

# PRODUCT INFORMATION

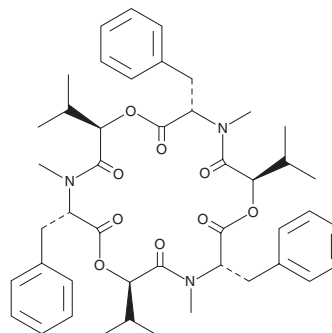


## Beauvericin

Item No. 11426

**CAS Registry No.:** 26048-05-5  
**Formal Name:** cyclo[(2R)-2-hydroxy-3-methylbutanoyl-N-methyl-L-phenylalanyl-(2R)-2-hydroxy-3-methylbutanoyl-N-methyl-L-phenylalanyl-(2R)-2-hydroxy-3-methylbutanoyl-N-methyl-L-phenylalanyl]

**MF:** C<sub>45</sub>H<sub>57</sub>N<sub>3</sub>O<sub>9</sub>  
**FW:** 784.0  
**Purity:** ≥95%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Beauvericin is supplied as a crystalline solid. A stock solution may be made by dissolving the beauvericin in the solvent of choice, which should be purged with an inert gas. Beauvericin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of beauvericin in these solvents is approximately 30 mg/ml.

Beauvericin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, beauvericin should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Beauvericin has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Beauvericin is a mycotoxin originally derived from *B. bassiana*, a common fungal parasite of arthropods. It is also produced by the fungus *Fusarium*, a pathogen of insects and plants.<sup>1</sup> Beauvericin is cytotoxic against insect cells, killing the insect cell line SF-9 with a 50% cytotoxic concentration of 2.5 μM.<sup>2</sup> In mammalian cells, beauvericin induces apoptosis with an IC<sub>50</sub> value of 4.5 μM.<sup>1,3</sup>

### References

1. Logrieco, A., Moretti, A., Castella, G., *et al.* Beauvericin production by *Fusarium* species. *Appl. Environ. Microbiol.* **64**(8), 3084-3088 (1998).
2. Calo, L., Fornelli, F., Nenna, S., *et al.* Beauvericin cytotoxicity to the invertebrate cell line SF-9. *J. Appl. Genet.* **44**(4), 515-520 (2003).
3. Tonshin, A.A., Teplova, V.V., Andersson, M.A., *et al.* The *Fusarium* mycotoxins enniatins and beauvericin cause mitochondrial dysfunction by affecting the mitochondrial volume regulation, oxidative phosphorylation and ion homeostasis. *Toxicology* **276**, 49-57 (2010).

#### 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, 10/06/2022

#### 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