# **PRODUCT** INFORMATION



## Padanamide A

Item No. 25085

CAS Registry No.:	1314881-80-5	NH <sub>2</sub>	$\frown$
Formal Name:	(S)-N-((2S,3S,4S)-5-(((S)-1-carbamoyl-2- oxopyrrolidin-3-yl)amino)-3-hydroxy-4-methyl-5- oxo-1-phenylpentan-2-yl)-2-((2R,3R)-3-hydroxy- 2-(2-methoxyacetamido)-4-methylpentanoyl)		
	hexahydropyridazine-3-carboxamide	✓ •N <sup>2</sup>	$\downarrow$ $\uparrow$ $H$ $\downarrow$ $\downarrow$ $\downarrow$
MF:	$C_{31}H_{47}N_7O_9$	н	
FW:	661.8		$\overline{N}$
Purity:	≥95%	Ì	
Supplied as:	A solid		0
Storage:	-20°C		
Stability:	≥4 years		
Item Origin:	Bacterium/Streptomyces sp.		
Information represents the product exectlications. Batch exectlic analytical results are provided on each cartificate of analysis			

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

#### Laboratory Procedures

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

#### Description

Padanamide A is a bacterial metabolite originally isolated from a marine Streptomyces species.<sup>1</sup> It is active against five strains of P. falciparum (EC<sub>50</sub>s = 160-340 nM), is not cytotoxic to HEK293 or HepG2 human cell lines, but is cytotoxic to Jurkat T lymphocytes with an IC<sub>50</sub> value of approximately 60  $\mu$ g/ml.<sup>2,3</sup>

#### References

- 1. Nam, S.J., Kauffman, C.A., Jensen, P.R., et al. Isolation and characterization of actinoramides A-C, highly modified peptides from a marine Streptomyces sp. Tetrahedron 67(35), 6707-6712 (2011).
- 2. Cheng, K.-C., Cao, S., Raveh, A., et al. Actinoramide A identified as a potent antimalarial from titration-based screening of marine natural product extracts. J. Nat. Prod. 78(10), 2411-2422 (2015).
- 3. Williams, D.E., Dalisay, D.S., Patrick, B.O., et al. Padanamides A and B, highly modified linear tetrapeptides produced in culture by a Streptomyces sp. isolated from a marine sediment. Org. Lett. 13(15), 3936-3939 (2011).

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

uver 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, 11/11/2022

### 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