

PRODUCT INFORMATION

Ochratoxin A

Item No. 11439

CAS Registry No.: 303-47-9

Formal Name: N-[[[(3R)-5-chloro-3,4-dihydro-8-hydroxy-3-methyl-1-oxo-1H-2-benzopyran-7-yl] carbonyl]-L-phenylalanine

Synonyms: NSC 201422, NSC 221991, OTA

MF: $C_{20}H_{18}ClNO_6$

FW: 403.8

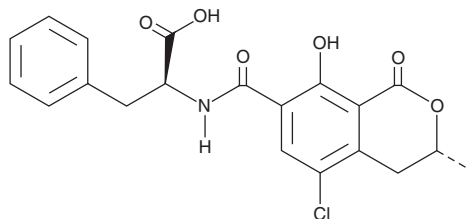
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 213, 331 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 2 years



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

Laboratory Procedures

Ochratoxin A (OTA) is supplied as a crystalline solid. A stock solution may be made by dissolving the OTA in the solvent of choice. OTA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of OTA in these solvents is approximately 50, 16, and 14 mg/ml, respectively.

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

Description

Ochratoxin A (OTA) is a mycotoxin that has been found in *Penicillium* and is an active metabolite of OTC (Item No. 20183).¹⁻³ It is formed from OTC *in vivo*, however, OTC can also be formed from OTA by gut microbiota.^{1,4} OTA (120 μ g/kg) increases renal lipid peroxide levels and the number of apoptotic cells, as well as reduces renal superoxide dismutase (SOD) activity in rats.² Topical application of OTA (80 μ g/animal) induces DNA damage, cell cycle arrest at the G₀/G₁ phase, and apoptosis in mouse skin cells.³ It also initiates tumor formation in a two-stage mouse skin tumorigenesis model. OTA has been found in food products and poultry feed.^{5,6}

References

1. Fuchs, R., Hult, K., Peraica, M., *et al.* Conversion of ochratoxin C into ochratoxin A *in vivo*. *Appl. Environ. Microbiol.* **48**(1), 41-42 (1984).
2. Petrik, J., Zanić-Grubišić, T., Barisić, K., *et al.* Apoptosis and oxidative stress induced by ochratoxin A in rat kidney. *Arch. Toxicol.* **77**(12), 685-693 (2003).
3. Kumar, R., Ansari, K.M., Chaudhari, B.P., *et al.* Topical application of ochratoxin A causes DNA damage and tumor initiation in mouse skin. *PLoS One* **7**(10), (2012).
4. Galtier, P. and Alvinerie, M. *In vitro* transformation of ochratoxin A by animal microbial floras. *Ann. Rech. Vet.* **7**(1), 91-98 (1976).
5. Al-Taher, F., Cappozzo, J., Zweigenbaum, J., *et al.* Detection and quantitation of mycotoxins in infant cereals in the U.S. market by LC-MS/MS using a stable isotope dilution assay. *Food Control* **72**(Part A), 27-35 (2017).
6. Ezekiel, C.N., Bandyopadhyay, R., Sulyok, M., *et al.* Fungal and bacterial metabolites in commercial poultry feed from Nigeria. *Food Addit. Contam. Part A Chem. Anal. Control Expo. Risk Assess.* **29**(8), 1288-1299 (2012).

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, 01/11/2024

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