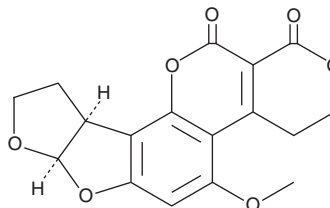


PRODUCT INFORMATION



Aflatoxin G₂ Item No. 11296

CAS Registry No.: 7241-98-7
Formal Name: (7aR,10aS)-3,4,7a,9,10,10a-hexahydro-5-methoxy-1H,12H-furo[3',2':4,5]furo[2,3-h]pyrano[3,4-c][1]benzopyran-1,12-dione
Synonym: Dihydroaflatoxin G₁
MF: C₁₇H₁₄O₇
FW: 330.3
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 G₂ is supplied as a solid. A stock solution may be made by dissolving the aflatoxin G₂ in the solvent of choice. Aflatoxin G₂ is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide, which should be purged with an inert gas.

Description

Aflatoxin G₂ is a food contaminant produced by various species of the common soil fungus, *Aspergillus* and is associated with toxicity and hepatocarcinogenicity in human and animal populations.^{1,2} This compound may be used as an internal standard when monitoring foodstuffs for aflatoxin contamination.

References

1. Bennett, J.W. and Klich, M. Mycotoxins. *Clin. Microbiol. Rev.* **16**(3), 497-516 (2003).
2. Wong, J.J. and Hsieh, D.P. Mutagenicity of aflatoxins related to their metabolism and carcinogenic potential. *Proc. Natl. Acad. Sci. USA* **73**(7), 2241-2244 (1976).

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, 06/30/2025

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