

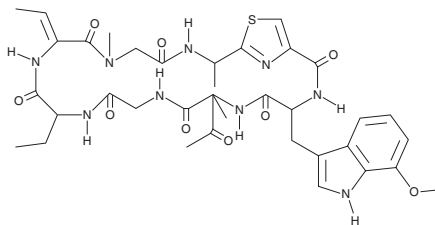
PRODUCT INFORMATION



Zelkovamycin

Item No. 27753

CAS Registry No.: 221197-33-7
MF: $C_{36}H_{45}N_9O_9S$
FW: 779.9
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 2 years
Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

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

Description

Zelkovamycin is a cyclic peptide antibiotic originally isolated from *Streptomyces*.¹ It inhibits growth of *X. oryzae*, *P. oryzae*, *S. aureus*, and *A. laidlawii* in a concentration-dependent manner when used at concentrations ranging from 0.01 to 300 $\mu g/ml$.

Reference

1. Zhang, H., Tomoda, H., Tabata, N., *et al.* Zelkovamycin, a new cyclic peptide antibiotic from *Streptomyces* sp. K96-0670. I. Production, isolation and biological properties. *J. Antibiot. (Tokyo)* **52(1)**, 29-33 (1999).

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, 06/13/2019

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