

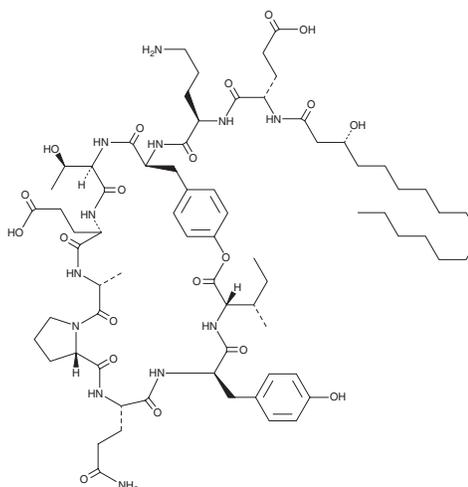
PRODUCT INFORMATION



Plipastatin A1

Item No. 38916

CAS Registry No.: 103651-09-8
Formal Name: (10→3)-lactone, N-[(3R)-3-hydroxy-1-oxohexadecyl]-L-α-glutamyl-D-ornithyl-L-tyrosyl-D-allothreonyl-L-α-glutamyl-prolyl-L-glutamyl-D-tyrosyl-L-isoleucine
Synonyms: Fengycin IX, SNA-60-367-3
MF: C₇₂H₁₁₀N₁₂O₂₀
FW: 1,463.7
Purity: ≥90%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Bacillus* sp.



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

Laboratory Procedures

Plipastatin A1 is supplied as a solid. A stock solution may be made by dissolving the plipastatin A1 in the solvent of choice. Plipastatin A1 is soluble in organic solvents such as methanol and DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Plipastatin A1 is a lipopeptide that has been found in *B. cereus* and has enzyme inhibitory and immunosuppressant activities.¹ It inhibits phospholipase A₂ (PLA₂), PLC, and PLD (IC₅₀s = 2.9, 1.3, and 1.4 μM, respectively). Plipastatin A1 (100 μg/ml) inhibits LPS-induced B cell blastogenesis and T cell blastogenesis induced by concanavalin A (Item No. 14951). *In vivo*, plipastatin A1 (0.1 mg/animal) inhibits delayed-type hypersensitivity in mice.

Reference

1. Umezawa, H., Ayoagi, T., Nishikiori, T., *et al.* Plipastatins: New inhibitors of phospholipase A₂, produced by *Bacillus cereus* BMG302-ff67. I. Taxonomy, production, isolation and preliminary characterization. *J. Antibiot. (Tokyo)* **39**(6), 737-744 (1986).

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, 10/11/2023

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