

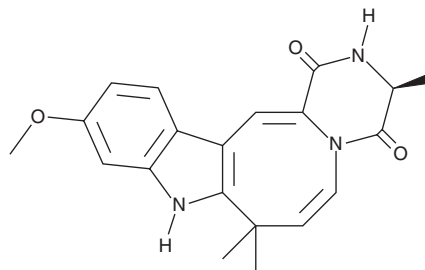
PRODUCT INFORMATION



Cycloechinulin

Item No. 25455

CAS Registry No.: 143086-29-7
Formal Name: (3S)-2,3,8,9-tetrahydro-11-methoxy-3,8,8-trimethyl-pyrazino[1':2':1,2]azocino[5,4-b]indole-1,4-dione
MF: C₂₀H₂₁N₃O₃
FW: 351.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Aspergillus ochraceus*



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

Laboratory Procedures

Cycloechinulin is supplied as a solid. A stock solution may be made by dissolving the cycloechinulin in the solvent of choice, which should be purged with an inert gas. Cycloechinulin is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Cycloechinulin is a diketopiperazine fungal metabolite originally isolated from *A. ochraceus*.¹ It reduces weight gain of corn earworms by 33% compared to controls when used at a dose of 100 ppm in the diet.

Reference

1. de Guzman, F.S., Gloer, J.B., Wicklow, D.T., *et al.* New diketopiperazine metabolites from the sclerotia of *Aspergillus ochraceus*. *J. Nat. Prod.* **55(7)**, 931-939 (1992).

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, 08/23/2023

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