

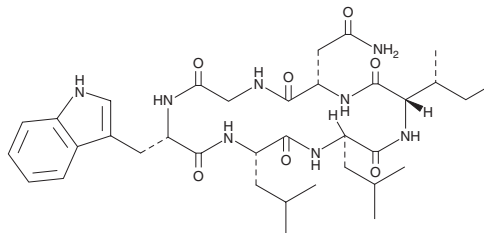
# PRODUCT INFORMATION



## Desotamide

Item No. 28129

**CAS Registry No.:** 194660-14-5  
**Formal Name:** cyclo(L-alloisoleucyl-L-asparaginylglycyl-L-tryptophyl-L-leucyl-D-leucyl)  
**Synonym:** Desotamide A  
**MF:** C<sub>35</sub>H<sub>52</sub>N<sub>8</sub>O<sub>7</sub>  
**FW:** 696.8  
**Purity:** ≥95%  
**Supplied as:** A lyophilisate  
**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

Desotamide is supplied as a lyophilisate. A stock solution may be made by dissolving the desotamide in the solvent of choice, which should be purged with an inert gas. Desotamide is soluble in ethanol, methanol, dimethyl formamide, and DMSO.

### Description

Desotamide is a cyclic hexapeptide antibiotic originally isolated from *Streptomyces*.<sup>1</sup> It is active against *S. aureus*, *S. pneumoniae*, and methicillin-resistant *S. epidermidis* (MRSE; MICs = 16, 12.5, and 32 µg/ml, respectively).<sup>2</sup>

### References

1. Miao, S., Anstee, M.R., LaMarco, K., *et al.* Inhibition of bacterial RNA polymerases. Peptide metabolites from the cultures of *Streptomyces* sp. *J. Nat. Prod.* **60(8)**, 858-861 (1997).
2. Song, Y., Li, Q., Liu, X., *et al.* Cyclic hexapeptides from the deep South China Sea-derived *Streptomyces scopuliridis* SCSIO ZJ46 active against pathogenic Gram-positive bacteria. *J. Nat. Prod.* **77(8)**, 1937 (2014).

#### 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, 10/11/2022

#### 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