

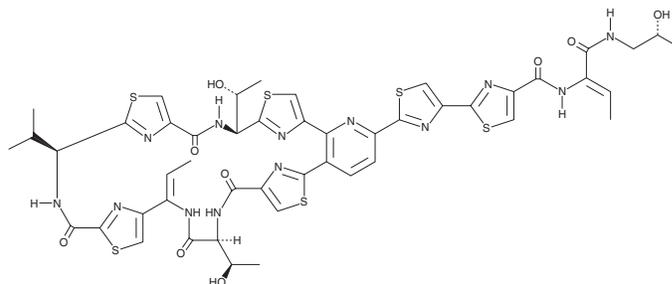
PRODUCT INFORMATION



Micrococcin P1

Item No. 17093

CAS Registry No.: 67401-56-3
Formal Name: 2'-[(11S,14Z,21S,28S)-14-ethylidene-9,10,11,12,13,14,20,21,27,28-decahydro-28-[(1R)-1-hydroxyethyl]-11-[(1R)-1-hydroxyethyl]-21-(1-methylethyl)-9,12,19,26-tetraoxo-19H,26H-8,5:18,15:25,22:32,29-tetranitrilo-5H,15H-pyrido[3,2-m][1,11,17,24,4,7,20,27]tetrathiatetraazacyclotriacontin-2-yl]-N-[(1Z)-1-[[[(2R)-2-hydroxypropyl]amino]carbonyl]-1-propen-1-yl]-[2,4'-bithiazole]-4-carboxamide
MF: C₄₈H₄₉N₁₃O₉S₆
FW: 1,144.4
Purity: ≥95%
Supplied as: A 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

Micrococcin P1 is supplied as a solid. A stock solution may be made by dissolving the micrococcin P1 in the solvent of choice. Micrococcin P1 is soluble in organic solvents such as ethanol, DMSO, dimethyl formamide, and methanol which should be purged with an inert gas.

Micrococcin P1 is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Micrococcin P1 is a macrocyclic peptide antibiotic that functions as an acceptor-site-specific inhibitor of ribosomal protein synthesis, effectively preventing the growth of Gram-positive bacteria without affecting that of Gram-negative bacteria.^{1,2}

References

1. Carnio, M.C., Hötzel, A., Rudolf, M., *et al.* The macrocyclic peptide antibiotic micrococcin P(1) is secreted by the food-borne bacterium *Staphylococcus equorum* WS 2733 and inhibits *Listeria monocytogenes* on soft cheese. *Appl. Environ. Microbiol.* **66(6)**, 2378-2384 (2000).
2. Mikolajka, A., Liu, H., Chen, Y., *et al.* Differential effects of thiopeptide and orthosomycin antibiotics on translational GTPases. *Chem. Biol.* **18(5)**, 589-600 (2011).

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, 12/14/2022

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