

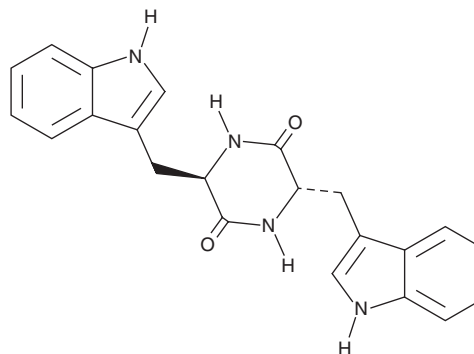
PRODUCT INFORMATION



Fellutanine A

Item No. 29320

CAS Registry No.: 175414-35-4
Formal Name: (3R,6S)-3,6-bis(1H-indol-3-ylmethyl)-2,5-piperazinedione
MF: C₂₂H₂₀N₄O₂
FW: 372.4
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/Actinomycete



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

Laboratory Procedures

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

Description

Fellutanine A is a diketopiperazine fungal metabolite that has been found in *A. candidus*.¹

Reference

1. Shan, T., Wang, Y., Wang, S., *et al.* A new *p*-terphenyl derivative From the insect-derived fungus *Aspergillus candidus* Bdf-2 and the synergistic effects of terphenyllin. *Peer J.* **8**, e8221 (2020).

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, 12/06/2022

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