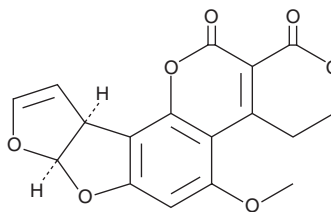


PRODUCT INFORMATION



Aflatoxin G₁ Item No. 11295

CAS Registry No.: 1165-39-5
Formal Name: (7aR,10aS)-3,4,7a,10a-tetrahydro-5-methoxy-1H,12H-furo[3',2':4,5]furo[2,3-h]pyrano[3,4-c][1]benzopyran-1,12-dione
MF: C₁₇H₁₂O₇
FW: 328.3
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

Aflatoxin G₁ is supplied as a powder. A stock solution may be made by dissolving the aflatoxin G₁ in the solvent of choice. Aflatoxin G₁ is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide, which should be purged with an inert gas.

Description

Aflatoxin G₁ is a mycotoxin that has been found in *A. terricola*.¹ *In vivo*, aflatoxin G₁ is lethal to ducklings (LD₅₀ = 1.18 mg/kg).² It induces hepatocellular carcinoma tumor formation and lethality in rats when administered at doses of 1.4 and 3 mg/animal, respectively. Aflatoxin G₁ also inhibits liver and kidney succinate dehydrogenase and fumarase, as well as kidney cytochrome oxidase, NADH oxidase, α-glycerophosphate dehydrogenase, isocitrate dehydrogenase, and malate dehydrogenase 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. 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).
3. Bai, N.J., Pai, M.R., and Venkitasubramanian, T.A. Mitochondrial function in aflatoxin toxicity. *Indian J. Biochem. Biophys.* **14**(4), 347-349 (1977).

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, 07/21/2025

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