

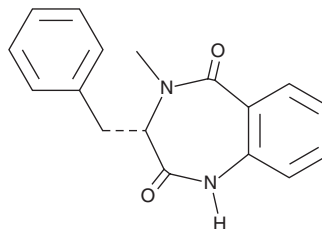
PRODUCT INFORMATION



Cyclopeptine

Item No. 23476

CAS Registry No.: 50886-63-0
Formal Name: (3S)-3,4-dihydro-4-methyl-3-(phenylmethyl)-1H-1,4-benzodiazepine-2,5-dione
MF: C₁₇H₁₆N₂O₂
FW: 280.3
Purity: ≥98%
UV/Vis.: λ_{max}: 217 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Penicillium* sp.



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

Laboratory Procedures

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

Description

Cyclopeptine is an intermediate in the synthesis of benzodiazepine alkaloids in *Penicillium*.¹ It is formed when cyclopeptine synthase catalyzes the activation of anthranilic acid (Item No. 18708) and phenylalanine, synthesizes two peptide bonds, and completes methylation using S-adenosylmethionine.

Reference

1. Luckner, M. Alkaloid biosynthesis in *Penicillium cyclopoum* - does it reflect general features of secondary metabolism? *J. Nat. Prod.* **43**(1), 21-40 (1980).

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, 08/09/2023

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