

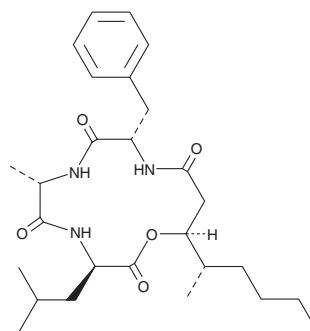
# PRODUCT INFORMATION



## Beauveriolide I

Item No. 31774

**CAS Registry No.:** 154491-55-1  
**Formal Name:** cyclo[L-alanyl-D-leucyl-(3S,4S)-3-hydroxy-4-methyloctanoyl-L-phenylalanyl]  
**MF:** C<sub>27</sub>H<sub>41</sub>N<sub>3</sub>O<sub>5</sub>  
**FW:** 487.6  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Fungus/*Beauveria bassiana*



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

### Laboratory Procedures

Beauveriolide I is supplied as a solid. A stock solution may be made by dissolving the beauveriolide I in the solvent of choice, which should be purged with an inert gas. Beauveriolide I is soluble in DMSO.

### Description

Beauveriolide I is a cyclodepsipeptide that has been found in *Beauveria* and an inhibitor of lipid droplet formation.<sup>1</sup> It inhibits lipid droplet formation when used at concentrations of 3 and 10 μM, as well as inhibits cholesterol synthesis (IC<sub>50</sub> = 0.78 μM), in primary mouse peritoneal macrophages.<sup>1,2</sup> Beauveriolide I also inhibits acyl-coenzyme A:cholesterol acyltransferase (ACAT) activity in mouse macrophage membranes (IC<sub>50</sub> = 6 μM).<sup>2</sup>

### References

1. Namatame, I., Tomoda, H., Si, S., et al. Beauveriolides, specific inhibitors of lipid droplet formation in mouse macrophages, produced by *Beauveria* sp. FO-6979. *J. Antibiot. (Tokyo)* **52(1)**, 1-6 (1999).
2. Namatame, I., Tomoda, H., Ishibashi, S., et al. Antiatherogenic activity of fungal beauveriolides, inhibitors of lipid droplet accumulation in macrophages. *Proc. Nat. Acad. Sci. USA* **101(3)**, 737-742 (2004).

#### 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, 08/17/2023

#### 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