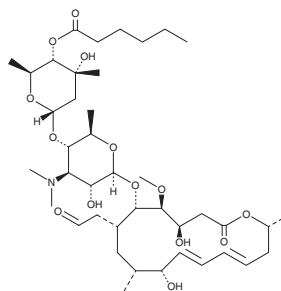


PRODUCT INFORMATION



Leucomycin A₁₃ Item No. 28121

CAS Registry No.: 78897-52-6
Formal Name: 4^B-hexanoate leucomycin V
Synonym: Kitasamycin A₁₃
MF: C₄₁H₆₉NO₁₄
FW: 800.0
Purity: ≥99%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces kitasatoensis*



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

Laboratory Procedures

Leucomycin A₁₃ is supplied as a solid. A stock solution may be made by dissolving the leucomycin A₁₃ in the solvent of choice. Leucomycin A₁₃ is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide, which should be purged with an inert gas.

Leucomycin A₁₃ is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Leucomycin A₁₃ is a macrolide antibiotic and a component of the leucomycin complex originally isolated from *S. kitasatoensis*.¹ It is active against *B. subtilis*, *S. aureus*, *M. luteus*, and *E. coli* with MIC values of 0.16, 0.16, 0.08 and >10 µg/ml, respectively.² It binds to ribosomes with an IC₅₀ value of 1.2 µM in a radioligand binding assay.

References

1. Hu, M. and Hu, C.-Q. Identification of the components of 16-membered macrolide antibiotics by LC/MS. *Anal. Chim. Acta* **535**(1-2), 89-99 (2005).
2. Tanaka, H., Moriuchi, I., Hirono, S., *et al.* Quantitative structure-activity relationships of O-acyl derivatives of leucomycin for antimicrobial and ribosome-binding activities. *Chem. Pharm. Bull. (Tokyo)* **33**(7), 2803-2808 (1985).

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, 10/11/2022

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