

PRODUCT INFORMATION



ARN24139

Item No. 31062

CAS Registry No.: 2699768-78-8
Formal Name: 6-hydroxy-4-oxo-1,3-diphenyl-2-thioxo-N-(3-(trifluoromethoxy)phenyl)-1,2,3,4-tetrahydropyrimidine-5-carboxamide

MF: C₂₄H₁₆F₃N₃O₄S

FW: 499.5

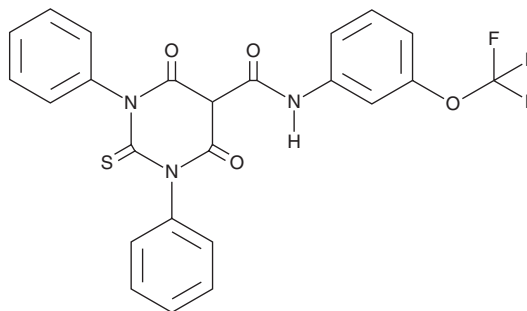
Purity: ≥98%

UV/Vis.: λ_{max}: 239, 321 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

ARN24139 is supplied as a crystalline solid. A stock solution may be made by dissolving the ARN24139 in the solvent of choice, which should be purged with an inert gas. ARN24139 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of ARN24139 in these solvents is approximately 10 and 30 mg/ml, respectively.

ARN24139 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ARN24139 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ARN24139 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ARN24139 is a topoisomerase II poison (IC₅₀ = 7.3 μM in a topoisomerase II decatenation assay).¹ It inhibits proliferation of DU145, HeLa, and A549 cells (IC₅₀s = 4.7, 3.8, and 3.1 μM, respectively).

Reference

1. Arencibia, J.M., Brindani, N., Franco-Ulloa, S., *et al.* Design, synthesis, dynamic docking, biochemical characterization, and *in vivo* pharmacokinetics studies of novel topoisomerase II poisons with promising antiproliferative activity. *J. Med. Chem.* **63(7)**, 3508-3521 (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, 01/07/2022

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