

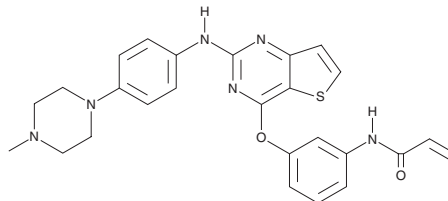
PRODUCT INFORMATION



HM61713

Item No. 19481

CAS Registry No.: 1353550-13-6
Formal Name: N-[3-[[2-[[4-(4-methyl-1-piperazinyl)phenyl]amino]thieno[3,2-d]pyrimidin-4-yl]oxy]phenyl]-2-propenamamide
Synonyms: BI-1482694, Olmutinib
MF: C₂₆H₂₆N₆O₂S
FW: 486.6
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

HM61713 is supplied as a crystalline solid. A stock solution may be made by dissolving the HM61713 in the solvent of choice, which should be purged with an inert gas. HM61713 is soluble in the organic solvent DMSO at a concentration of approximately 44 mg/ml.

Description

HM61713 is an inhibitor of mutant EGFR (IC₅₀ = 0.01 μM for EGFR^{T790M/L858R}).¹ It is selective for EGFR^{T790M/L858R} over PI3Kα (IC₅₀ = >10 μM). HM61713 is cytotoxic to A549, H1975, and H460 lung and MCF-7 breast cancer cells (IC₅₀s = 4.29, 0.52, 5.29, and 26.9 μM, respectively). It reduces tumor growth in an ETS1-overexpressing doxorubicin-resistant K562 leukemia mouse xenograft model when administered at a dose of 30 mg/kg in combination with doxorubicin.²

References

1. Hu, X., Tang, S., Yang, F., *et al.* Design, synthesis, and antitumor activity of olmutinib derivatives containing acrylamide moiety. *Molecules* **26(10)**, 3041 (2021).
2. Zhong, J., Zhang, J., Yu, X., *et al.* Olmutinib reverses doxorubicin resistance in ETS1-overexpressing leukemia cells. *Med. Sci. Monit.* **26**, e924922 (2020).

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, 10/06/2022

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