

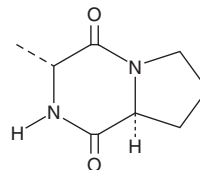
PRODUCT INFORMATION



Cyclo(D-Ala-L-Pro)

Item No. 24940

CAS Registry No.: 36238-64-9
Formal Name: (3R,8aS)-hexahydro-3-methyl-
pyrrolo[1,2-a]pyrazine-1,4-dione
MF: C₈H₁₂N₂O₂
FW: 168.2
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.

Laboratory Procedures

Cyclo(D-Ala-L-Pro) is supplied as a solid. A stock solution may be made by dissolving the cyclo(D-Ala-L-Pro) in the solvent of choice. Cyclo(D-Ala-L-Pro) is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide, which should be purged with an inert gas.

Description

Cyclo(D-Ala-L-Pro) is a diketopiperazine fungal metabolite originally isolated from *P. terrestre*.¹

Reference

1. Lui, H., Chen, J., Deng, Z.-W., *et al.* Secondary metabolites from marine derived fungus *Penicillium terrestre*. *J. Chin. Pharma. Sci.* **19(6)**, 482-486 (2010).

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, 12/19/2022

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