

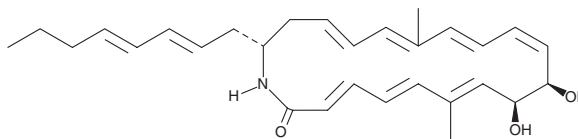
PRODUCT INFORMATION



Heronamide C

Item No. 23644

CAS Registry No.: 1257083-94-5
Formal Name: 9S,10R-dihydroxy-7,15-dimethyl-20R-(2E,4E)-2,4-octadien-1-yl-azacycloeicosa-3E,5E,7E,11Z,13E,15E,17E-heptaen-2-one
MF: $C_{29}H_{39}NO_3$
FW: 449.6
Purity: $\geq 90\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
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

Heronamide C is supplied as a crystalline solid. A stock solution may be made by dissolving the heronamide C in the solvent of choice, which should be purged with an inert gas. Heronamide C is soluble in the organic solvent DMSO.

Description

Heronamide C is a polyketide macrolactam antifungal agent produced by marine *Actinomycetes*.¹ It inhibits the growth of and induces abnormal accumulation of cell wall material in wild-type fission yeast cells (MIC = 0.13-0.28 μM). Heronamide C lacks activity against Gram-positive bacteria and human HeLa and MDA-MB-231 cancer cell lines.² However, it induces reversible morphological changes, including formation of large intracellular structures, in HeLa cells at a concentration of 20 μM .

References

1. Fujita, K., Sugiyama, R., Nishimura, S., *et al.* Stereochemical assignment and biological evaluation of BE-14106 unveils the importance of one acetate unit for the antifungal activity of polyene macrolactams. *J. Nat. Prod.* **79**(7), 1877-1880 (2016).
2. Paju, R., Piggott, A.M., Conte, M.M., *et al.* Heronamides A-C, new polyketide macrolactams from an Australian marine-derived *Streptomyces* sp. A biosynthetic case for synchronized tandem electrocyclization. *Org. Biomol. Chem.* **8**(20), 4682-4689 (2010).

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, 12/22/2022

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