

PRODUCT INFORMATION

Herquiline A

Item No. 27731

CAS Registry No.: 71812-08-3

Formal Name: (2R,5R,11S,11aS,14aS,14bR)-rel-(-)-1,2,5,6,8,9,11,11a,13,14,14a,14b-dodecahydro-16-methyl-4H-5,2-(iminomethano)-11,7-methenopyrrolo[3,2,1-no][1]benzazacycloundecine-10,12-dione

MF: $C_{19}H_{26}N_2O_2$

FW: 314.4

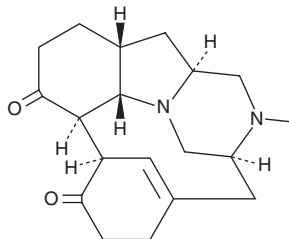
Purity: $\geq 98\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Fungus/*Penicillium herquei*



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

Laboratory Procedures

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

Description

Herquiline A is an alkaloid fungal metabolite originally isolated from *P. herquei*.¹ It inhibits cell death induced by influenza A strain A/PR/8/34 in MDCK cells with an IC_{50} value of $10\text{ }\mu\text{g/ml}$ and inhibits viral replication in a plaque assay in a concentration-dependent manner.² Herquiline A also inhibits platelet aggregation induced by ADP (Item No. 16778) and platelet-activating factor (PAF) *in vitro* (IC_{50} s = 180 and $240\text{ }\mu\text{M}$, respectively).¹

References

- Enomoto, Y., Shiomi, K., Hayashi, M., *et al.* Herquiline B, a new platelet aggregation inhibitor produced by *Penicillium herquei* Fg-372. *J. Antibiot. (Tokyo)* **49**(1), 50-53 (1996).
- Chiba, T., Asami, Y., Suga, T., *et al.* Herquiline A, produced by *Penicillium herquei* FKI-7215, exhibits anti-influenza virus properties. *Biosci. Biotechnol. Biochem.* **81**(1), 59-62 (2017).

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, 08/10/2023

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