

PRODUCT INFORMATION



Tryptoquivaline D

Item No. 33531

CAS Registry No.: 60676-56-4
Formal Name: (2S,2'S,4R,9'aS)-4-[2-[(1S)-1-(acetyloxy)-2-methylpropyl]-4-oxo-3(4H)-quinazoliny]-1',3,4,9'a-tetrahydro-1'-hydroxy-2'-methyl-spiro[furan-2(5H),9'-[9H]imidazo[1,2-a]indole]-3',5(2'H)-dione
Synonyms: Nortryptoquivaline, NSC 292204

MF: C₂₈H₂₈N₄O₇

FW: 532.5

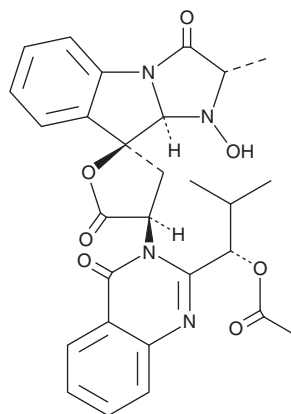
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungi/*Aspergillus fumigatus*



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

Laboratory Procedures

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

Description

Tryptoquivaline D is a fungal metabolite that has been found in *Neosartorya siamensis* and has anticancer activity.^{1,2} It induces nuclear chromatin condensation, a marker of apoptosis, in HCT116 colon and HepG2 liver cancer cells when used at a concentration of 150 μM.¹ Tryptoquivaline D (1-100 μM), alone or in combination with doxorubicin (Item No. 15007), reduces the viability of A549 lung cancer cells.²

References

1. Prata-Sena, M., Ramos, A.A., Buttachon, S., *et al.* Cytotoxic activity of secondary metabolites from marine-derived fungus *Neosartorya siamensis* in human cancer cells. *Phytother. Res.* **30(11)**, 1862-1871 (2016).
2. Ramos, A.A., Castro-Carvalho, B., Prata-Sena, M., *et al.* Can marine-derived fungus *Neosartorya siamensis* KUFA 0017 extract and its secondary metabolites enhance antitumor activity of doxorubicin? An in vitro survey unveils interactions against lung cancer cells. *Environ. Toxicol.* **35(4)**, 507-517 (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, 11/15/2022

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