

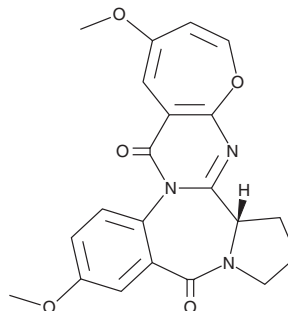
PRODUCT INFORMATION



Circumdatin A

Item No. 31436

CAS Registry No.: 223130-52-7
Formal Name: (16aS)-14,15,16,16a-tetrahydro-4,10-dimethoxy-6H,12H-oxepino[2',3':4,5]pyrimido[1,2-a]pyrrolo[2,1-c][1,4]benzodiazepine-6,12-dione
MF: C₂₁H₁₉N₃O₅
FW: 393.4
Purity: ≥95%
UV/Vis.: λ_{max}: 238 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Aspergillus* sp.



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

Laboratory Procedures

Circumdatin A is supplied as a solid. A stock solution may be made by dissolving the circumdatin A in the solvent of choice, which should be purged with an inert gas. Circumdatin A is soluble in ethanol.

Description

Circumdatin A is a benzodiazepine alkaloid fungal metabolite originally isolated from *A. ochraceus*.¹

Reference

1. Rahbaek, L., Breinholt, J., Frisvad, J.C., et al. Circumdatin A, B, and C: Three new benzodiazepine alkaloids isolated from a culture of the fungus *Aspergillus ochraceus*. *J. Org. Chem.* **64**(5), 1689-1692 (1999).

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, 08/10/2023

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