

PRODUCT INFORMATION

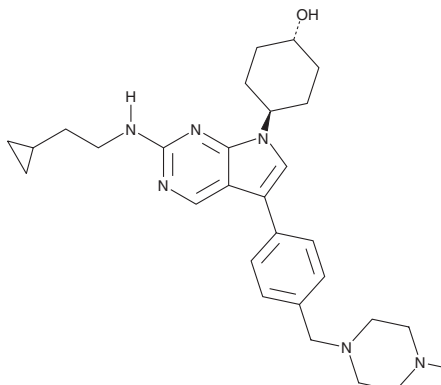


MRX-2843

Item No. 27923

CAS Registry No.: 1429882-07-4
Formal Name: *trans*-4-[2-[(2-cyclopropylethyl)amino]-5-[4-[(4-methyl-1-piperazinyl)methyl]phenyl]-7H-pyrrolo[2,3-d]pyrimidin-7-yl]-cyclohexanol

MF: C₂₉H₄₀N₆O
FW: 488.7
Purity: ≥95%
UV/Vis.: λ_{max}: 223, 255, 290 nm
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

MRX-2843 is supplied as a crystalline solid. A stock solution may be made by dissolving the MRX-2843 in the solvent of choice. MRX-2843 is soluble in the organic solvent methanol, which should be purged with an inert gas.

Description

MRX-2843 is an inhibitor of Mer and FMS-like tyrosine kinase 3 (FLT3; IC₅₀s = 1.3 and 1 nM, respectively).¹ It also inhibits Axl and Tyro3 (IC₅₀s = 15 and 17 nM, respectively). MRX-2843 (10-300 nM) inhibits Mer phosphorylation in MOLM-14 and MV4-11 acute myeloid leukemia (AML) cells and reduces clonal expansion of Kasumi-1 AML cells (IC₅₀ = 143.5 nM).² *In vivo*, MRX-2843 increases survival in NOMO-1 and MOLM-14 AML mouse xenograft models when administered at doses of 65 and 50 mg/kg, respectively.

References

1. Zhang, W., DeRyckere, D., Hunter, D., *et al.* UNC2025, a potent and orally bioavailable MER/FLT3 dual inhibitor. *J. Med. Chem.* **57(16)**, 7031-7041 (2014).
2. Minson, K.A., Smith, C.C., DeRyckere, D., *et al.* The MERTK/FLT3 inhibitor MRX-2843 overcomes resistance-conferring FLT3 mutations in acute myeloid leukemia. *JCI Insight* **1(3)**, e85630 (2016).

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, 12/19/2022

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