

# PRODUCT INFORMATION



## Asukamycin

Item No. 20133

**CAS Registry No.:** 61116-33-4  
**Formal Name:** (2E,4E,6E)-7-cyclohexyl-N-[(1S,5S,6R)-5-hydroxy-5-[(1E,3E,5E)-7-[(2-hydroxy-5-oxo-1-cyclopenten-1-yl)amino]-7-oxo-1,3,5-heptatrien-1-yl]-2-oxo-7-oxabicyclo[4.1.0]hept-3-en-3-yl]-2,4,6-heptatrienamide

**Synonyms:** AM1042, Asukamycin A

**MF:** C<sub>31</sub>H<sub>34</sub>N<sub>2</sub>O<sub>7</sub>

**FW:** 546.6

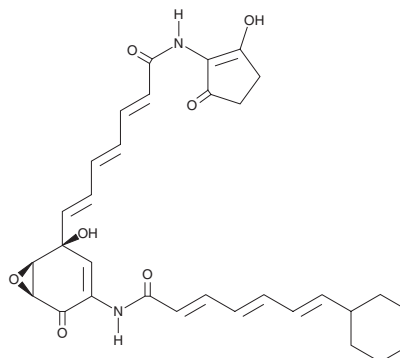
**Purity:** ≥95%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** *Streptomyces* sp.



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

### Laboratory Procedures

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

### Description

Asukamycin is polyketide isolated from the *S. nodosus* subspecies *asukaensis* that demonstrates a broad range of antibiotic functions. It has been shown to inhibit growth of various tumor cell lines (IC<sub>50</sub>S = 1-5 μM) by activating caspases 8 and 3.<sup>1</sup>

### Reference

1. Shipley, P. R., Donnelly, C. C. A., Le, C. H., *et al.* Antitumor activity of asukamycin, a secondary metabolite from the actinomycete bacterium *Streptomyces nodosus* subspecies *asukaensis*. *Int. J. Mol. Med.* **24**(5) 711-715 (2009).

#### 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, 11/03/2022

#### 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