

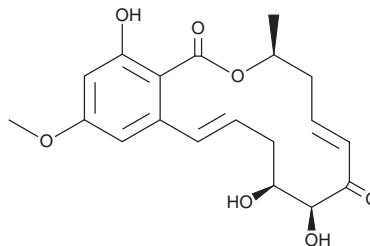
PRODUCT INFORMATION



(5E)-7-Oxozeaenol

Item No. 29840

CAS Registry No.: 1198574-97-8
Formal Name: (3S,5E,8S,9S,11E)-3,4,9,10-tetrahydro-8,9,16-trihydroxy-14-methoxy-3-methyl-1H-2-benzoxacyclotetradecin-1,7(8H)-dione
Synonym: LL-Z 1640-2
MF: C₁₉H₂₂O₇
FW: 362.4
Purity: ≥95%
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

(5E)-7-Oxozeaenol is supplied as a solid. A stock solution may be made by dissolving the (5E)-7-oxozeaenol in the solvent of choice, which should be purged with an inert gas. (5E)-7-Oxozeaenol is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Description

(5E)-7-Oxozeaenol is a resorcylic acid lactone that has been found in the fungus MSX 63935 and has enzyme inhibitory and anticancer activities.^{1,2} It inhibits TGF- β -activated kinase 1 (TAK-1; IC₅₀ = 1.3 μ M).¹ (5E)-7-Oxozeaenol inhibits proliferation of MCF-7, H460, SF-268, HT-29, and MDA-MB-435 human cancer cells with IC₅₀ values of 4.9, 1.2, 5.6, 4.4, and 5.5 μ M, respectively.²

References

1. Fakhouri, L., El-Elimat, T., Hurst, D.P., *et al.* Isolation, semisynthesis, covalent docking and transforming growth factor beta-activated kinase 1 (TAK1)-inhibitory activities of (5Z)-7-oxozeaenol analogues. *Bioorg. Med. Chem.* **23(21)**, 6993-6999 (2015).
2. Ayers, S., Graf, T.N., Adcock, A.F., *et al.* Resorcylic acid lactones with cytotoxic and NF- κ B inhibitory activities and their structure-activity relationships. *J. Nat. Prod.* **74(5)**, 1126-1131 (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, 10/06/2022

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