

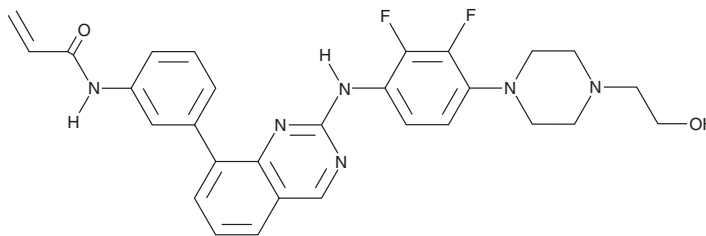
PRODUCT INFORMATION



RX-518

Item No. 40922

CAS Registry No.: 1660963-42-7
Formal Name: N-[3-[2-[[2,3-difluoro-4-[4-(2-hydroxyethyl)-1-piperazinyl]phenyl]amino]-8-quinazoliny]phenyl]-2-propenamide
Synonyms: CK-101, Olafertinib
MF: C₂₉H₂₈F₂N₆O₂
FW: 530.6
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

RX-518 is supplied as a solid. A stock solution may be made by dissolving the RX-518 in the solvent of choice, which should be purged with an inert gas. RX-518 is slightly soluble (0.1-1 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in methanol.

Description

RX-518 is an inhibitor of EGFR.¹ It inhibits the growth of A431, H1975, and HCC827 cancer cells expressing wild-type EGFR, EGFR L858R/T790M, and EGFR del19, respectively (GI₅₀s = 5, 10, and 689 nM, respectively).

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, 06/26/2024

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