

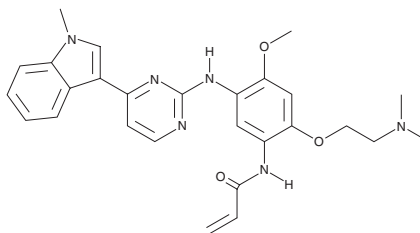
PRODUCT INFORMATION



Rezivertinib

Item No. 42537

CAS Registry No.: 1835667-12-3
Formal Name: N-[2-[2-(dimethylamino)ethoxy]-4-methoxy-5-[[4-(1-methyl-1H-indol-3-yl)-2-pyrimidinyl]amino]phenyl]-2-propenamide
Synonym: BPI-7711
MF: C₂₇H₃₀N₆O₃
FW: 486.6
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

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

Description

Rezivertinib is an inhibitor of mutant EGFR.¹ It selectively inhibits the proliferation of mutant-EGFR-containing PC-9, HCC827, and H1975 cells (GI₅₀s = 13.3, 6.8, and 22 nM, respectively) over wild-type EGFR-containing A431 cells (GI₅₀ = >1,000 nM). Rezivertinib increases survival in an H1975 brain orthotopic rat model.

Reference

1. Nagasaka, M., Zhu, V.W., Lim, S.M., *et al.* Beyond osimertinib: The development of third-generation EGFR tyrosine kinase inhibitors for advanced EGFR+ NSCLC. *J. Thorac. Oncol.* **16**(5), 740-763 (2021).

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, 12/17/2024

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