

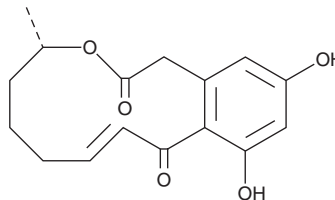
PRODUCT INFORMATION



10,11-dehydro Curvularin

Item No. 21801

CAS Registry No.: 21178-57-4
Formal Name: (4S,8E)-4,5,6,7-tetrahydro-11,13-dihydroxy-4-methyl-2H-3-benzoxacyclododecin-2,10(1H)-dione
Synonym: α,β -dehydro Curvularin
MF: C₁₆H₁₈O₅
FW: 290.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 226, 298 nm
Supplied as: A crystalline 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

10,11-dehydro Curvularin is supplied as a crystalline solid. A stock solution may be made by dissolving the 10,11-dehydro curvularin in the solvent of choice. 10,11-dehydro Curvularin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of 10,11-dehydro curvularin in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

10,11-dehydro Curvularin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 10,11-dehydro curvularin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 10,11-dehydro Curvularin has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

10,11-dehydro Curvularin is a natural mycotoxin that has cytotoxic activity against select cancer cell lines.^{1,2} It has been shown to activate the heat shock response, block TGF- β signaling, and scavenge superoxide anions.^{1,3,4}

References

1. Kumar, C.G., Mongolla, P., Sujitha, P., *et al.* *Springerplus* **2**(1), 122 (2013).
2. Zhan, J., Wijeratne, E.M., Seliga, C.J., *et al.* *J. Antibiot. (Tokyo)* **57**(5), 341-344 (2004).
3. Rudolph, K., Serwe, A., and Erkel, G. *Nat. Rev. Drug Discov.* **8**(1), 23-33 (2009).
4. Xu, Y., Espinosa-Artiles, P., Schubert, V., *et al.* *Appl. Environ. Microbiol.* **79**(6), 2038-2047 (2013).

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, 12/22/2022

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