

PRODUCT INFORMATION

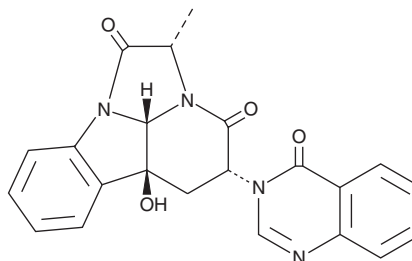


(-)-Chaetominine

Item No. 13135

CAS Registry No.: 918659-56-0
Formal Name: (2S,4R,5aS,9cS)-4,5,5a,9c-tetrahydro-5a-hydroxy-2-methyl-4-(4-oxo-3(4H)-quinazoliny)-3H-2a,9b-diazacyclopenta[jk]fluorene-1,3(2H)-dione

MF: C₂₂H₁₈N₄O₄
FW: 402.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Description

(-)-Chaetominine is a cytotoxic alkaloid originally isolated from *Chaetomium* sp. IFB-E015.¹ It inhibits the growth of K562 leukemia and SW1116 colon cancer cells (IC₅₀s = 20 and 28 nM, respectively). (-)-Chaetominine induces apoptosis of K562 cells via upregulation of the Bax/Bcl-2 ratio, decreasing mitochondrial membrane potential, inducing mitochondrial cytochrome C release, and activation of caspase-3 and caspase-9.² It also decreases doxorubicin efflux mediated by multidrug resistance-associated protein 1 (MRP1) and restores sensitivity to doxorubicin (Item No. 15007) in resistant K562 cells.³

References

1. Jiao, R.H., Xu, S., Liu, J.Y., *et al.* Chaetominine, a cytotoxic alkaloid produced by endophytic *Chaetomium* sp. IFB-E015. *Org. Lett.* **8(25)**, 5709-5712 (2006).
2. Yao, J., Jiao, R.H., Liu, C., *et al.* Assessment of the cytotoxic and apoptotic effects of chaetominine in a human leukemia cell line. *Biomol. Ther. (Seoul)* **24(2)**, 147-155 (2016).
3. Yao, J., Wei, X., and Lu, Y. Chaetominine reduces MRP1-mediated drug resistance via inhibiting PI3K/Akt/Nrf2 signaling pathway in K562/Adr human leukemia cells. *Biochem. Biophys. Res. Commun.* **473(4)**, 867-873 (2016).

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, 12/13/2022

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