

PRODUCT INFORMATION



Lariatins A

Item No. 28165

CAS Registry No.: 732286-09-8

Formal Name: (8→1)-lactam glycyl-L-seryl-L-glutaminy-L-leucyl-L-valyl-L-tyrosyl-L-arginyl-L-α-glutamyl-L-tryptophyl-L-valylglycyl-L-histidyl-L-seryl-L-asparaginy-L-valyl-L-isoleucyl-L-lysyl-L-proline

MF: C₉₄H₁₄₃N₂₇O₂₅

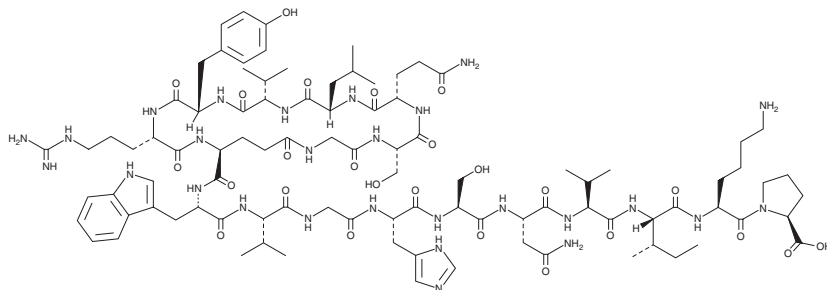
FW: 2,051.3

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

Lariatins A is supplied as a solid. A stock solution may be made by dissolving the lariatins A in the solvent of choice. Lariatins A is soluble in organic solvents such as methanol, which should be purged with an inert gas.

Lariatins A 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

Lariatins A is an antimycobacterial lasso peptide originally isolated from *R. jostii*.¹ It is active against *M. tuberculosis* and *M. smegmatis* *in vitro* (MICs = 0.39 and 0.1 µg/ml, respectively).^{1,2} Lariatins A increases survival in a silkworm model of *M. smegmatis* infection (ED₅₀ = 0.5 µg/g).

References

1. Iwatsuki, M., Uchida, R., Takakusagi, Y., *et al.* Lariatins, novel anti-mycobacterial peptides with a lasso structure, produced by *Rhodococcus jostii* K01-B0171. *J. Antibiot. (Tokyo)* **60(6)**, 357-363 (2007).
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, 12/08/2022

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