

PRODUCT INFORMATION

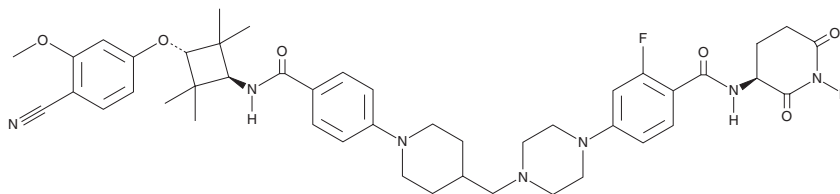


ARV-766

Item No. 41985

CAS Registry No.: 2750830-09-0
Formal Name: 4-[4-[[1-[4-[[[trans-3-(4-cyano-3-methoxyphenoxy)-2,2,4,4-tetramethylcyclobutyl]amino]carbonyl]phenyl]-4-piperidinyl]methyl]-1-piperazinyl]-N-[(3S)-2,6-dioxo-3-piperidinyl]-2-fluorobenzamide

Synonym: Luxdegalutamide
MF: C₄₅H₅₄FN₇O₆
FW: 808.0
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-766 is supplied as a solid. A stock solution may be made by dissolving the ARV-766 in the solvent of choice, which should be purged with an inert gas. ARV-766 is soluble (≥10 mg/ml) in DMSO.

Description

ARV-766 is a proteolysis-targeting chimera (PROTAC) that drives androgen receptor degradation.¹ It induces degradation of the androgen receptor in VCaP cells with a half-maximal degradation concentration (DC₅₀) value of less than 1 nM. ARV-766 inhibits tumor growth in various LNCaP and VCaP prostate cancer mouse xenograft models, including an enzalutamide-insensitive, non-castrated VCaP model.

References

1. Snyder, L., Lee, S.H., Neklesa, T.K., *et al.* Discovery of ARV-766, an androgen receptor degrading PROTAC® for the treatment of men with metastatic castration resistant prostate cancer. *Oral presentation: Proceedings of the American Association for Cancer Research Annual Meeting Part 1 (Regular and Invited Abstracts) (2023).*

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, 10/16/2024

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