

PRODUCT INFORMATION



(S)-Sunvozertinib

Item No. 38900

CAS Registry No.: 2370013-49-1

Formal Name: N-[5-[[4-[[5-chloro-4-fluoro-2-(1-hydroxy-1-methylethyl)phenyl]amino]-2-pyrimidinyl]amino]-2-[(3S)-3-(dimethylamino)-1-pyrrolidinyl]-4-methoxyphenyl]-2-propenamide

Synonym: T9304

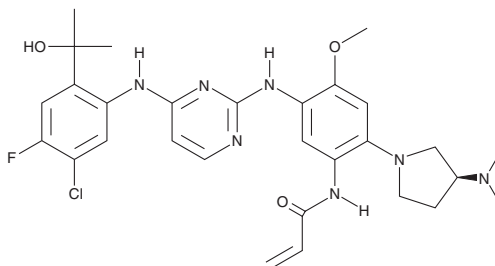
MF: $C_{29}H_{35}ClFN_7O_3$

FW: 584.1

Purity: $\geq 95\%$

Storage: -20°C

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

(S)-Sunvozertinib is supplied as a solid. A stock solution may be made by dissolving the (S)-sunvozertinib in the solvent of choice, which should be purged with an inert gas. (S)-Sunvozertinib is soluble in acetonitrile.

Description

(S)-Sunvozertinib is an inhibitor of mutant EGFR and the enantiomer of sunvozertinib (Item No. 46085).¹ It selectively inhibits the phosphorylation of EGFR^{L858R/T790M} and EGFR containing exon 20 insertion (ex20ins) mutations, including EGFR ex20ins ASV and EGFR ex20ins NPH, over wild-type EGFR (IC_{50} s = 1, 52, 51, and 362 nM, respectively). (S)-Sunvozertinib also inhibits the proliferation of PaTu 8988t pancreatic cancer cells (IC_{50} = 5.6 μM).²

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).
2. Deng, Y., Li, S., Wang, L., *et al.* Discovery of a covalent small-molecule eEF1A1 inhibitor via structure-based virtual screening. *J. Chem. Inf. Model.* **66**(2), 1083-1096 (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, 07/20/2026

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