

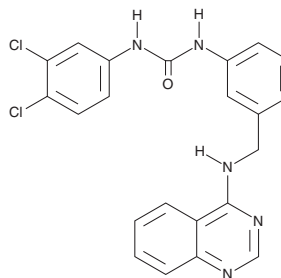
PRODUCT INFORMATION



CAY10773

Item No. 31806

CAS Registry No.: 1648546-79-5
Formal Name: N-(3,4-dichlorophenyl)-N'-[3-[(4-quinazolinylamino)methyl]phenyl]-urea
MF: C₂₂H₁₇Cl₂N₅O
FW: 438.3
Purity: ≥98%
UV/Vis.: λ_{max}: 264 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

CAY10773 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10773 in the solvent of choice, which should be purged with an inert gas. CAY10773 is soluble in DMSO and dimethyl formamide at a concentration of approximately 1 mg/ml.

Description

CAY10773 is a derivative of the ferroptosis inducer sorafenib (Item No. 10009644).¹ It selectively inhibits proliferation of BEL-7402, MGC803, and T24 bladder cancer cells (IC₅₀s = 5.77, 5.21 and 3.97 μM, respectively) over non-cancerous HCV-29 cells (IC₅₀ = 23.19 μM). CAY10773 induces apoptosis in T24 cells when used at concentrations ranging from 2 to 6 μM but induces ferroptosis at concentrations greater than or equal to 6 μM with 10 hour or longer incubation times. It increases the production of reactive oxygen species (ROS) and decreases the mitochondrial membrane potential in T24 cells. CAY10773 (≥6 μM) also induces autophagy with incubation times longer than eight hours.

Reference

1. Chen, J.-N., Li, T., Cheng, L., *et al.* Synthesis and *in vitro* anti-bladder cancer activity evaluation of quinazoliny-arylurea derivatives. *Eur. J. Med. Chem.* **205**, 112661 (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, 11/18/2020

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