

PRODUCT INFORMATION

Ochratoxin B-¹³C₂₀

Item No. 40238

Formal Name: ((R)-8-hydroxy-3-(methyl-¹³C)-1-oxoisochromane-7-carbonyl-¹³C₉)-L-phenylalanine-¹³C₉

Synonym: OTB-¹³C₂₀

MF: [¹³C]₂₀H₁₉NO₆

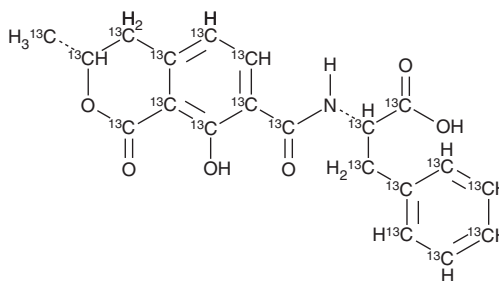
FW: 389.2

Purity: ≥98%

Supplied as: A 10 µg/ml solution in methanol

Storage: -20°C

Stability: ≥4 years



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

Description

Ochratoxin B-¹³C₂₀ (OTB-¹³C₂₀) is intended for use as an internal standard for the quantification of ochratoxin B (Item No. 16167) by GC- or LC-MS. OTB is a mycotoxin that has been found in *Aspergillus* and is a non-chlorinated derivative of OTA (Item No. 11439).¹ Unlike OTA, OTB is not genotoxic to HepG2 cells but does inhibit cell division when used at concentrations ranging from 1 to 25 µg/ml.² It induces mortality (LC₅₀ = 700 nM) and craniofacial malformations (EC₅₀ = 200 nM) in *Xenopus* embryos.³ OTB has been found in poultry feed.⁴

References

1. Varga, J., Kevei, E., Rinyu, E., *et al.* Ochratoxin production by *Aspergillus* species. *Appl. Environ. Microbiol.* **62**(12), 4461-4464 (1996).
2. Knasmüller, S., Cavin, C., Chakraborty, A., *et al.* Structurally related mycotoxins ochratoxin A, ochratoxin B, and citrinin differ in their genotoxic activities and in their mode of action in human-derived liver (HepG2) cells: Implications for risk assessment. *Nutr. Cancer* **50**(2), 190-197 (2004).
3. O'Brien, E., Prietz, A., and Dietrich, D.R. Investigation of the teratogenic potential of ochratoxin A and B using the FETAX system. *Birth Defects Res. B Dev. Reprod. Toxicol.* **74**(5), 417-423 (2005).
4. 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, 05/08/2026

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