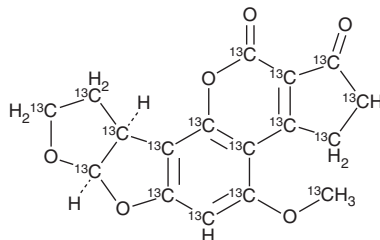


PRODUCT INFORMATION



Aflatoxin B₂-¹³C₁₇ Item No. 31261

CAS Registry No.: 1217470-98-8
Formal Name: (6aR,9aS)-4-(methoxy-¹³C)-2,3,6a,8,9,9a-hexahydrocyclopenta[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
Synonyms: AFB₂-¹³C₁₇, Dihydroaflatoxin B₁-¹³C₁₇
MF: [¹³C]₁₇H₁₄O₆
FW: 331.2
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 B₂-¹³C₁₇ (AFB₂-¹³C₁₇) is intended for use as an internal standard for the quantification of AFB₂ (Item No. 11294) by GC- or LC-MS. AFB₂ is a mycotoxin that has been found in *A. terricola*.¹ It induces hepatic autophagy and apoptosis in broiler chickens when administered at doses of 0.2, 0.4, and 0.8 mg/kg.² AFB₂ (0.5 and 1 mg/animal) also induces parenchymal cell hyperplasia in rats.³

References

1. Moubasher, A.H., el-Kady, I.A., and Shoriet, A. Toxigenic *Aspergilli* isolated from different sources in Egypt. *Ann. Nutr. Aliment.* **31**(4-6), 607-615 (1977).
2. Chen, B., Li, D., Li, M., *et al.* Induction of mitochondria-mediated apoptosis and PI3K/Akt/mTOR-mediated autophagy by aflatoxin B₂ in hepatocytes of broilers. *Oncotarget* **7**(51), 84989-84998 (2016).
3. 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