# **PRODUCT** INFORMATION



## 14α-hydroxy Paspalinine

Item No. 25081

CAS Registry No.: Formal Name:	151341-77-4 (3R,5bS,6S,7aR,13bS,13cR,15aS)- 2,3,5b,6,7,7a,8,13,13b,13c,14,15- dodecahydro-5b,6-dihydroxy- 2,2,13b,13c-tetramethyl-4H- 3,15a-epoxy-1-benzoxepino[6- ',7':6,7]indeno[1,2-b]indol-4-one	
Synonym:	14(S)-hydroxy Paspalinine	
MF:	$C_{27}H_{31}NO_5$	
FW:	449.5	
Purity:	≥95%	H VII
Supplied as:	A solid	
Storage:	-20°C	
Stability:	≥4 years	
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.		

ire p

### Laboratory Procedures

 $14\alpha$ -hydroxy Paspalinine is supplied as a solid. A stock solution may be made by dissolving the  $14\alpha$ -hydroxy paspalinine in the solvent of choice, which should be purged with an inert gas.  $14\alpha$ -hydroxy Paspalinine is soluble in organic solvents such as methanol, ethanol, DMSO, and dimethyl formamide.

#### Description

14 $\alpha$ -hydroxy Paspalinine is an indole diterpene fungal metabolite originally isolated from A. nomius.<sup>1</sup> Dietary administration of  $14\alpha$ -hydroxy paspalinine (100 ppm) reduces weight gain of *H. zea* first instar larvae by 91%.

#### Reference

1. Staub, G.M., Gloer, K.B., and Gloer, J.B. New paspalinine derivatives with antiinsectan activity from the sclerotia of Aspergillus nominus. Tetrahedron Lett. 34(16), 2569-2572 (1993).

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.

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, 12/19/2022

## 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