

PRODUCT INFORMATION



ARD-266

Item No. 45321

CAS Registry No.: 2666951-70-6

Formal Name: (2S,4R)-N-[(1S)-3-[4-[2-[4-[[[trans-3-(3-chloro-4-cyanophenoxy)-2,2,4,4-tetramethylcyclobutyl]amino]carbonyl]phenyl]ethynyl]-1-piperidinyl]-3-oxo-1-phenylpropyl]-4-hydroxy-1-[(2R)-3-methyl-2-(3-methyl-5-isoxazolyl)-1-oxobutyl]-2-pyrrolidinecarboxamide

MF: C₅₂H₅₉ClN₆O₇

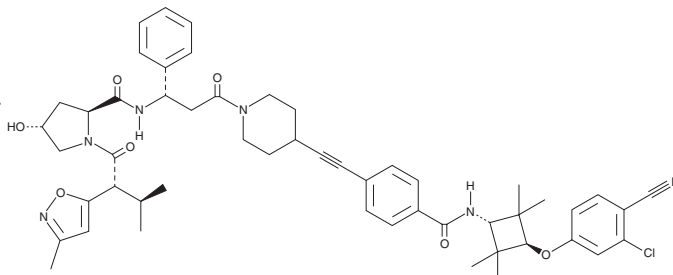
FW: 915.5

Purity: ≥95%

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

ARD-266 is supplied as a solid. A stock solution may be made by dissolving the ARD-266 in the solvent of choice, which should be purged with an inert gas. ARD-266 is slightly soluble (0.1-1 mg/ml) in acetonitrile and methanol.

Description

ARD-266 is a proteolysis-targeting chimera (PROTAC) containing an androgen receptor antagonist conjugated to a VHL ligand via a piperidine-alkyne linker.¹ It induces degradation of the androgen receptor in LNCaP, VCaP, and 22Rv1 prostate cancer cells with half-maximal degradation concentration (DC₅₀) values of 0.54, 0.97, and 0.15 nM, respectively. ARD-266 inhibits the growth of LNCaP and VCaP cells (IC₅₀s = 2 and 6 nM, respectively).

Reference

1. Han, X., Zhao, L., Xiang, W., *et al.* Discovery of highly potent and efficient PROTAC degraders of androgen receptor (AR) by employing weak binding affinity VHL E3 ligase ligands. *J. Med. Chem.* **62**(24), 11218-11231 (2019).

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, 04/21/2026

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