

PRODUCT INFORMATION



Bactobolin A (hydrochloride)

Item No. 38912

CAS Registry No.: 73543-68-7
Formal Name: 2S-amino-N-[(3S,4R,4aR,5R,6R)-3-(dichloromethyl)-3,4,4a,5,6,7-hexahydro-5,6,8-trihydroxy-3-methyl-1-oxo-1H-2-benzopyran-4-yl]-propanamide, monohydrochloride

MF: C₁₄H₂₀Cl₂N₂O₆ • HCl

FW: 419.7

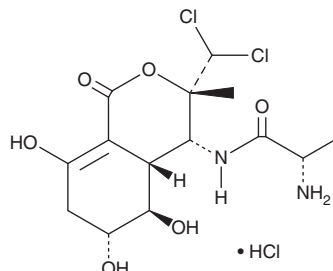
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Pseudomonas* sp.



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

Laboratory Procedures

Bactobolin A (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the bactobolin A (hydrochloride) in the solvent of choice. Bactobolin A (hydrochloride) is soluble in the organic solvent DMSO, which should be purged with an inert gas. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Bactobolin A is a microbial metabolite originally isolated from *Pseudomonas* and has antibiotic and anticancer activities.¹ It is active against *S. aureus*, *S. epidermidis*, *S. faecalis*, *B. anthracis*, *B. subtilis*, *E. coli*, *S. typhi*, and *S. dysenteriae* (MICs = 0.1-12.5 µg/ml). *In vivo*, bactobolin A (0.25-4 mg/kg) increases survival in an L1210 murine leukemia model.

Reference

1. Ezaki, N., Miyadoh, S., Hisamatsu, T., *et al.* BN-183B, a new antitumor antibiotic produced by *Pseudomonas*. Taxonomy, isolation, physico-chemical and biological properties. *J. Antibiot. (Tokyo)* **33(2)**, 213-220 (1980).

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