹³C₁₆
Synonym: AFM₁-¹³C₁₇
MF: [¹³C]₁₇H₁₂O₇
FW: 345.1
Purity: ≥98%
Supplied as: A 0.5 µg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥4 years



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

Description

AFM₁-¹³C₁₇ is intended for use as an internal standard for the quantification of AFM₁ (Item No. 11297) by GC- or LC-MS. AFM₁ is an active metabolite of aflatoxin B₁ (Item No. 11293).¹ It induces hepatocyte hyperplasia and hepatocellular carcinoma tumor formation in rats when administered at a dose of 1 mg/animal. AFM₁ (6 mg/kg) increases hepatic total lipid levels and inhibits hepatic mitochondrial and cytoplasmic protein synthesis in rats.²

References

1. Wogan, G.N. and Paglialunga, S. Carcinogenicity of synthetic aflatoxin M₁ in rats. *Food Cosmet. Toxicol.* **12**(3), 381-385 (1974).
2. Pai, M.R., Bai, N.J., and Venkitasubramanian, T.A. Aflatoxin induced inhibition of protein synthesis. *Toxicon* **16**(3), 283-287 (1978).

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