

# PRODUCT INFORMATION



## Urdamycin A

Item No. 29510

CAS Registry No.: 98474-21-6

Formal Name: (3R,4aR,12bS)-9-[2,6-dideoxy-3-O-[(2S,5S,6S)-5-[(2,6-dideoxy-β-D-arabino-hexopyranosyl)oxy]tetrahydro-6-methyl-2H-pyran-2-yl]-β-D-arabino-hexopyranosyl]-3,4,4a,12b-tetrahydro-3,4a,8-trihydroxy-3-methyl-12b-[[[(2S,5S,6S)-tetrahydro-5-hydroxy-6-methyl-2H-pyran-2-yl]oxy]-benz[a]anthracene-1,7,12(2H)-trione

Synonyms: Antibiotic A 32Y2, Kerriamycin B, NSC 372208, NSC 606397

MF: C<sub>43</sub>H<sub>56</sub>O<sub>17</sub>

FW: 844.9

Purity: ≥70%

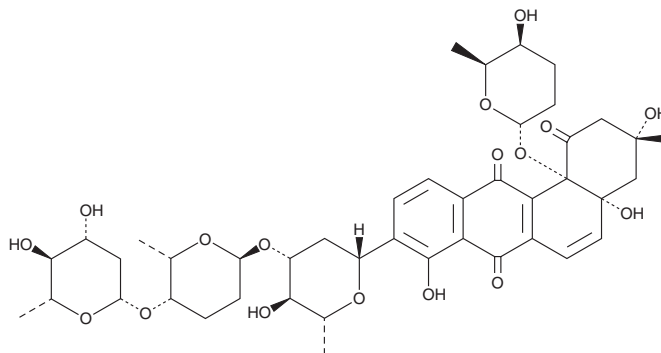
Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Streptomyces* sp.

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



### Laboratory Procedures

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

### Description

Urdamycin A is a bacterial metabolite originally isolated from *S. fradiae* that has antibacterial and anticancer activities.<sup>1,2</sup> It is active against a variety of Gram-positive and Gram-negative bacteria, including *B. subtilis* and strains of *Arthrobacter* and *Streptomyces*, but not the fungus *S. cerevisiae*, in a disc assay when used at a concentration of 1 mg/ml.<sup>2</sup> Urdamycin A is cytotoxic to L1210 and HT-29, but not A549, cancer cells (IC<sub>50</sub>s = 7.5, 5, and >10 µg/ml, respectively).<sup>1</sup>

### References

1. Henkel, T., Ciesiolka, T., Rohr, J., et al. Urdamycins, new angucycline antibiotics from *Streptomyces fradiae*. V. Derivatives of urdamycin A. *J. Antibiot. (Tokyo)* **42**(2), 299-311 (1989).
2. Drautz, H., Zähler, H., Rohr, J., et al. Metabolic products of microorganisms. 234. Urdamycins, new angucycline antibiotics from *Streptomyces fradiae*. I. Isolation, characterization and biological properties. *J. Antibiot. (Tokyo)* **39**(12), 1657-1669 (1986).

#### 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, 11/10/2022

#### 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