

PRODUCT INFORMATION

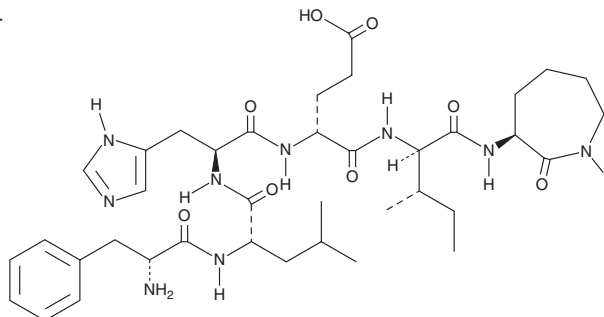


Calpinactam

Item No. 27735

CAS Registry No.: 1205538-83-5
Formal Name: D-phenylalanyl-L-leucyl-L-histidyl-D- α -glutamyl-N-[(3S)-hexahydro-2-oxo-1H-azepin-3-yl]-D-alloisoleucinamide

Synonym: FKI-4905
MF: C₃₈H₅₇N₉O₈
FW: 767.9
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Fungus/*Mortierella alpina*



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

Laboratory Procedures

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

Description

Calpinactam is a fungal metabolite originally isolated from *M. alpina* that has antimycobacterial activity.¹ It is active against *M. smegmatis* (MIC = 0.78 μ g/ml) but not a panel of 13 other bacteria and fungi. Calpinactam increases survival in a silkworm model of *M. smegmatis* infection (ED₅₀ = 5 μ g/larva).²

References

1. Koyama, N., Kojima, S., Nonaka, K., et al. Calpinactam, a new anti-mycobacterial agent, produced by *Mortierella alpina* FKI-4905. *J. Antibiot. (Tokyo)* **63**(4), 183-186 (2010).
2. Yagi, A., Uchida, R., Hamamoto, H., et al. Anti-*Mycobacterium* activity of microbial peptides in a silkworm infection model with *Mycobacterium smegmatis*. *J. Antibiot. (Tokyo)* **70**(5), 685-690 (2017).

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, 08/10/2023

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