

PRODUCT INFORMATION

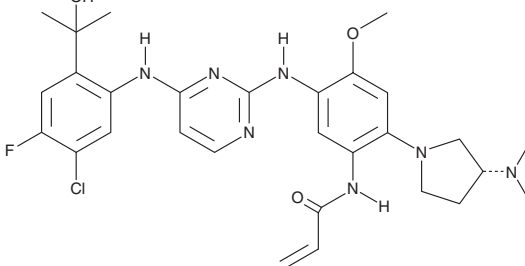


Sunvozertinib

Item No. 46085

CAS Registry No.: 2370013-12-8
Formal Name: N-[5-[[4-[[5-chloro-4-fluoro-2-(1-hydroxy-1-methylethyl)phenyl]amino]-2-pyrimidinyl]amino]-2-[(3R)-3-(dimethylamino)-1-pyrrolidinyl]-4-methoxyphenyl]-2-propenamide

Synonym: DZD9008
MF: C₂₉H₃₅ClFN₇O₃
FW: 584.1
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

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

Description

Sunvozertinib is an irreversible inhibitor of EGFR containing exon 20 insertion (ex20ins) mutations (IC₅₀ = 2.1 nM for EGFR ex20ins 770_NPG).¹ It is selective for EGFR ex20ins and EGFR^{L858R/T790M} (IC₅₀ = 13 nM) over wild-type EGFR, HER2, and HER4 (IC₅₀s = 226, 392, and 704 nM, respectively), a panel of 106 kinases at 1 μM, and other kinases, including BTK, BRK, ALK, Yes, and TrkA (IC₅₀s = 213, 529, 1,172, 2,895, and >10,000 nM, respectively). Sunvozertinib selectively inhibits the proliferation of mutant EGFR-expressing cancer cells (GI₅₀s = 6-88 nM) over those containing wild-type EGFR (GI₅₀ = 113 nM). Sunvozertinib (25 and 50 mg/kg twice per day) induces tumor regression in a patient-derived xenograft (PDX) mouse model of non-small cell lung cancer (NSCLC) carrying EGFR ex20ins insNPH. Formulations containing sunvozertinib have been used in the treatment of NSCLC with EGFR ex20ins mutations.

Reference

1. Wang, M., Yang, J.C., Mitchell, P.L., *et al.* Sunvozertinib, a selective EGFR inhibitor for previously treated non-small cell lung cancer with EGFR Exon 20 insertion mutations. *Cancer Discov.* **12**(7), 1676-1689 (2022).

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, 07/15/2026

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