

PRODUCT INFORMATION



Paraherquamide E

Item No. 26680

CAS Registry No.: 125600-53-5
Formal Name: (1'S,5'aS,7'R,8'aS,9'aR)-2',3',8'a,9'-tetrahydro-1',4,4,8',11'-hexamethyl-spiro[4H,8H-[1,4]dioxepino[2,3-g]indole-8,7'(8'H)-[5H,6H-5a,9a](iminomethano)[1H]cyclopent[f]indolizine]-9,10'(10H)-dione

Synonym: VM 54159

MF: $C_{28}H_{35}N_3O_4$

FW: 477.6

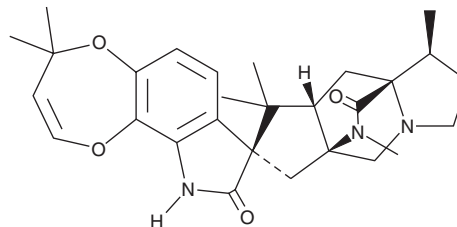
Purity: $\geq 85\%$

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

Paraherquamide E is supplied as a solid. A stock solution may be made by dissolving the paraherquamide E in the solvent of choice, which should be purged with an inert gas. Paraherquamide E is slightly soluble (0.1-1 mg/ml) in methanol.

Description

Paraherquamide E is a fungal metabolite originally isolated from *P. charlesii* with anthelmintic and insecticidal activities.¹⁻³ It is lethal to *C. elegans* ($\text{LD}_{50} = 6 \mu\text{g/ml}$).¹ Paraherquamide E is also lethal to *O. fasciatus* ($\text{LD}_{50} = 0.089 \mu\text{g/nymph}$).² Oral administration of paraherquamide E (0.5-4 mg/kg) reduces *T. colubriformis* fecal egg count in gerbils.³

References

1. Ondeyka, J.G., Goegelman, R.T., Schaeffer, J.M., et al. Novel antinematodal and antiparasitic agents from *Penicillium charlesii*. *J. Antibiot. (Tokyo)* **43(11)**, 1375-1379 (1990).
2. López-Gresa, M.P., González, M.C., Ciavatta, L., et al. Insecticidal activity of Paraherquamides, including paraherquamide H and paraherquamide I, two new alkaloids isolated from *Penicillium cluniae*. *J. Agric. Food Chem.* **54(8)**, 2921-2925 (2006).
3. Banks, R.M., Blanchflower, S.E., Everett, J.R., et al. Novel anthelmintic metabolites from an *Aspergillus* species; the aspergillimides. *J. Antibiot. (Tokyo)* **50(10)**, 840-846 (1997).

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, 03/21/2024

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