

PRODUCT INFORMATION



AZD 9750

Item No. 45751

CAS Registry No.: 3056515-10-4

Formal Name: 4-chloro-7-[(3S)-3-[4-[4-[1-methyl-6-(tetrahydro-2,4-dioxo-1(2H)-pyrimidinyl)-1H-indol-2-yl]-1-piperidinyl]methyl]-1-piperidinyl]phenyl]-1-piperidinyl]-1H-indazole-3-carbonitrile

MF: $C_{43}H_{48}ClN_9O_2$

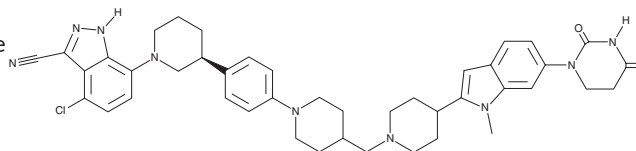
FW: 758.4

Purity: $\geq 98\%$

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

AZD 9750 is supplied as a solid. A stock solution may be made by dissolving the AZD 9750 in the solvent of choice, which should be purged with an inert gas. AZD 9750 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

AZD 9750 is a proteolysis-targeting chimera (PROTAC) composed of an androgen receptor ligand and a lenalidomide derivative.¹ It induces the degradation of wild-type androgen receptor (AR) and drug-resistant mutant AR containing a leucine-to-histidine substitution at position 702 (AR^{L702H}) with half-maximal degradation concentration (DC₅₀) values of 10 and 12 nM, respectively. AZD 9750 (30 mg/kg) decreases tumor volume in a wild-type AR-expressing patient-derived xenograft (PDX) mouse model of prostate cancer.

Reference

1. Scott, J.S., Evans, L., Astles, P.C., *et al.* Discovery of AZD9750, an orally bioavailable androgen receptor degrader for the treatment of prostate cancer. *J. Med. Chem.* **69**(3), 3209-3232 (2026).

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, 05/29/2026

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