

PRODUCT INFORMATION



Bacitracin A (technical grade)

Item No. 36252

CAS Registry No.: 22601-59-8

Formal Name: N-[[[(4R)-2-[(1S,2S)-1-amino-2-methylbutyl]-4,5-dihydro-4-thiazolyl]carbonyl]-L-leucyl-D- α -glutamyl-L-isoleucyl-L-lysyl-D-ornithyl-L-isoleucyl-D-phenylalanyl-L-histidyl-D- α -aspartyl-L-asparagine, (10 $\rightarrow$ 4)-lactam

Synonym:

NSC 45737

MF: $C_{66}H_{103}N_{17}O_{16}S$

FW: 1,422.7

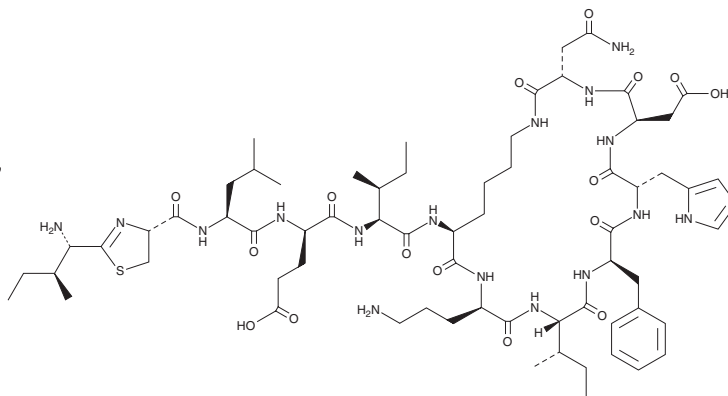
Purity: $\geq 60\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years

Item Origin: Bacterium/*Bacillus licheniformis*



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

Laboratory Procedures

Bacitracin A (technical grade) is supplied as a solid. A stock solution may be made by dissolving the bacitracin A (technical grade) in the solvent of choice, which should be purged with an inert gas. Bacitracin A (technical grade) is slightly soluble in methanol (sonicated) and DMSO.

Description

Bacitracin A is a polypeptide antibiotic that has been found in *B. licheniformis* and the main active component of the bacitracin family of polypeptide antibiotics.¹ It is active against *M. lysodeikticus* when used at a concentration of 0.1 μM and forms a complex with divalent cations and the cell membrane component C55-isoprenyl pyrophosphate.²

References

1. Newton, G.G.F. and Abraham, E.P. Observations on the nature of bacitracin A. *Biochem. J.* **53**(4), 604-613 (1953).
2. Storm, D.R. and Strominger, J.L. Complex formation between bacitracin peptides and isoprenyl pyrophosphates. The specificity of lipid-peptide interactions. *J. Biol. Chem.* **248**(11), 3940-3945 (1973).

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, 08/25/2023

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