

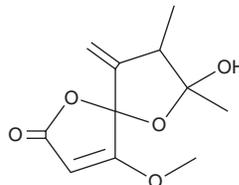
PRODUCT INFORMATION



Papyracillic Acid

Item No. 29324

CAS Registry No.: 179308-49-7
Formal Name: 7-hydroxy-4-methoxy-7,8-dimethyl-9-methylene-1,6-dioxaspiro[4.4]non-3-en-2-one
MF: C₁₁H₁₄O₅
FW: 226.2
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

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

Description

Papyracillic acid is a fungal metabolite and a derivative of penicillic acid (Item No. 11440) originally isolated from *L. papyraceum* and has antibiotic, antifungal, and phytotoxic activities.^{1,2} It is active against the bacteria *X. campestris* and *B. subtilis* and the fungus *C. tropicalis* in a disc assay when used at a concentration of 5 µg/disc.² Papyracillic acid (1 mg/ml) induces necrotic lesion formation in a panel of 10 plants.

References

1. Shan, R., Anke, H., Stadler, M., et al. Papyracillic acid, a new penicillic acid analogue from the Ascomycete *Lachnum papyraceum*. *Tetrahedron* **52(30)**, 10249-10254 (1996).
2. Evidente, A., Berestetskiy, A., Cimmino, A., et al. Papyracillic acid, a phytotoxic 1,6-dioxaspiro[4,4]nonene produced by *Ascochyta agropyrina* Var. *nana*, a potential mycoherbicide for *Elytrigia repens* biocontrol. *J. Agric. Food Chem.* **57(23)**, 11168-11173 (2009).

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, 09/14/2023

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