

PRODUCT INFORMATION



Kigamicin C

Item No. 19445

CAS Registry No.: 680571-51-1

Formal Name: (1S,3R,4S,8aR,11aR)-3-[[[(2S,5S,6R)-5-[(2,6-dideoxy-3-O-methyl-b-D-arabinohexopyranosyl)oxy]tetrahydro-6-methyl-2H-pyran-2-yl]oxy]-1,2,3,4,8a,9,11,11a,13,14-decahydro-1,4,17,18-tetrahydroxy-11a-methyl-oxazolo[3,2-b]xantheno[4',3':2',4,5][1,3]benzodioxino[7,6-g]isoquinoline-16,19-dione

MF: C₄₁H₄₇NO₁₆

FW: 809.8

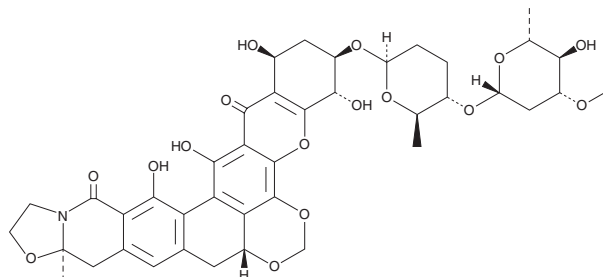
Purity: ≥85%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Amycolatopsis* sp.



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

Laboratory Procedures

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

Description

Kigamicin C is an antitumor antibiotic that has been found in *A. regifaucium*.^{1,2} It is active against various strains of *S. aureus*, including methicillin-resistant *S. aureus* (MRSA; MICs = 0.05-0.2 µg/ml), as well as *M. luteus* and *B. subtilis* (MICs = 0.05-0.2 µg/ml).¹ Kigamicin C is cytotoxic to PANC-1 pancreatic cancer cells in nutrient-deprived media at a 100-fold lower concentration than that required for cells cultured in nutrient-rich media.

References

1. Kunitomo, S., Lu, J., Esumi, H., *et al.* Kigamicins, novel antitumor antibiotics. I. Taxonomy, isolation, physico-chemical properties and biological activities. *J. Antibiot. (Tokyo)* **56(12)**, 1004-1011 (2003).
2. Tan, G.Y.A., Robinson, S., Lacey, E., *et al.* *Amycolatopsis regifaucium* sp. nov., a novel actinomycete that produces kigamicins. *Int. J. Syst. Evol. Microbiol.* **57(Pt 110)**, 2562-2567 (2007).

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, 01/04/2024

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