

PRODUCT INFORMATION

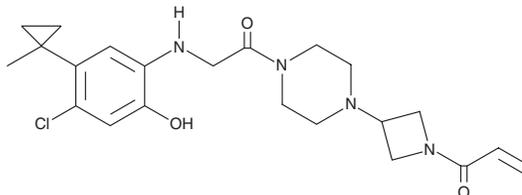


ARS853

Item No. 19137

CAS Registry No.: 1629268-00-3
Formal Name: 1-[3-[4-[2-[[4-chloro-2-hydroxy-5-(1-methylcyclopropyl)phenyl]amino]acetyl]-1-piperazinyl]-1-azetidiny]-2-propen-1-one

MF: C₂₂H₂₉ClN₄O₃
FW: 432.9
Purity: ≥95%
UV/Vis.: λ_{max}: 211, 249 nm
Supplied as: A crystalline 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

ARS853 is supplied as a crystalline solid. A stock solution may be made by dissolving the ARS853 in the solvent of choice, which should be purged with an inert gas. ARS853 is soluble in the organic solvent DMSO at a concentration of approximately 50 mg/ml.

Description

ARS853 is an inhibitor of K-RAS^{G12C}, a mutant form of K-RAS that accumulates in the active GTP-bound state in certain cancer cells.¹ ARS853 binds to the inactive GDP-bound state of K-RAS^{G12C}. It decreases the level of active GTP-bound K-RAS by 95% when used at a concentration of 10 μM in K-RAS^{G12C} mutant lung cancer cells. It also decreases phosphorylation of c-RAF, ERK, and Akt. ARS853 inhibits proliferation of six lung cancer cell lines containing the K-RAS^{G12C} mutation, including H358 cells (IC₅₀ = 2.5 μM), but does not inhibit proliferation of PC-9 cells, which contain wild-type K-RAS.

Reference

1. Lito, P., Solomon, M., Li, L.-S., *et al.* Allele-specific inhibitors inactivate mutant KRAS G12C by a trapping mechanism. *Science* **351**(6273), 604-608 (2016).

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, 11/04/2022

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