

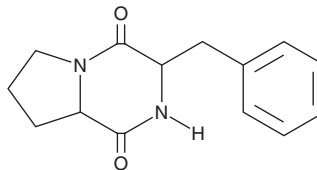
PRODUCT INFORMATION



Cyclo(Phe-Pro)

Item No. 34758

CAS Registry No.: 14705-60-3
Formal Name: hexahydro-3-(phenylmethyl)-pyrrolo[1,2-a]pyrazine-1,4-dione
Synonyms: A-64863, Cyclo(phenylalanylprolyl), cFP
MF: C₁₄H₁₆N₂O₂
FW: 244.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

Cyclo(Phe-Pro) is supplied as a solid. A stock solution may be made by dissolving the cyclo(Phe-Pro) in the solvent of choice, which should be purged with an inert gas. Cyclo(Phe-Pro) is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of cyclo(Phe-Pro) in these solvents is approximately 1 mg/ml.

Description

Cyclo(Phe-Pro) is a bacterial quorum-sensing molecule that has been found in *V. cholerae*.¹ It enhances ToxR-dependent transcription of *leuO*, a DNA binding protein, and inhibits virulence factor production in *V. cholerae*. Cyclo(Phe-Pro) is an inhibitor of the interaction between retinoic acid-inducible gene I (RIG-I) and the ubiquitin ligase TRIM25.² It increases the viral load in Huh7 cells infected with hepatitis C virus (HCV) when used at concentrations of 1, 2.5, and 5 mM. Cyclo(Phe-Pro) (50 mg/kg) reduces the expression of IFN-β induced by poly(I:C) and increases serum HCV titers in animal models of HCV infection.

References

1. Bina, X.R., Taylor, D.L., Vikram, A., et al. *Vibrio cholerae* ToxR downregulates virulence factor production in response to cyclo(Phe-Pro). *mBio* **4**(5), e00366-13 (2013).
2. Lee, W., Lee, S.-H., Kim, M., et al. *Vibrio vulnificus* quorum-sensing molecule cyclo(Phe-Pro) inhibits RIG-I-mediated antiviral innate immunity. *Nat. Commun.* **9**(1), 1606 (2018).

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, 01/03/2023

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