

PRODUCT INFORMATION

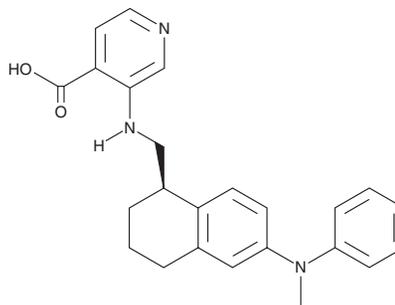


QC6352

Item No. 26231

CAS Registry No.: 1851373-36-8
Formal Name: 3-[[[(1R)-1,2,3,4-tetrahydro-6-(methylphenylamino)-1-naphthalenyl]methyl]amino]-4-pyridinecarboxylic acid

MF: C₂₄H₂₅N₃O₂
FW: 387.5
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

QC6352 is supplied as a solid. A stock solution may be made by dissolving the QC6352 in the solvent of choice, which should be purged with an inert gas. QC6352 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

QC6352 is an inhibitor of lysine demethylase 4 (KDM4).¹ It is selective for the KDM4 isoforms KDM4A, KDM4B, KDM4C, and KDM4D (IC₅₀s = 104, 56, 35, and 104 nM, respectively) over KDM2B, KDM3A, KDM3B, KDM6B, and PHD finger protein 8 (PHF8; IC₅₀s = >4,000 nM for all) but does inhibit KDM5B (IC₅₀ = 750 nM). QC6352 increases trimethylation of lysine 36 on histone H3 (H3K36Me3) in KYSE-150 esophageal carcinoma cells overexpressing KDM4C (EC₅₀ = 1.3 nM). It reduces the proliferation of KYSE-150 cells (EC₅₀ = 3.5 nM) but not IMR-90 fibroblasts (IC₅₀ = >30 μM). QC6352 (10, 25, and 50 mg/kg) reduces tumor volume in a BR0859f patient-derived xenograft (PDX) mouse model of HER2-positive breast cancer.

Reference

1. Chen, Y.K., Bonaldi, T., Cuomo, A., *et al.* Design of KDM4 inhibitors with antiproliferative effects in cancer models. *ACS Med. Chem. Lett.* **8**(8), 869-874 (2017).

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, 05/15/2025

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