

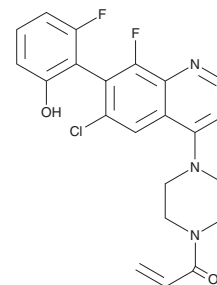
PRODUCT INFORMATION



ARS1620

Item No. 27915

CAS Registry No.: 1698055-85-4
Formal Name: 1-[4-[(7S)-6-chloro-8-fluoro-7-(2-fluoro-6-hydroxyphenyl)-4-quinazoliny]-1-piperazinyl]-2-propen-1-one
MF: C₂₁H₁₇ClF₂N₄O₂
FW: 430.8
Purity: ≥98%
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

ARS1620 is supplied as a solid. A stock solution may be made by dissolving the ARS1620 in the solvent of choice, which should be purged with an inert gas. ARS1620 is soluble in DMSO.

Description

ARS1620 is a covalent inhibitor of K-RAS^{G12C} (IC₅₀ = <3 μM).¹ It selectively inhibits growth of H358, MIA-PaCa2, and LU65 cancer cells that express K-RAS^{G12C} (IC₅₀s = 150 nM) over H441, A549, and HCT116 cells expressing wild-type K-RAS (IC₅₀s = >10 μM). ARS1620 (200 mg/kg) induces tumor regression in a MIA-PaCa2, but not an H441, mouse xenograft model. It also decreases tumor volume in patient-derived xenograft (PDX) mouse models expressing K-RAS^{G12C}, but not K-RAS^{G12D} or wild-type K-RAS.

Reference

1. Janes, M.R., Zhang, J., Li, L.-S., *et al.* Targeting KRAS mutant cancer cells with a covalent G12C-specific inhibitor. *Cell* **172**(3), 578-589 (2018).

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, 12/14/2022

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