

PRODUCT INFORMATION



Quarfloxin

Item No. 36552

CAS Registry No.: 865311-47-3
Formal Name: 5-fluoro-N-[2-[(2S)-1-methyl-2-pyrrolidinyl]ethyl]-3-oxo-6-[3-(2-pyrazinyl)-1-pyrrolidinyl]-3H-benzo[b]pyrido[3,2,1-k]phenoxazine-2-carboxamide

Synonyms: CX-3543, Itarnaflloxin, Quarfloxacin

MF: C₃₅H₃₃FN₆O₃

FW: 604.7

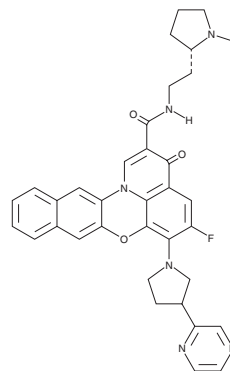
Purity: ≥98%

UV/Vis.: λ_{max}: 218, 319 nm

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

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

Description

Quarfloxin is an inhibitor of RNA polymerase I-mediated transcription.¹ It binds to G-quadruplex DNA and inhibits the interaction between ribosomal DNA G-quadruplexes and nucleolin, a nucleolar protein required for RNA polymerase I-mediated transcription. Quarfloxin (10 μM) induces nucleolin mislocalization from the nucleoli to the nucleoplasm in A549 cells. It induces apoptosis in A549 and HL-60 cells when used at a concentration of 5 μM. Quarfloxin (5 mg/kg) reduces tumor growth in MDA-MB-231 breast and MiaPaCa-2 pancreatic cancer mouse xenograft models.

Reference

1. Drygin, D., Siddiqui-Jain, A., O'Brien, S., *et al.* Anticancer activity of CX-3543: A direct inhibitor of rRNA biogenesis. *Cancer Res.* **69**(19), 7653-7661 (2009).

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, 06/12/2023

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