

PRODUCT INFORMATION



ARV-110

Item No. 39662

CAS Registry No.: 2222112-77-6
Formal Name: N-[*trans*-4-(3-chloro-4-cyanophenoxy)cyclohexyl]-6-[4-[[4-[2-(2,6-dioxo-3-piperidinyl)-6-fluoro-2,3-dihydro-1,3-dioxo-1H-isoindol-5-yl]-1-piperazinyl]methyl]-1-piperidinyl]-3-pyridazinecarboxamide

Synonym: Bavdegalutamide

MF: C₄₁H₄₃ClFN₉O₆

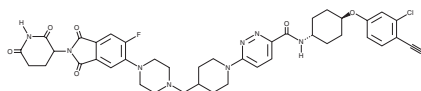
FW: 812.3

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

ARV-110 is supplied as a solid. A stock solution may be made by dissolving the ARV-110 in the solvent of choice, which should be purged with an inert gas. ARV-110 is soluble in DMSO.

Description

ARV-110 is a proteolysis-targeting chimera (PROTAC) that contains a cereblon ligand 1-derivative connected to an androgen receptor (AR) ligand by a linker region.¹ It reduces AR levels in VCaP prostate cancer cells in a time- and concentration-dependent manner. ARV-110 (1 μM) decreases VCaP cancer cell proliferation. *In vivo*, ARV-110 reduces tumor volume in a VCaP prostate cancer xenograft model that is resistant to the non-steroidal AR antagonist enzalutamide (Item No. 11596).

Reference

1. Crew, A.P., Snyder, L.B., Wang, J., *et al.* Methods of treating prostate cancer *Arvinas Operation Inc.* US20210113557A1. (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, 06/28/2024

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