

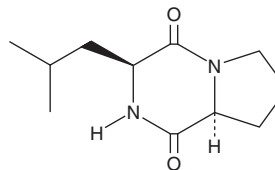
PRODUCT INFORMATION



Cyclo(L-Leu-L-Pro)

Item No. 24941

CAS Registry No.: 2873-36-1
Formal Name: (3S,8aS)-hexahydro-3-(2-methylpropyl)-pyrrolo[1,2-a]pyrazine-1,4-dione
Synonym: Maculosin 6
MF: C₁₁H₁₈N₂O₂
FW: 210.3
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(L-Leu-L-Pro) is supplied as a solid. A stock solution may be made by dissolving the cyclo(L-Leu-L-Pro) in the solvent of choice. Cyclo(L-Leu-L-Pro) is soluble in the organic solvent methanol, which should be purged with an inert gas, at a concentration of approximately 50 mg/ml.

Description

Cyclo(L-Leu-L-Pro) is a diketopiperazine metabolite that has been isolated from various bacterial and fungal species including *Streptomyces*.¹ It is active against twelve strains of vancomycin-resistant enterococci (VRE) with MIC values of 12.5 µg/ml for *E. faecalis* strains K-99-34, K-00-184, and K-00-221.¹ It also inhibits growth of K562, HL-60, and U937 leukemia cells in a concentration-dependent manner when used at concentrations of 1-500 µg/ml. Cyclo(L-Leu-L-Pro) also has antifouling activity, inhibiting attachment of *B. amphitrite* larva with an EC₅₀ value of 0.15 mM.²

References

1. Rhee, K.H. Isolation and characterization of *Streptomyces* sp KH-614 producing anti-VRE (vancomycin-resistant enterococci) antibiotics. *J. Gen. Appl. Microbiol.* **48(6)**, 321-327 (2002).
2. Li, X., Dobretsov, S., Xu, Y., et al. Antifouling diketopiperazines produced by a deep-sea bacterium, *Streptomyces fungicidicus*. *Biofouling* **22(3-4)**, 201-208 (2006).

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, 11/02/2022

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