

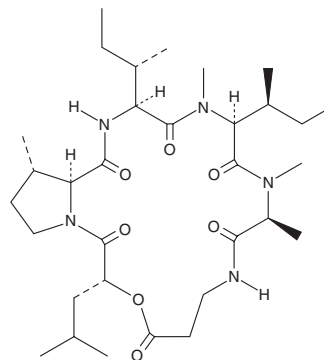
PRODUCT INFORMATION



Homodestcardin

Item No. 30052

CAS Registry No.: 917382-84-4
Formal Name: cyclo[N-methyl-L-alanyl-β-alanyl-(2R)-2-hydroxy-4-methylpentanoyl-(3S)-3-methyl-L-prolyl-L-isoleucyl-N-methyl-L-isoleucyl]
MF: C₃₂H₅₅N₅O₇
FW: 621.8
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Trichothecium* sp.



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

Laboratory Procedures

Homodestcardin is supplied as a solid. A stock solution may be made by dissolving the homodestcardin in the solvent of choice, which should be purged with an inert gas. Homodestcardin is soluble in organic solvents such as ethanol, methanol, DMSO, and dichloromethane.

Description

Homodestcardin is a cyclic depsipeptide fungal metabolite that has been found in *T. roseum* and has immunosuppressant activity.¹ It inhibits concanavalin A-induced proliferation of isolated mouse T cells (IC₅₀ = 0.86 μM).

Reference

1. Zhang, A.H., Wang, X.Q., Han, W.B., *et al.* Discovery of a new class of immunosuppressants from *Trichothecium roseum* co-inspired by cross-kingdom similarity in innate immunity and pharmacophore motif. *Chem. Asian J.* **8(12)**, 3101-3107 (2013).

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, 08/10/2023

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