

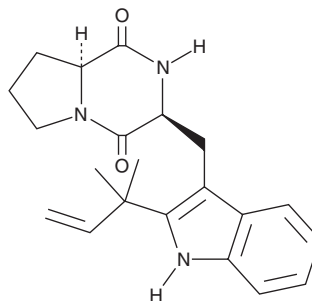
PRODUCT INFORMATION



Deoxybrevianamide E

Item No. 26662

CAS Registry No.: 34610-68-9
Formal Name: (3S,8aS)-3-[[2-(1,1-dimethyl-2-propen-1-yl)-1H-indol-3-yl]methyl]hexahydro-pyrrolo[1,2-a]pyrazine-1,4-dione
MF: C₂₁H₂₅N₃O₂
FW: 351.4
Purity: ≥99%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Penicillium* sp.



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

Laboratory Procedures

Deoxybrevianamide E is supplied as a solid. A stock solution may be made by dissolving the deoxybrevianamide E in the solvent of choice, which should be purged with an inert gas. Deoxybrevianamide E is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Deoxybrevianamide E is an alkaloid fungal metabolite that has been found in *Aspergillus*.¹ It is a precursor in the biosynthesis of notoamides.

Reference

1. Kato, H., Nakamura, Y., Finefield, J.M., *et al.* Study on the biosynthesis of the notoamides: Pinacol-type rearrangement of the isoprenyl unit in deoxybrevianamide E and 6-hydroxydeoxybrevianamide E. *Tetrahedron Lett.* **52(51)**, 6923-6926 (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, 08/23/2023

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