

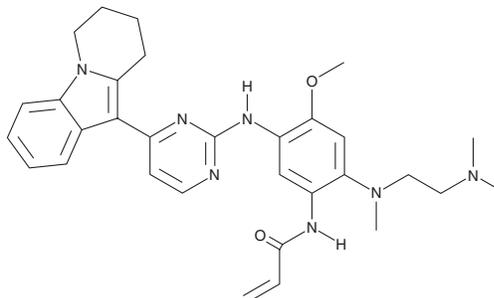
PRODUCT INFORMATION



Oritinib

Item No. 45042

CAS Registry No.: 2035089-28-0
Formal Name: N-[2-[[2-(dimethylamino)ethyl]methylamino]-4-methoxy-5-[[4-(6,7,8,9-tetrahydropyrido[1,2-a]indol-10-yl)-2-pyrimidinyl]amino]phenyl]-2-propenamide
Synonym: SH-1028
MF: C₃₁H₃₇N₇O₂
FW: 539.7
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

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

Description

Oritinib is an inhibitor of mutant EGFR.¹ It selectively inhibits EGFR^{L858R}, EGFR^{L858R/T790M}, EGFR^{Δ746-750}, and EGFR^{Δ746-750/T790M} (IC₅₀s = 0.7, 0.1, 1.4, and 0.89 nM, respectively) over wild-type EGFR (IC₅₀ = 18 nM). *In vivo*, oritinib (2.5, 5, and 15 mg/kg) inhibits tumor growth in PC-9 and H1975 mouse xenograft models expressing EGFR^{Δ746-750} and EGFR^{L858R/T790M}, respectively.

Reference

1. Han, L., Zhang, X., Wang, Z., *et al.* SH-1028, an irreversible third-generation EGFR TKI, overcomes T790M-mediated resistance in non-small cell lung cancer. *Front. Pharmacol.* **12**, (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, 02/03/2026

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