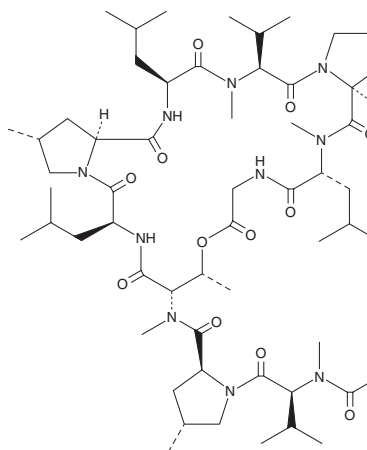


PRODUCT INFORMATION



Griselimycin Item No. 39306

CAS Registry No.: 26034-16-2
Formal Name: N-acetyl-N-methyl-L-valyl-(4R)-4-methyl-L-prolyl-N-methyl-L-threonyl-L-leucyl-(4R)-4-methyl-L-prolyl-L-leucyl-N-methyl-L-valyl-L-prolyl-N-methyl-D-leucyl-glycine, (10→3)-lactone
Synonym: NSC 246122
MF: C₅₇H₉₆N₁₀O₁₂
FW: 1,113.4
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces muensis*



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

Description

Griselimycin is a cyclic depsipeptide that has been found in *Streptomyces* and has antibacterial activity.^{1,2} It is active against *M. tuberculosis* strain H37Rv (MICs = 1-1.2 and 6.2 µg/ml in liquid culture and RAW 264.7 macrophages, respectively) but associated with resistance.² Griselimycin inhibits the protein-protein interaction between the *M. smegmatis* DNA polymerase sliding clamp, DnaN, and *M. smegmatis* DNA polymerase, leading to inhibition of DNA replication.

References

1. Terlain, B. and Thomas, J.P. Structure of griselimycin, a polypeptide antibiotic extracted from streptomyces cultures. *C.R. Acad. Sci. C Chim.* **269(24)**, 1546-1549 (1969).
2. Kling, A., Lukat, P., Almeida, D.V., *et al.* Targeting DnaN for tuberculosis therapy using novel griselimycins. *Science* **348(6239)**, 1106-1112 (2015).

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, 02/24/2025

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