

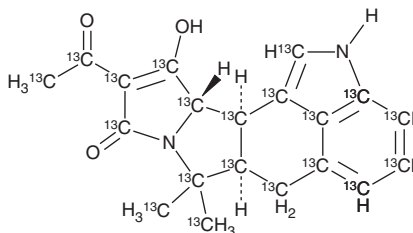
PRODUCT INFORMATION



Cyclopiazonic Acid-¹³C₂₀ Item No. 31297

Formal Name: (6aR,11aS,11bR)-10-(acetyl-¹³C₂)-11-hydroxy-7,7-di(methyl-¹³C)-2,6,6a,7,11a,11b-hexahydro-9H-pyrrolo[1',2':2,3]isoindolo[4,5,6-cd]indol-9-one-1,2a,2a1,3,4,5,5a,6,6a,7,9,10,11,11a,11b,11c-¹³C₁₆

Synonym: CPA-¹³C₂₀
MF: [¹³C]₂₀H₂₀N₂O₃
FW: 356.2
Purity: ≥98%
Supplied as: A 10 µg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥2 years



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

Description

Cyclopiazonic acid-¹³C₂₀ is intended for use as an internal standard for the quantification of cyclopiazonic acid (Item No. 11326) by GC- or LC-MS. Cyclopiazonic acid is a mycotoxin that has been found in *Penicillium* and *Aspergillus* and an inhibitor of sarcoplasmic/endoplasmic reticulum Ca²⁺-ATPase (SERCA; IC₅₀ = 0.6 µM).^{1,2} *In vivo*, cyclopiazonic acid (11-14 mg/kg) induces ptosis, hypothermia, tremor, and cachexia, as well as induces lethality (LD₅₀ = 13 mg/kg), in rats.³ It induces tachycardia and sedation in rabbits when administered at a dose of 10 mg/kg.

References

1. Seidler, N.W., Jona, I., Vegh, M., et al. Cyclopiazonic acid is a specific inhibitor of the Ca²⁺-ATPase of sarcoplasmic reticulum. *J. Biol. Chem.* **264**(30), 17816-17823 (1989).
2. Uyama, Y., Imaizumi, Y., and Watanabe, M. Effects of cyclopiazonic acid, a novel Ca²⁺-ATPase inhibitor, on contractile responses in skinned ileal smooth muscle. *Brit. J. Pharmacol.* **106**(1), 208-214 (1992).
3. Nishie, K., Cole, R.J., and Dorner, J.W. Toxicity and neuropharmacology of cyclopiazonic acid. *Food Chem. Toxicol.* **23**(9), 831-839 (1985).

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, 08/30/2023

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