

PRODUCT INFORMATION

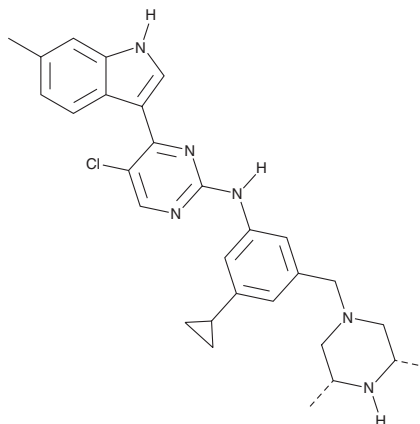


HM43239

Item No. 37737

CAS Registry No.: 2294874-49-8
Formal Name: *rel*-5-chloro-N-[3-cyclopropyl-5-[[[(3*R*,5*S*)-3,5-dimethyl-1-piperazinyl]methyl]phenyl]-4-(6-methyl-1*H*-indol-3-yl)]-2-pyrimidinamine

Synonym: Tuspetinib
MF: C₂₉H₃₃ClN₆
FW: 501.1
Purity: ≥98%
UV/Vis.: λ_{max}: 215, 283, 330 nm
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

HM43239 is supplied as a solid. A stock solution may be made by dissolving the HM43239 in the solvent of choice, which should be purged with an inert gas. HM43239 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of HM43239 in these solvents is approximately 10 mg/ml.

Description

HM43239 is an inhibitor of wild-type (WT) and internal tandem duplication (ITD) mutant FMS-related tyrosine kinase 3 (FLT3; IC₅₀s = 1.1 and 1.8 nM, respectively).¹ It also inhibits VEGFR2 and spleen tyrosine kinase (Syk) by 100 and 95%, respectively, when used at a concentration of 100 nM. HM43239 inhibits the growth of MOLM-13 and MV4-11 leukemia cells (GI₅₀s = 5.2 and 1.3 nM, respectively).

References

1. Bae, I.H., Kim, J.S., Choi, J.Y., *et al.* Pyrimidine compounds and pharmaceutical compositions for preventing or treating cancers including the same. *Hanmi Pharm. Co., LTD US 10,870,639 B2* (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, 06/12/2023

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