

PRODUCT INFORMATION

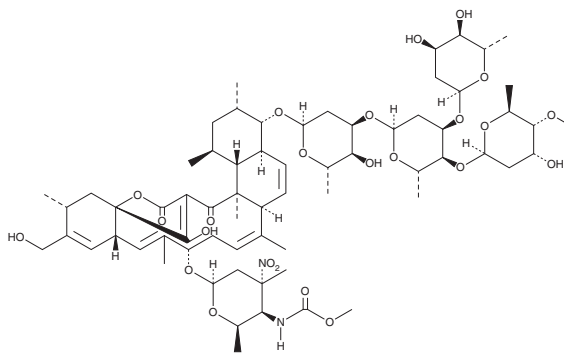


Kijanimitin

Item No. 20586

CAS Registry No.: 78798-08-0

Formal Name: 4-[[O-2,6-dideoxy- α -L-ribo-hexopyranosyl-(1 \rightarrow 3)-O-[2,6-dideoxy-4-O-methyl- β -L-ribo-hexopyranosyl-(1 \rightarrow 4)]-O-2,6-dideoxy- α -L-ribo-hexopyranosyl-(1 \rightarrow 3)-2,6-dideoxy- α -L-ribo-hexopyranosyl]oxy]-2,3S,4S,4



aS,6aS,9,10S,12aS,15R,16,20a,20bR-dodecahydro-21-hydroxy-14-(hydroxymethyl)-1S,3,7E,11E,15,20aS-hexamethyl-10-[[2,3,4,6-tetra-deoxy-4-[(methoxycarbonyl)amino]-3-C-methyl-3-nitro- β -D-xylo-hexopyranosyl]oxy]-18H-16aS,19-Metheno-16aH-benzo[b]naphth[2,1-j]oxacyclotetradecin-18,20(1H)-dione

Synonym: NSC 329515

MF: $C_{67}H_{100}N_2O_{24}$

FW: 1,317.5

Purity: $\geq 95\%$

Supplied as: A solid

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

Stability: ≥ 2 years

Item Origin: Bacteria/*Actinomadura* sp.

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

Laboratory Procedures

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

Description

Kijanimitin is an antibiotic first isolated from the fermentation broth of *A. kijaniata* SCC 1256.¹ It is active against a broad spectrum of microorganisms *in vitro* including *P. acnes* (MIC = 0.86 μ g/ml), *B. subtilis* (MIC < 0.13 μ g/ml), *Enterobacter* sp. (MIC = 64 μ g/ml), *Trichophyton* sp. (MIC = 17.5 μ g/ml), and *Microsporium* sp. (MIC = 17.5 μ g/ml).¹ At 250 mg/kg, kijanimitin was also shown to be effective in mice against *P. berghei*, a protozoan parasite that causes malaria in certain rodents.¹

Reference

1. Waitz, J. A., Horan, A. C., Kalyanpur, M., *et al.* Kijanimitin (Sch 25663), a novel antibiotic produced by *Actinomadura kijaniata* SCC 1256. Fermentation, isolation, characterization and biological properties. *J. Antibiot. (Tokyo)*. **34(9)**, 1101-1106 (1981).

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, 06/27/2022

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