

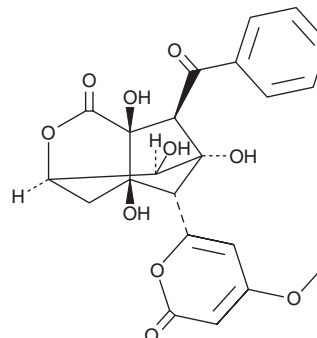
PRODUCT INFORMATION



Enterocin

Item No. 20592

CAS Registry No.: 59678-46-5
Formal Name: (3R,4aR,5S,6S,7S,7aS,8R)-7-benzoylhexahydro-4a,6,7a,8-tetrahydroxy-5-(4-methoxy-2-oxo-2H-pyran-6-yl)-3,6-methanocyclopenta[c]pyran-1(3H)-one
Synonym: Vulgamycin
MF: C₂₂H₂₀O₁₀
FW: 444.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥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

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

Enterocin 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

Enterocin is an antibiotic originally isolated from *Streptomyces* sp. It has static activities against a broad range of Gram-positive and Gram-negative bacteria but has no activity against fungi, including yeasts.¹ This natural product was named for its strong activity against *Enterobacteriaceae* and is not related to the large group of bacterial proteins isolated from *Enterococci* that are classified as bacteriocins.

Reference

1. Miyairi, N., Saki, H.-I., Konomi, T., *et al.* Enterocin, a new antibiotic taxonomy, isolation and characterization. *J. Antibiot. (Tokyo)* **29**(3), 227-235 (1976).

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, 11/11/2022

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