

# PRODUCT INFORMATION



## Turonicin A

Item No. 45511

**CAS Registry No.:** 3059457-35-8  
**Formal Name:** (2E,4E,6E,8E,12E,14S,15S,16S,17R,18E,20E,22E,24E,26E,29S,30S,33R,35S,37S,38E,41R,43S,45R,47S,48E,51S,52E,55R)-58-guanidino-15,17,29,33,35,37,41,43,45,47,51,55-dodecahydroxy-2,14,16,30-tetramethyl-31-oxooctapentacontane-2,4,6,8,12,18,20,22,24,26,38,48,52-tridecaenoic acid

**MF:** C<sub>63</sub>H<sub>99</sub>N<sub>3</sub>O<sub>15</sub>

**FW:** 1,138.5

**Purity:** ≥95%

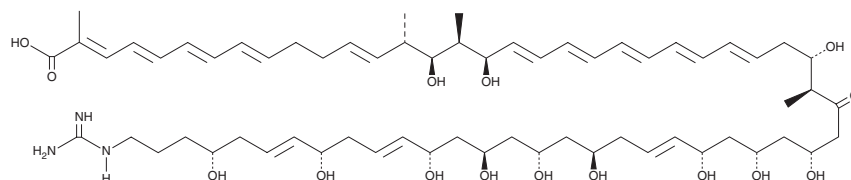
**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Bacterium/*Streptomyces* sp.

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



### Laboratory Procedures

Turonicin A is supplied as a solid. A stock solution may be made by dissolving the turonicin A in the solvent of choice, which should be purged with an inert gas. Turonicin A is soluble (≥10 mg/ml) in methanol and DMSO.

### Description

Turonicin A is a polyketide synthase-derived linear polyene originally isolated from *Streptomyces* that has antifungal and antibacterial activities.<sup>1</sup> It is active against *C. albicans*, *S. cerevisiae*, *B. subtilis*, and *S. aureus* (MICs = 2.7, 0.7, 85, and 340 nM, respectively).

### Reference

1. Chen, R., Minns, S.A., Kalaitzis, J.A., et al. Turonicin A, an antifungal linear polyene polyketide from an Australian *Streptomyces* sp. *J. Nat. Prod.* **86**(8), 2054-2058 (2023).

#### 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, 05/19/2026

#### 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