

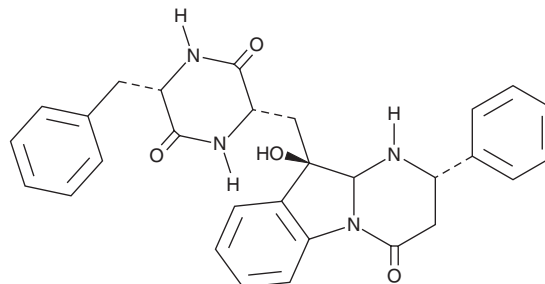
PRODUCT INFORMATION



Citreoindole

Item No. 22069

CAS Registry No.: 138655-14-8
Formal Name: (3S,6S)-3-[[[(2S,10R,10aR)-1,2,3,4,10,10a-hexahydro-10-hydroxy-4-oxo-2-phenylpyrimido[1,2-a]indol-10-yl]methyl]-6-(phenylmethyl)-2,5-piperazinedione
MF: C₂₉H₂₈N₄O₄
FW: 496.6
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

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

Description

Citreoindole is a diketopiperazine metabolite isolated from a hybrid cell fusion of two strains of *P. citreovirde* that is cytotoxic *in vitro* against HeLa cells at 8.4 μM.^{1,2}

References

1. Matsunaga, K., Shizuri, Y., Yamamura, S., *et al.* Isolation and structure of citreoindole, a new metabolite of hybrid strain KO 0052 derived from *Penicillium citreo-viride* B. IFO 6200 and 4692. *Tetrahedron Lett.* **32(47)**, 6883-6884 (1991).
2. Song, F., He, H., Ma, R., *et al.* Structure revision of the *Penicillium* alkaloids haenamindole and citreoindole. *Tetrahedron Lett.* **57(24)**, 3851-3852 (2016).

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, 11/14/2022

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