

# PRODUCT INFORMATION

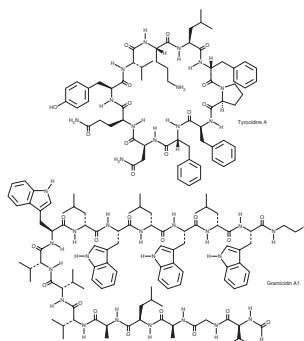
## Tyrothricin

Item No. 22177

**CAS Registry No.:** 1404-88-2  
**MF:**  $C_{66}H_{87}N_{13}O_{13}$  (for Tyrocidine A)  
**FW:** 1,270.5  
**Composite Purity:**  $\geq 95\%$  composite purity (based on 80% tyrocidine and 20% gramicidin complexes)

**Supplied as:** A solid  
**Storage:**  $-20^{\circ}\text{C}$   
**Stability:**  $\geq 4$  years

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



### Laboratory Procedures

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

### Description

Tyrothricin is a polypeptide antibiotic mixture produced by *B. brevis*.<sup>1,2</sup> Tyrothricin is a mixture of tyrocidine and gramicidin that has rapid *in vitro* bactericidal activity against *T. microdentium*, *T. macrodentium*, *T. vincentii*, *T. buccalis*, *T. mucosum*, *T. vaginalis*, *T. cruzi*, and *L. tropica*.<sup>2</sup> Preclinical studies demonstrate that intravenous use of tyrothricin leads to hemolysis-induced lethality ( $\text{LD}_{50}\text{s} = 0.4$  and  $0.3$  mg/kg, in dogs and mice, respectively) and damages liver, kidney, and olfactory tissues. Formulations containing tyrothricin are used topically to treat bacterial skin infections.<sup>1,2</sup>

### References

1. Dubos, R.J. and Hotchkiss, R.D. The production of bactericidal substances by aerobic sporulating bacilli. *J. Exp. Med.* **73**(5), 629-640 (1941).
2. Lask, S.A. Tyrothricin as an antibiotic. *Arch. Surg.* **56**(4), 475-483 (1948).

#### 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, 12/02/2022

#### 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