

# PRODUCT INFORMATION



## Unguisin A

Item No. 23878

CAS Registry No.: 226956-06-5

Formal Name: cyclo(D-alanyl-D-tryptophyl-4-aminobutanoyl-D-alanyl-D-valyl-L-phenylalanyl-D-valyl)

MF:  $C_{40}H_{54}N_8O_7$

FW: 758.9

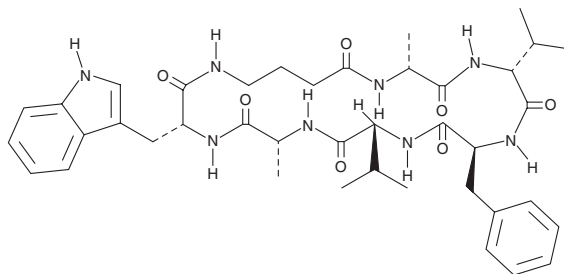
Purity:  $\geq 95\%$

Supplied as: A solid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 4$  years

Item Origin: Fungus/*Aspergillus unguis*



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

### Laboratory Procedures

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

### Description

Unguisin A is a cyclic heptapeptide originally isolated from the marine fungus *E. unguis* whose structure is comprised of amino acids and GABA.<sup>1</sup> It binds to phosphate, pyrophosphate, and chloride but has no effect on chloride transport in a liposome-based assay.<sup>2</sup> It exhibited moderate antibacterial activity against *S. aureus* and *V. parahaemolyticus* in one assay but no activity against a variety of bacteria, including *S. aureus*, in another.<sup>1,2</sup>

### References

1. Malmstrom, J. Unguisins A and B: New cyclic peptides from the marine-derived fungus *Emericella unguis* *J. Nat. Prod.* **62**(5), 787-789 (1999).
2. Ariawan, A.D., Webb, J.E.A., Howe, E.N.W., *et al.* Cyclic peptide unguisin A is an anion receptor with high affinity for phosphate and pyrophosphate. *Org. Biomol. Chem.* **15**(14), 2962-2967 (2017).

#### 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, 08/23/2023

#### 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