

PRODUCT INFORMATION



Aflatoxin R₀ Item No. 20437

CAS Registry No.: 29611-03-8
Formal Name: 2,3,6aR,9aS-tetrahydro-1S-hydroxy-4-methoxy-cyclopenta[c]furo[3',2':4,5]furo[2,3-h][1]benzopyran-11(1H)-one

Synonym: Aflatoxicol

MF: C₁₇H₁₄O₆

FW: 314.3

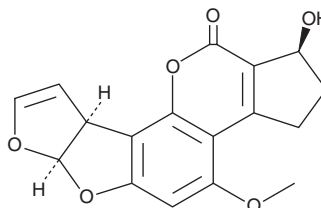
Purity: ≥98%

Supplied as: A powder

Storage: -20°C

Stability: ≥2 years

Item Origin: Fungi/*Aspergillus flavus*



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Aflatoxin R₀ is supplied as a powder. A stock solution may be made by dissolving the aflatoxin R₀ in the solvent of choice. Aflatoxin R₀ is soluble in methanol at a concentration of approximately 1 mg/ml.

Description

Aflatoxin R₀ is a metabolite of the carcinogenic mycotoxin aflatoxin B₁ (Item No. 11293). Aflatoxin R₀ is produced by metabolism in animals as well as in fungi and can be reverted to the parent compound in fungi.^{1,2} Aflatoxin R₀ is highly mutagenic when tested in the Ames' *in vitro* microbial detection system for carcinogens.²

References

1. Nakazato, Morozumi, M., S., Saito, K., *et al.* Interconversion of aflatoxin B₁ and aflatoxicol by several fungi. *Appl. Environ. Microbiol* **56(5)**, 1465-1470 (1990).
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, 09/04/2019

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