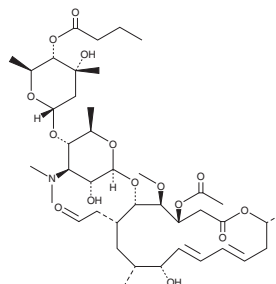


PRODUCT INFORMATION



Leucomycin A₄ Item No. 28066

CAS Registry No.: 18361-46-1
Formal Name: 3-acetate 4^B-butanoate leucomycin V
Synonym: Kitasamycin A₄
MF: C₄₁H₆₇NO₁₅
FW: 814.0
Purity: ≥85%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces kitatoensis*



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

Laboratory Procedures

Leucomycin A₄ is supplied as a solid. A stock solution may be made by dissolving the leucomycin A₄ in the solvent of choice, which should be purged with an inert gas. Leucomycin A₄ is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Leucomycin A₄ 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 A₄ is a macrolide antibiotic that has been found in *S. kitatoensis*.¹ It is active against a variety of bacteria, including *S. aureus*, *B. subtilis*, *C. diphtheriae*, *N. gonorrhoeae*, and *H. influenzae* (MICs = 0.15, 1.25, 0.15, 0.6, and 0.15 µg/ml, respectively).

Reference

1. Omura, S., Katagiri, M., Umezawa, I., et al. Structure-biological activities relationships among the leucomycins and their derivatives. *J. Antibiot. (Tokyo)* **21(9)**, 532-538 (1968).

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, 09/10/2024

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