

# PRODUCT INFORMATION



## A-54556B

Item No. 17090

**CAS Registry No.:** 1629166-56-8  
**Formal Name:** N-[(2E,4E,6E)-1-oxo-2,4,6-octatrien-1-yl]-L-phenylalanyl-L-seryl-L-prolyl-N-methyl-L-alanyl-L-alanyl-L-proline, (6→2)-lactone

**Synonym:** A-54556 Factor B

**MF:** C<sub>37</sub>H<sub>48</sub>N<sub>6</sub>O<sub>8</sub>

**FW:** 704.8

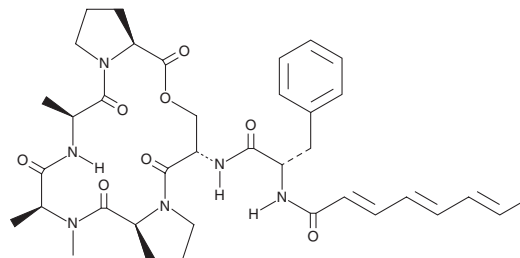
**Purity:** ≥95%

**Supplied as:** A powder

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** *Streptomyces* sp.



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

### Laboratory Procedures

A-54556B is supplied as a powder. A stock solution may be made by dissolving the A-54556B in the solvent of choice. A-54556B is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide, which should be purged with an inert gas.

### Description

A-54556B is a natural acyldepsipeptide (ADEP) antibiotic isolated from the fermentation broth of *S. hawaiiensis*.<sup>1</sup> Like the closely related ADEP, A-54556A (Item No. 15305), A-54556B is effective against Gram-positive bacteria, including penicillin-resistant *S. aureus*.<sup>1</sup> A-54556B is also effective against many strains of anaerobic bacteria.<sup>1</sup>

### Reference

1. Michel, K.H. and Kastner, R.E. A54556 antibiotics and process for production thereof. *Eli Lilly and Company*, 4,492,650 (1985).

#### 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, 10/18/2022

#### 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