

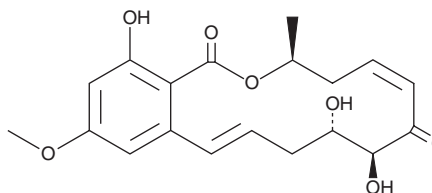
PRODUCT INFORMATION



(5Z)-7-Oxozeaenol

Item No. 17459

CAS Registry No.: 253863-19-3
Formal Name: (3S,5Z,8S,9S,11E)-3,4,9,10-tetrahydro-8,9,16-trihydroxy-14-methoxy-3-methyl-1H-2-benzoxacyclotetradecin-1,7(8H)-dione
Synonyms: FR148083, L-783,279, LL-Z 1640-2
MF: $C_{19}H_{22}O_7$
FW: 362.4
Purity: $\geq 99\%$
UV/Vis.: λ_{max} : 232, 273, 317 nm
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

(5Z)-7-Oxozeaenol is supplied as a solid. A stock solution may be made by dissolving the (5Z)-7-oxozeaenol in the solvent of choice, which should be purged with an inert gas. (5Z)-7-Oxozeaenol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (5Z)-7-oxozeaenol in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

(5Z)-7-Oxozeaenol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (5Z)-7-oxozeaenol should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. (5Z)-7-Oxozeaenol has a solubility of approximately 0.05 mg/ml in a 1:20 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(5Z)-7-Oxozeaenol is a resorcylic acid lactone fungal metabolite and an inhibitor of TGF- β activated kinase 1 (TAK1; IC_{50} = 8.1 nM in the presence of TAK1-binding protein 1 (TAB1)).¹ It is selective for TAK1 over MEK1 and MEKK1 (IC_{50} s = 411 and 268 nM, respectively), as well as apoptosis signal-regulating kinase 1 (ASK1) and I κ B kinase β (IKK β) at 300 nM. (5Z)-7-Oxozeaenol inhibits NF- κ B activity in a reporter assay using HEK293 cells expressing TAK1 and TAB1 (IC_{50} = 83 nM) and reduces IL-1-induced increases in COX-2 levels in mouse embryonic fibroblasts (MEFs). It sensitizes HeLa cells to apoptosis induced by TNF-related apoptosis-inducing ligand (TRAIL) when used at a concentration of 300 nM and induces necroptosis when used in combination with TNF and zVAD in primary mouse bone marrow-derived macrophages (BMDMs) at 100 nM.^{2,3} Topical application of (5Z)-7-oxozeaenol (10 μ l of a 1 mg/ml solution) prevents picryl chloride-induced increases in ear thickness in mice.¹

References

1. Ninomiya-Tsuji, J., Kajino, T., Ono, K., *et al.* A resorcylic acid lactone, 5Z-7-oxozeaenol, prevents inflammation by inhibiting the catalytic activity of TAK1 MAPK kinase kinase. *J. Biol. Chem.* **278**(20), 18485-18490 (2003).
2. Choo, M.-K., Kawasaki, N., Singhirunnusorn, P., *et al.* Blockade of transforming growth factor- β -activated kinase 1 activity enhances TRAIL-induced apoptosis through activation of a caspase cascade. *Mol. Cancer Ther.* **5**(12), 2970-2976 (2006).
3. Place, D.E., Samir, P., Malireddi, R.K.S., *et al.* Integrated stress response restricts macrophage necroptosis. *Life Sci. Alliance* **5**(1), e202101260 (2021).

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, 08/08/2023

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