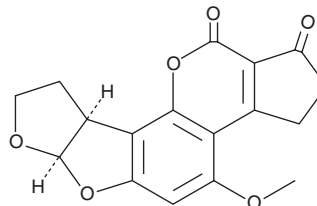


PRODUCT INFORMATION



Aflatoxin B₂ Item No. 11294

CAS Registry No.: 7220-81-7
Formal Name: 2,3,6aR,8,9,9aS-hexahydro-4-methoxy-cyclopenta[c]furo[3',2':4,5]furo[2,3-h][1]benzopyran-1,11-dione
Synonyms: AFB₂, Dihydroaflatoxin B₁
MF: C₁₇H₁₄O₆
FW: 314.3
Purity: ≥98%
UV/Vis.: λ_{max}: 202, 220, 265, 363 nm
Supplied as: A crystalline 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 B₂ is supplied as a crystalline solid. A stock solution may be made by dissolving the aflatoxin B₂ in the solvent of choice which should be purged with an inert gas. Aflatoxin B₂ is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of aflatoxin B₂ in DMSO is approximately 12 mg/ml and approximately 20 mg/ml in DMF.

Aflatoxin B₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, aflatoxin B₂ should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Aflatoxin B₂ has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Aflatoxin B₂ (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

Reference

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, 12/16/2022

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