

# PRODUCT INFORMATION



## Kanchanamycin A

Item No. 44778

CAS Registry No.: 181637-98-9

Formal Name: propanedioic acid, mono[(10E,12E,18E,20E)-15-[(5E)-9-[(aminocarbonyl)amino]-1-methyl-5-nonenyl]-5,7,9,23,25,27,31,33,34,35-decahydroxy-10,14,22,26,30-pentamethyl-17-oxo-16,37-dioxabicyclo[31.3.1]heptatriaconta-10,12,18,20-tetraen-3-yl] ester

MF:  $C_{54}H_{90}N_2O_{18}$

FW: 1,055.3

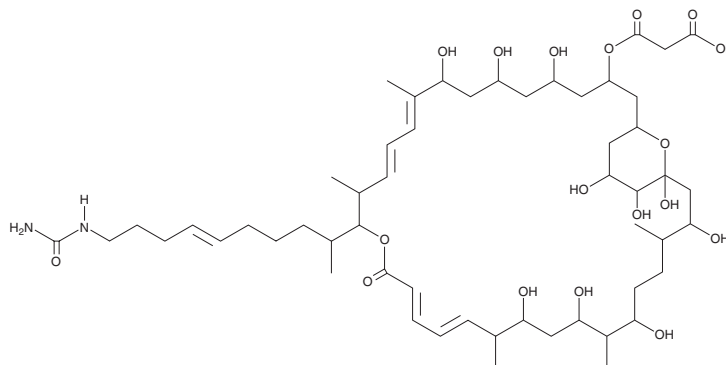
Purity:  $\geq 95\%$

Supplied as: A solid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 4$  years

Item Origin: Bacterium/*Streptomyces* sp.



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

### Laboratory Procedures

Kanchanamycin A is supplied as a solid. A stock solution may be made by dissolving the kanchanamycin A in the solvent of choice, which should be purged with an inert gas. Kanchanamycin A is soluble in methanol and DMSO.

### Description

Kanchanamycin A is a polyol macrolide antibiotic that has been found in *S. olivaceus*.<sup>1</sup> It is active against *S. aureus*, *P. fluorescens*, *A. viridinutans*, and *P. notatum* (MICs = 30, 3, 100, and 30  $\mu\text{g/ml}$ , respectively).

### Reference

1. Fiedler, H.P., Nega, M., Pfefferle, C., *et al.* Kanchanamycins, new polyol macrolide antibiotics produced by *Streptomyces olivaceus* Tü 4018. I. Taxonomy, fermentation, isolation and biological activities. *J. Antibiot. (Tokyo)* **49(8)**, 758-764 (1996).

#### 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, 11/25/2025

#### 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