

PRODUCT INFORMATION



Onychocin B

Item No. 33134

Formal Name: 6S,12R-dibenzyl-3S-((S)-sec-butyl)-9S-isopropyl-1,7-dimethyl-1,4,7,10-tetraazacyclododecane-2,5,8,11-tetraone

MF: C₃₁H₄₂N₄O₄

FW: 534.7

Purity: ≥95%

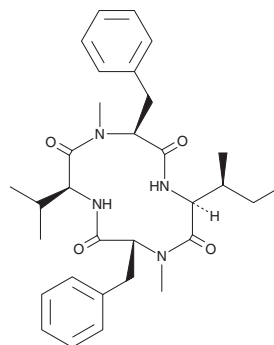
Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungi/*Aspergillus* sp.

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



Laboratory Procedures

Onychocin B is supplied as a solid. A stock solution may be made by dissolving the onychocin B in the solvent of choice, which should be purged with an inert gas. Onychocin B is soluble in methanol and DMSO.

Description

Onychocin B is a cyclic tetrapeptide originally isolated from *Onychocola sclerotica* that has calcium channel inhibitory activity.¹ It inhibits the voltage-gated calcium (Ca_v) channel Ca_v1.2 with an IC₅₀ value of 7.1 μM.

Reference

1. Pérez-Victoria, I., Martín, J., González-Menéndez, V., *et al.* Isolation and structural elucidation of cyclic tetrapeptides from *Onychocola sclerotica*. *J. Nat. Prod.* **75**(6), 1210-1214 (2012).

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, 11/14/2022

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