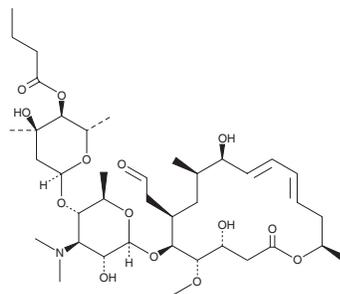


# PRODUCT INFORMATION



## Leucomycin A<sub>5</sub> Item No. 24177

**CAS Registry No.:** 18361-45-0  
**Formal Name:** 4<sup>B</sup>-butanoate leucomycin V  
**Synonyms:** Kitasamycin A<sub>5</sub>, Leukomycin A<sub>5</sub>,  
Turimycin H<sub>4</sub>  
**MF:** C<sub>39</sub>H<sub>65</sub>NO<sub>14</sub>  
**FW:** 771.9  
**Purity:** ≥85%  
**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>5</sub> is supplied as a solid. A stock solution may be made by dissolving the leucomycin A<sub>5</sub> in the solvent of choice, which should be purged with an inert gas. Leucomycin A<sub>5</sub> is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

### Description

Leucomycin A<sub>5</sub> is a metabolite from the leucomycin complex, which was originally isolated from *S. kitasatoensis*.<sup>1,2</sup> It is active against a variety of Gram-positive and Gram-negative bacteria (MICs = 0.04-0.8 µg/ml) but not against *K. pneumoniae*, *S. typhimurium*, or *E. coli* with MICs ranging from 5 to >10 µg/ml.<sup>1</sup> Leucomycin A<sub>5</sub> is also active against penicillin-susceptible and -resistant *S. pyogenes* (MICs = 0.8 and 3.2 µg/ml, respectively) and *S. faecalis* (MIC = 0.8 µg/ml).<sup>3</sup>

### References

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).
2. 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).
3. Singh, K. and Rakhit, S. Microbial transformation of leucomycin A<sub>5</sub>. *J. Antibiot. (Tokyo)* **32(1)**, 78-80 (1979).

#### 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, 11/14/2022

#### 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