

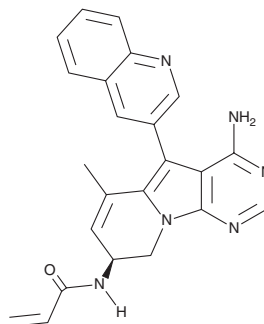
PRODUCT INFORMATION



TAS 6417

Item No. 27911

CAS Registry No.: 1661854-97-2
Formal Name: N-[(8S)-4-amino-8,9-dihydro-6-methyl-5-(3-quinolinyl)pyrimido[5,4-b]indolizin-8-yl]-2-propenamide
Synonyms: CLN-081, Zipalertinib
MF: C₂₃H₂₀N₆O
FW: 396.4
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

TAS 6417 is supplied as a solid. A stock solution may be made by dissolving the TAS 6417 in the solvent of choice, which should be purged with an inert gas. TAS 6417 is soluble in acetonitrile and DMSO.

Description

TAS 6417 is a pan-EGFR inhibitor.¹ It inhibits wild-type EGFR (IC₅₀ = 8 nM), as well as the EGFR point mutants EGFR^{L858R}, EGFR^{L861Q}, EGFR^{T790M}, EGFR^{T790M/L858R} (IC₅₀s = 1.9, 5.5, 4.1, and 2 nM, respectively) and the insertion or deletion mutants EGFR D770_N771insNPG, EGFR d746-750, and EGFR d746-770/T790M (IC₅₀s = 7.4, 1.4, and 1.1 nM, respectively). It is selective for these kinases over a panel of 20 additional kinases (IC₅₀s = 44-850 nM) but does inhibit TXK, BMX, HER4, Tec, and BTK (IC₅₀s = 1.1-22 nM). TAS 6417 inhibits the proliferation of LXF 2478L, H1975, HCC827, and PC-9 lung cancer cells (GI₅₀s = 86.5, 45.4, 1.92, and 2.96 nM, respectively). *In vivo*, TAS 6417 (50, 100, and 200 mg/kg) reduces tumor volume and increases survival in an H1975 EGFR D770-N771insSVD mouse

Reference

1. Hasako, S., Terasaka, M., Abe, N., *et al.* TAS6417, a novel EGFR inhibitor targeting exon 20 insertion mutations. *Mol. Cancer Ther.* **17**(8), 1648-1658 (2018).

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, 02/13/2024

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