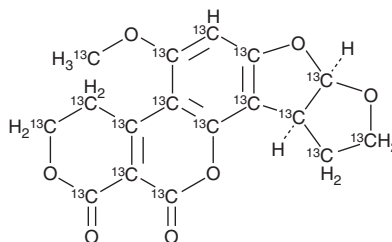


PRODUCT INFORMATION



Aflatoxin G₂-¹³C₁₇
Item No. 31263

CAS Registry No.: 1217462-49-1
Formal Name: (7aR,10aS)-5-(methoxy-¹³C)-3,4,7a,9,10,10a-hexahydro-1H,12H-furo[3',2':4,5]furo[2,3-h]pyrano[3,4-c]chromene-1,12-dione-1,3,4,4a,4b,5,6,6a,7a,9,10,10a,10b,10c,12,12a-¹³C₁₆
Synonym: Dihydroaflatoxin G₁-¹³C₁₇
MF: [¹³C]₁₇H₁₄O₇
FW: 347.2
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

Aflatoxin G₂-¹³C₁₇ is intended for use as an internal standard for the quantification of aflatoxin G₂ (Item No. 11296) by GC- or LC-MS. Aflatoxin G₂ is a mycotoxin that has been found in *Aspergillus*.¹ It is lethal to ducklings (LD₅₀ = 2.83 mg/kg) but is non-toxic to rats when administered at a dose of 200 mg/kg.²

References

1. Bennett, J.W. and Klich, M. Mycotoxins. *Clin. Microbiol. Rev.* **16**(3), 497-516 (2003).
2. Wogan, G.N., Edwards, G.S., and Newberne, P.M. Structure-activity relationships in toxicity and carcinogenicity of aflatoxins and analogs. *Cancer Res.* **31**(12), 1936-1942 (1971).

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