

PRODUCT INFORMATION



GKK1032B

Item No. 28855

CAS Registry No.: 358375-11-8

Formal Name: (3aR,9aR,9bS,10R,12S,13aR,13bR,13cR,16R,16aS,17aS)-rel-(+)-16-ethenyl-3a,9b,10,11,12,13,13a,13b,13c,16,16a,17a-dodecahydro-10,12,13a,14,16-pentamethyl-5,8-ethenofluoreno[9',1':2,3,4]oxacyclododecino[6,7-c]pyrrole-1,3,17(2H,4H,9aH)-trione

MF: $C_{32}H_{39}NO_4$

FW: 501.7

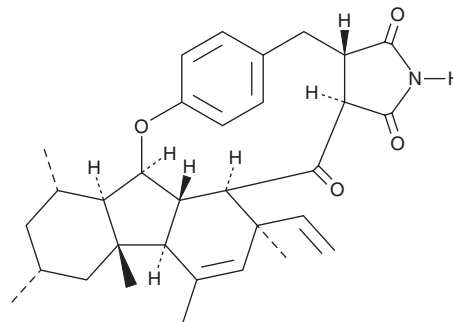
Purity: $\geq 70\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

GKK1032B is supplied as a solid. A stock solution may be made by dissolving the GKK1032B in the solvent of choice, which should be purged with an inert gas. GKK1032B is soluble in organic solvents such as ethanol, DMSO, and dichloromethane.

Description

GKK1032B is a fungal metabolite originally isolated from *Penicillium* sp. GKK1032 that has antiproliferative and antibacterial activities.^{1,2} GKK1032B inhibits the growth of HeLa S3 cervical and MCF-7 breast cancer cells (IC_{50} s = 17.7 and 14.71 μM , respectively) and Vero cells (IC_{50} = 29.55 μM). It also inhibits the growth of *B. subtilis* (MIC = 20.8 $\mu\text{g/ml}$) and *M. tuberculosis* (MIC = 48.35 μM).

References

1. Li, X.-W., Ear, A., and Nay, B. Hirsutellones and beyond: Figuring out the biological and synthetic logics toward chemical complexity in fungal PKS-NRPS compounds. *Nat. Prod. Rep.* **30**(6), 765-782 (2013).
2. Rukachaisirikul, V., Satpradit, S., Klaiklay, S., et al. Polyketide anthraquinone, diphenyl ether, and xanthone derivatives from the soil fungus *Penicillium* sp. PSU-RSPG99. *Tetrahedron* **70**(34), 5148-5152 (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, 08/16/2023

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