

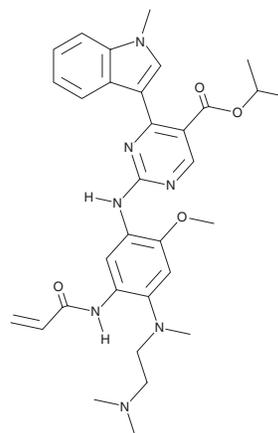
PRODUCT INFORMATION



Mobocertinib

Item No. 31522

CAS Registry No.: 1847461-43-1
Formal Name: 2-[[4-[[2-(dimethylamino)ethyl]methylamino]-2-methoxy-5-[(1-oxo-2-propen-1-yl)amino]phenyl]amino]-4-(1-methyl-1H-indol-3-yl)-5-pyrimidinecarboxylic acid, 1-methylethyl ester
Synonyms: AP32788, TAK-788
MF: C₃₂H₃₉N₇O₄
FW: 585.7
Purity: ≥98%
Supplied as: A crystalline 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

Mobocertinib is supplied as a crystalline solid. A stock solution may be made by dissolving the mobocertinib in the solvent of choice, which should be purged with an inert gas. Mobocertinib is soluble in the organic solvent dimethyl formamide (DMF) at a concentration of approximately 10 mg/ml.

Mobocertinib is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, mobocertinib should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Mobocertinib has a solubility of approximately 0.11 mg/ml in a 1:8 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Mobocertinib is an inhibitor of mutant EGFR and HER2 receptors.^{1,2} It is selective for the HER2 exon 20 mutants HER2^{V659E}, HER2^{G660D}, HER2^{G309E}, and HER2^{S310F} over wild-type EGFR in Ba/F3 cells.¹

References

1. Chouitar, J., Vincent, S., Brake, R., *et al.* TAK-788 is a novel and potent tyrosine kinase inhibitor with selective activity against EGFR/HER2. *Abstracts* **S811**, P2(13-32) (2018).
2. Neal, J., Doebele, R., Riely, G., *et al.* Safety, PK, and preliminary antitumor activity of the oral EGFR/HER2 exon 20 inhibitor TAK-788 in NSCLC. *Abstracts* **S599**, P1(13-44) (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, 10/31/2022

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