

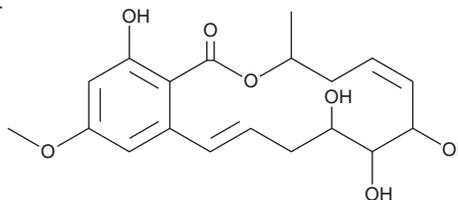
PRODUCT INFORMATION



LL-Z 1640-4

Item No. 19159

CAS Registry No.: 66018-41-5
Formal Name: 3,4,7,8,9,10-hexahydro-7,8,9,16-tetrahydroxy-14-methoxy-3-methyl-1H-2-benzoxacyclotetradecin-1-one
Synonyms: Antibiotic LL-Z 1640-4, (5Z)-Zeaenol
MF: C₁₉H₂₄O₇
FW: 364.4
Purity: ≥99%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Curvularia* sp.



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

Laboratory Procedures

LL-Z 1640-4 is supplied as a solid. A stock solution may be made by dissolving the LL-Z 1640-4 in the solvent of choice, which should be purged with an inert gas. LL-Z 1640-4 is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

LL-Z 1640-4 is a *cis*-enol resorcylic acid lactone that has been shown to exhibit both antiviral and antiprotozoan activity.¹ While LL-Z 1640-4 is inactive against JNK/p38 kinase signaling, its companion antibiotic, (5Z)-7-oxo zeaenol (Item No. 17459), has been identified as an irreversible inhibitor of the MAP3K TAK1, blocking T cell activation.² Thus, LL-Z 1640-4 is useful as a negative control to help dissect the selectivity of this MAP3K inhibitor.

References

1. Ellestad, G.A., Lovell, F.M., Perkinson, N.A., *et al.* New zearalenone related macrolides and isocoumarins from an unidentified fungus. *J. Org. Chem.* **43(12)**, 2339-2343 (1978).
2. Sellès, P. and Lett, R. Convergent stereospecific synthesis of LL-Z1640-2 (or C292), hypothemycin and related macrolides. Part 2. *Tetrahedron Lett.* **43(26)**, 4627-4631 (2002).

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, 08/09/2023

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