

PRODUCT INFORMATION



Bactenecin (bovine) (trifluoroacetate salt)

Item No. 34534

Formal Name: L-arginyl-L-leucyl-L-cysteinyl-L-arginyl-L-isoleucyl-L-valyl-L-valyl-L-isoleucyl-L-arginyl-L-valyl-L-cysteinyl-L-arginine, cyclic (3→11)-disulfide, trifluoroacetate salt

Synonyms: H-Arg-Leu-Cys-Arg-Ile-Val-Val-Ile-Arg-Val-Cys-Arg-OH, RLCRIVVIRVCR-OH

MF: C₆₃H₁₁₈N₂₄O₁₃S₂ • XCF₃COOH

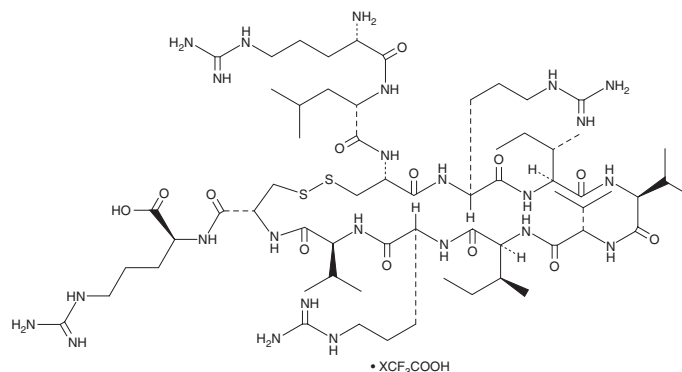
FW: 1,483.9

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

Bactenecin (bovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the bactenecin (bovine) (trifluoroacetate salt) in water. The solubility of bactenecin (bovine) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Bactenecin is a cyclic cationic peptide with antimicrobial activity.¹ It is found in bovine neutrophils. Bactenecin is active against the bacteria *E. coli* and *S. aureus* (IC₅₀s = 8 and 32 mg/ml, respectively) and the fungus *C. albicans* (MIC = 64 mg/ml).²

References

1. Romeo, D., Skerlavaj, B., Bolognesi, M., *et al.* Structure and bactericidal activity of an antibiotic dodecapeptide purified from bovine neutrophils. *J. Biol. Chem.* **263**(20), 9573-9575 (1988).
2. Hancock, R.E.W. Cationic peptides: Effectors in innate immunity and novel antimicrobials. *Lancet Infect. Dis.* **1**(3), 156-164 (2001).

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, 11/22/2022

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