

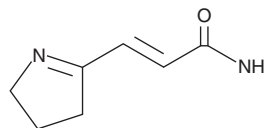
PRODUCT INFORMATION



Cyclamidomycin

Item No. 36387

CAS Registry No.: 35663-85-5
Formal Name: (2E)-3-(3,4-dihydro-2H-pyrrol-5-yl)-2-propenamide
Synonyms: Desdanine, NSC 121234, Pyracrimycin A
MF: C₇H₁₀N₂O
FW: 138.2
Purity: ≥90%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces* sp. MA130-A1



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

Laboratory Procedures

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

Description

Cyclamidomycin is a bacterial metabolite that has been found in *Streptomyces* and has antibiotic activity.¹ It is active against *S. aureus*, *M. flavus*, *S. lutea*, *B. subtilis*, *E. coli*, *S. flexneri*, *S. typhosa*, *P. vulgaris*, and *K. pneumoniae* (MICs = 3.12-25 mg/ml). Cyclamidomycin inhibits *E. coli* nucleoside diphosphokinase when used at concentrations ranging from 20 to 160 µg/ml.²

References

1. Takahashi, S., Nakajima, M., Ikeda, Y., et al. A new antibiotic, cyclamidomycin. *J. Antibiot. (Tokyo)* **24**(12), 902-903 (1971).
2. Saeki, T., Hori, M., and Umezawa, H. Cyclamidomycin (desdanine), an inhibitor of nucleoside diphosphokinase of *Escherichia coli*. *J. Antibiot. (Tokyo)* **25**(6), 343-349 (1972).

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