

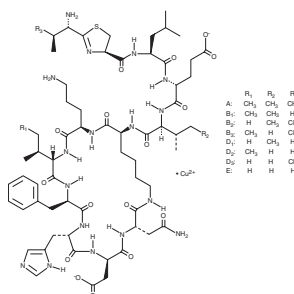
PRODUCT INFORMATION



Bacitracin Complex

Item No. 30189

MF: $C_{66}H_{101}N_{17}O_{16}S \cdot Cu$ (for A)
FW: 1,484.2
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years
Item Origin: Bacterium/*Bacillus subtilis*



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

Laboratory Procedures

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

Description

Bacitracin complex is a bacitracin (Item No. 21212) antibiotic complex composed of cyclic dipeptides that has been found in *Bacillus*.¹ It is metal-dependent and is complexed to a divalent metal cation, such as Cu^{2+} , Ni^{2+} , Co^{2+} , Zn^{2+} , or Mn^{2+} , for its activity.^{1,2} Bacitracin complex is a mixture of bacitracin polypeptides in complex with copper.

References

1. Piacham, T., Isarankura-Na-Ayudhya, C., Nantasenamat, C., *et al.* Metalloantibiotic Mn(II)-bacitracin complex mimicking manganese superoxide dismutase. *Biochem. Biophys. Res. Commun.* **341**(4), 925-930 (2006).
2. Rodríguez-Urbe, N.A., Pérez-González, R., Domínguez-Mendoza, B.E., *et al.* Metallobacitracins: Affinity and structural study in aqueous solution. *Inorg. Chem. Commun.* **104**, 61-70 (2019).

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, 10/27/2022

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