

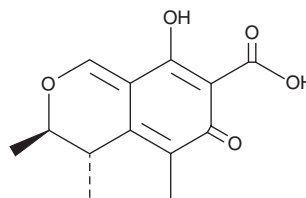
PRODUCT INFORMATION



Citrinin

Item No. 11320

CAS Registry No.: 518-75-2
Formal Name: (3R,4S)-4,6-dihydro-8-hydroxy-3,4,5-trimethyl-6-oxo-3H-2-benzopyran-7-carboxylic acid
Synonyms: NSC 186, (-)-Citrinin, CTN
MF: C₁₃H₁₄O₅
FW: 250.3
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 251, 330 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Penicillium citrinum*



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

Laboratory Procedures

Citrinin is supplied as a crystalline solid. A stock solution may be made by dissolving the citrinin in the solvent of choice, which should be purged with an inert gas. Citrinin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of citrinin in ethanol is approximately 2 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Citrinin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, citrinin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Citrinin 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

Citrinin is a mycotoxin that has been found in *Monascus* and has diverse biological activities.¹⁻⁴ It is active against *S. aureus*, methicillin-resistant *S. aureus* (MRSA), rifampicin-resistant *S. aureus*, and vancomycin-resistant *E. faecium* (MICs = 1.95, 3.9, 0.97, and 7.81 μg/ml, respectively), as well as the pathogenic yeast *C. neoformans* (MIC = 3.9 μg/ml).² It is cytotoxic to a variety of cells *in vitro*, including bovine kidney cells and mice embryonic stem cells.⁴ Citrinin (30 μM) induces reactive oxygen species (ROS) production, mitochondrial membrane potential loss, and apoptosis in HepG2 cells, effects that can be blocked by the antioxidant resveratrol.³ In contrast, citrinin reduces glutamate-induced excitotoxicity in primary rat cortical neurons at concentrations ranging from 0.1 to 1,000 nM and inhibits LPS-induced production of nitric oxide (NO) in RAW 264.7 cells at 0.625 to 40 μM.⁴ It is toxic to brine shrimp larvae (LD₅₀ = 96 μg/ml), as well as to rats and mice with oral LD₅₀ values of 50 and 87-105 mg/kg, respectively.^{2,4} It induces reproductive abnormalities in male mice and toxic effects in the liver, kidney, heart, and gastrointestinal tracts of various animals.⁴ Citrinin has been found in stored cereal grains, as well as beans, fruit, and herbs.

References

1. Blanc, P.J., Laussac, J.P., Le Bars, P., et al. *Int. J. Food Microbiol.* **27(2-3)**, 201-213 (1995).
2. Subramani, R., Kumar, R., Prasad, P., et al. *Asian Pac. J. Allergy Immunol.* **3(4)**, 291-296 (2013).
3. Chen, C.-C. and Chan, W.-H. *Int. J. Mol. Sci.* **10(8)**, 3338-3357 (2009).
4. Filho, A.R.d.S., Islam, M.T., Ali, E.S., et al. *Food Chem. Toxicol.* **110**, 130-141 (2017).

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, 08/11/2023

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