

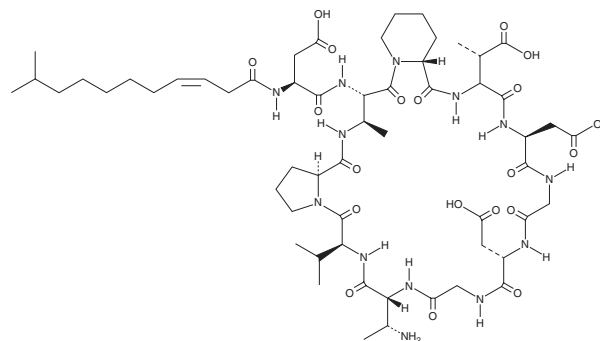
PRODUCT INFORMATION



Aspartocin D

Item No. 23143

CAS Registry No.: 1211953-97-7
Formal Name: N-[(3Z)-10-methyl-1-oxo-3-undecen-1-yl]-L- α -aspartyl-(2S,3R)-2,3-diaminobutanoyl-(2R)-2-piperidinecarbonyl-(3S)-3-methyl-L- α -aspartyl-L- α -aspartylglycyl-L- α -aspartylglycyl-(2R,3R)-2,3-diaminobutanoyl-L-valyl-L-proline (11 \rightarrow 2)-lactam
MF: C₅₇H₈₉N₁₃O₂₀
FW: 1,276.4
Purity: \geq 85%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

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

Description

Aspartocin D is a lipopeptide antibiotic originally isolated from *S. canus* FIM0916.¹ It has activity against Gram-positive bacteria, including *B. subtilis* and *S. aureus* (MICs = 0.125 and 0.5 μ g/ml, respectively). The antibiotic activity is calcium-dependent, with MICs ranging from 0.06-4 and 0.25-8 μ g/ml for *B. subtilis* and *S. aureus*, respectively, in the presence of higher (2.5 mM) to lower (0.125 mM) calcium concentrations.

Reference

1. Yang, H.J., Huang, X.Z., Zhang, Z.L., *et al.* Two novel amphomycin analogues from *Streptomyces canus* strain FIM-0916. *Nat. Prod. Res.* **28(12)**, 861-867 (2014).

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, 12/05/2022

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