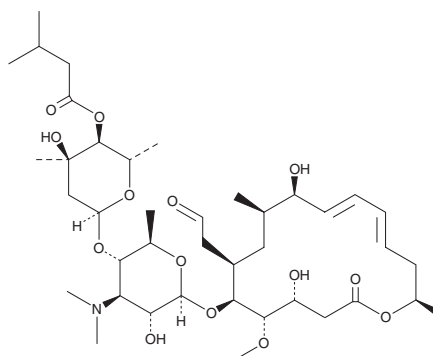


# PRODUCT INFORMATION



## Leucomycin A<sub>1</sub> Item No. 20585

**CAS Registry No.:** 16846-34-7  
**Formal Name:** leucomycin V, 4<sup>B</sup>-(3-methylbutanoate)  
**Synonyms:** 3-deacetyl Josamycin, Kitasamycin A<sub>1</sub>, Leucomycin V 4-isovalerate, 9-dihydro Niddamycin, Turimycin H5  
**MF:** C<sub>40</sub>H<sub>67</sub>NO<sub>14</sub>  
**FW:** 786.0  
**Purity:** ≥85%  
**UV/Vis.:** λ<sub>max</sub>: 233 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Leucomycin A<sub>1</sub> is supplied as a solid. A stock solution may be made by dissolving the leucomycin A<sub>1</sub> in the solvent of choice, which should be purged with an inert gas. Leucomycin A<sub>1</sub> is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

### Description

Leucomycin A<sub>1</sub> 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.<sup>1</sup> Leucomycin A<sub>1</sub> 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, 12/15/2022

#### 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