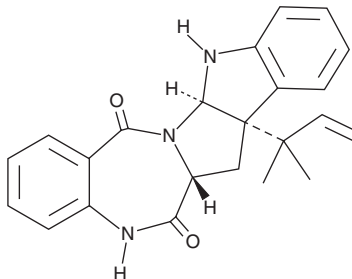


PRODUCT INFORMATION



Aszonalenin Item No. 25018

CAS Registry No.: 81797-27-5
Formal Name: (5aR,13aR,14aS)-14a-(1,1-dimethyl-2-propenyl)-5a,13a,14,14a-tetrahydro-indolo[3',2':4,5]pyrrolo[2,1-c][1,4]benzodiazepine-7,13(5H,12H)-dione
Synonym: NSC 374337
MF: C₂₃H₂₃N₃O₂
FW: 373.5
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Neosartorya spinosa*



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

Laboratory Procedures

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

Description

Aszonalenin is a hexahydropyrrolo[2,3-b]indole-containing fungal metabolite produced by *A. zonatus* and *N. tatenoi*.^{1,2}

References

1. Yasuo, K., Takashi, H., Hiromitsu, N., *et al.* Structure of aszonalenin, a new metabolite of *Aspergillus zonatus*. *Tetrahedron Lett.* **23(2)**, 225-228 (1982).
2. Yim, T., Kanokmedhakul, K., Kanokmedhakul, S., *et al.* A new meroterpenoid tatenolic acid from the fungus *Neosartorya tatenoi* KKU-2NK23. *Nat. Prod. Res.* **28(21)**, 1847-1852 (2014).

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, 10/05/2023

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