

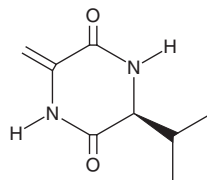
PRODUCT INFORMATION



Cyclo(Δ -Ala-L-Val)

Item No. 25739

CAS Registry No.: 25516-00-1
Formal Name: 3-methylene-6S-(1-methylethyl)-
2,5-piperazinedione
MF: C₈H₁₂N₂O₂
FW: 168.2
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

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

Cyclo(Δ -Ala-L-Val) is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Cyclo(Δ -Ala-L-Val) is a bacterial cyclic dipeptide characterized as a diketopiperazine.¹ It activates a LuxR-based N-acylhomoserine lactone (AHL) *E. coli* biosensor and blocks activation of the biosensor by the quorum sensing signal molecule N-(β -ketocaproyl)-L-homoserine lactone (3-oxo-C6-HSL; Item No. 10011207) with an IC₅₀ value of 0.8 mM. Cyclo(Δ -Ala-L-Val) (15 μ M) reduces *S. liquefaciens* colony expansion by 21%, indicating inhibition of swarming motility. It also inhibits the interaction of the kinases Ras and Raf-1 in a yeast two-hybrid assay in a concentration-dependent manner.²

References

1. Holden, M.T.G., Chhabra, S.R., de Nys, R., *et al.* Quorum-sensing cross talk: Isolation and chemical characterization of cyclic dipeptides from *Pseudomonas aeruginosa* and other gram-negative bacteria. *Mol. Microbiol.* **33**(6), 1254-1266 (1999).
2. Cheenpracha, S., Borris, R.P., Tran, T.T., *et al.* Three new amides from *Streptomyces* sp. H7372. *J. Braz. Chem. Soc.* **22**(2), 223-229 (2011).

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, 10/19/2022

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