

# PRODUCT INFORMATION



## Phepropeptin D

Item No. 10008111

**Formal Name:** (3S,6R,9S,12S,15R,20aS)-3,15-dibenzyl-9-((S)-sec-butyl)-6,12-diisobutyltetradecahydropyrrolo [1,2-a][1,4,7,10,13,16]hexaazacyclooctadecine-1,4,7,10,13,16-hexaone

**Synonym:** Cyclic(Leu-D-Phe-D-Pro-Phe-D-Leu- $\xi$ -Ile)

**Peptide Sequence:** C(LfpFIX) (X =  $\xi$ -Isoleucine)

**MF:** C<sub>41</sub>H<sub>58</sub>N<sub>6</sub>O<sub>6</sub>

**FW:** 730.9

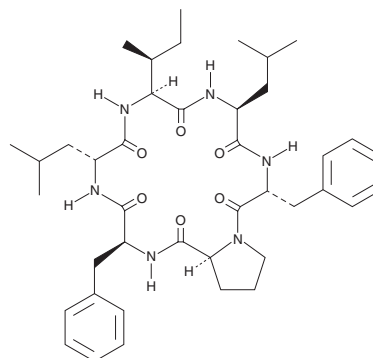
**Purity:**  $\geq 98\%$

**Supplied as:** A solid

**Storage:** -20°C

**Stability:**  $\geq 4$  years

**Item Origin:** Bacterium/*Streptomyces* sp.



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

### Laboratory Procedures

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

### Description

Phepropeptin D is a cyclic peptide originally isolated from *Streptomyces* that has proteasome inhibitory activity.<sup>1</sup> It inhibits isolated mouse liver proteasomes (IC<sub>50</sub> = 7.8  $\mu$ g/ml).

### Reference

1. Sekizawa, R., Momose, I., Kinoshita, N., *et al.* Isolation and structural determination of phepropeptins A, B, C, and D, new proteasome inhibitors, produced by *Streptomyces* sp. *J. Antibiot. (Tokyo)* **54(11)**, 874-881 (2001).

#### 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, 07/02/2025

#### 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](http://WWW.CAYMANCHEM.COM)