

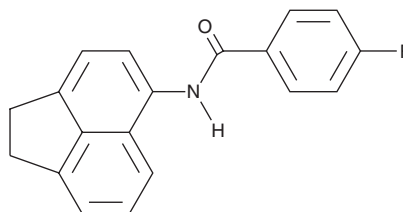
PRODUCT INFORMATION



Questin

Item No. 40904

CAS Registry No.: 3774-64-9
Formal Name: 1,6-dihydroxy-8-methoxy-3-methyl-9,10-anthracenedione
MF: $C_{16}H_{12}O_5$
FW: 284.3
Purity: $\geq 70\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungus/*Aspergillus terreus*



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

Laboratory Procedures

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

Description

Questin is an anthraquinone that has been found in *A. fumigatus* and has diverse biological activities.¹⁻⁴ It inhibits cell division cycle 25B (Cdc25B) phosphatase ($IC_{50} = 34 \mu\text{g/ml}$) and reduces the proliferation of SW620 colon cancer cells ($GI_{50} = 0.9 \mu\text{g/ml}$).¹ Questin (50 or 100 μM) reduces the viability of A549 lung cancer cells.² It reduces viral replication in MDCK cells infected with influenza A H1N1 or H3N2 (IC_{50} s = 1.92 and 9.62 μM , respectively).³ Questin decreases proliferation of the protozoa *T. fetus* ($MIC = 12.5 \mu\text{g/ml}$).⁴

References

1. Choi, S.-G. and Kim, J. Anthraquinones, Cdc25B phosphatase inhibitors, isolated from the roots of *Polygonum multiflorum* Thunb. *Nat. Prod. Res.* **21**(6), 487-493 (2007).
2. Gauthier, T., Wang, X., Dos Santos, J.S., et al. Trypacidin, a spore-borne toxin from *Aspergillus fumigatus*, is cytotoxic to lung cells. *PLoS One* **7**(2), e29906 (2012).
3. Liu, N., Peng, S., Yang, J., et al. Structurally diverse sesquiterpenoids and polyketides from a sponge-associated fungus *Aspergillus sydowii* SCSIO41301. *Fitoterapia* **135**, 27-32 (2019).
4. Chaudhary, N.K., Pitt, J.I., Lacey, E., et al. Banksialactones and banksiamarins: Isochromanones and isocoumarins from an Australian fungus, *Aspergillus banksianus*. *J. Nat. Prod.* **81**(7), 1517-1526 (2018).

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, 05/16/2024

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