

PRODUCT INFORMATION

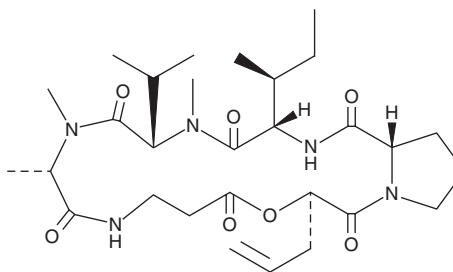


Destruxin A

Item No. 20428

CAS Registry No.: 6686-70-0
Formal Name: cyclo[N-methyl-L-alanyl-β-alanyl-(2R)-2-hydroxy-4-pentenoyl-L-prolyl-L-isoleucyl-N-methyl-L-valyl]

Synonym: SB 242533
MF: $C_{29}H_{47}N_5O_7$
FW: 577.7
Purity: ≥96%
Supplied as: A powder
Storage: -20°C
Stability: ≥4 years



Special Conditions: Protect from moisture; avoid direct light

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

Laboratory Procedures

Destruxin A is supplied as a powder. A stock solution may be made by dissolving the destruxin A in the solvent of choice. Destruxin A is soluble in organic solvents such as acetonitrile, DMSO, and dimethyl formamide, which should be purged with an inert gas.

Description

Destruxin A is a cyclic hexadepsipeptide mycotoxin first isolated from *O. destructor* and commonly found in entomopathogenic fungi.¹ It has insecticidal activity by affecting diverse pathways, including those related to immune response, calcium signaling, and development.^{2,3}

References

1. Pedras, M., Zaharia, L. I. and Ward, D. E. The destruxins: Synthesis, biosynthesis, biotransformation, and biological activity. *Phytochemistry* **59**(6), 579-596 (2002).
2. Han, P., Jin, F., Dong, X., et al. Transcript and protein profiling analysis of the Destruxin A-induced response in larvae of *Plutella xylostella*. *PLoS One* **8**(4), (2013).
3. Pal, S., St.Leger, R. J. and Wu, L. P. Fungal peptide Destruxin A plays a specific role in suppressing the innate immune response in *Drosophila melanogaster*. *J. Biol. Chem.* **282**(12), 8969-8977 (2007).

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, 02/25/2025

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