

PRODUCT INFORMATION

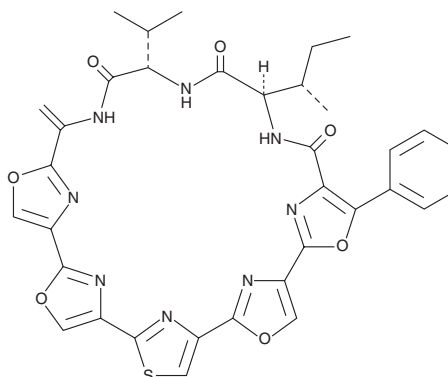


Mechercharmycin A

Item No. 40868

CAS Registry No.: 822520-96-7
Formal Name: cyclo[D-alloisoleucyl-L-valyl-2'-[2'-(1-aminoethenyl)[2,4'-bioxazol]-4-yl]-4-thiazolyl]-5-phenyl[2,4'-bioxazole]-4-carbonyl]

Synonym: IB-01211
MF: C₃₅H₃₂N₈O₇S
FW: 708.7
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Microbial



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

Laboratory Procedures

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

Description

Mechercharmycin A is a cyclic peptide originally isolated from *Thermoactinomyces* that has anticancer activity.¹ It induces cytotoxicity in A549 lung cancer and Jurkat leukemia cells (IC₅₀s = 40 and 46 nM, respectively).

Reference

1. Kanoh, K., Matsuo, Y., Adachi, K., *et al.* Mechercharmycins A and B, cytotoxic substances from marine-derived *Thermoactinomyces* sp. YM3-251. *J. Antibiot. (Tokyo)* **58(4)**, 289-292 (2005).

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, 04/24/2024

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