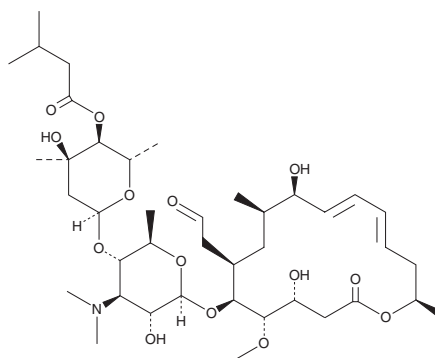


PRODUCT INFORMATION



Leucomycin A₁ Item No. 20585

CAS Registry No.: 16846-34-7
Formal Name: leucomycin V, 4^B-(3-methylbutanoate)
Synonyms: 3-deacetyl Josamycin, Kitasamycin A₁, Leucomycin V 4-isovalerate, 9-dihydro Niddamycin, Turimycin H5
MF: C₄₀H₆₇NO₁₄
FW: 786.0
Purity: ≥85%
UV/Vis.: λ_{max}: 233 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Leucomycin A₁ is supplied as a solid. A stock solution may be made by dissolving the leucomycin A₁ in the solvent of choice, which should be purged with an inert gas. Leucomycin A₁ is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Leucomycin A₁ is a major metabolite from the leucomycin complex, a family of macrolide antibiotics produced by *S. kitasatoensis* that are effective against Gram-positive bacteria, Gram-negative cocci, *Leptospira*, and mycoplasma.¹ Leucomycin A₁ is one of the more potent members of the complex.

Reference

1. Ishii, T., Hida, T., Iinuma, S., *et al.* TAN-1323 C and D, new concanamycin-group antibiotics; detection of the angiostatic activity with a wide range of macrolide antibiotics. *J. Antibiot. (Tokyo)* **48(1)**, 12-20 (1995).

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, 06/23/2021

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