

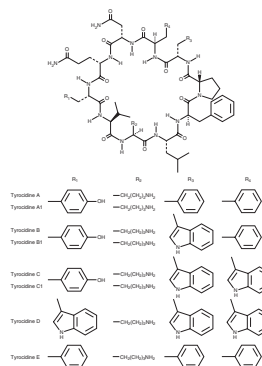
PRODUCT INFORMATION



Tyrocidine Complex

Item No. 25655

CAS Registry No.: 8011-61-8
MF: $C_{66}H_{87}N_{13}O_{13}$ (for A)
FW: 1,269.7
Purity: $\geq 95\%$
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

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

Description

Tyrocidine complex is a mixture of cyclic decapeptides originally isolated from *B. brevis*.^{1,2} It is active against Gram-positive and Gram-negative bacteria.¹ It forms a complex with single- and double-stranded DNA to inhibit RNA synthesis.³

References

1. Dubos, R.J. Studies on bactericidal agent extracted from a soil bacillus: I. Preparation of the agent. Its activity *in vitro*. *J. Exp. Med.* **70**(1), 1-10 (1939).
2. Hotchkiss, R.D. The chemical nature of gramicidin and tyrocidine. *J. Biol. Chem.* **141**, 171-185 (1941).
3. Schazschneider, B., Ristow, H., and Kleinkauf, H. Interaction between the antibiotic tyrocidine and DNA *in vitro*. *Nature* **249**(459), 757-759 (1974).

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, 01/03/2023

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