

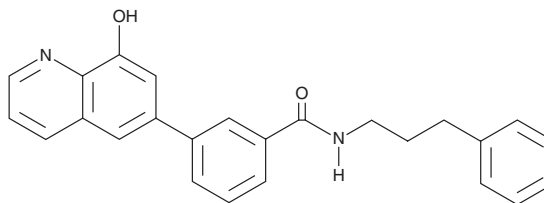
PRODUCT INFORMATION



NCGC00244536

Item No. 18478

CAS Registry No.: 2003260-55-5
Formal Name: 3-(8-hydroxy-6-quinoliny)-N-(3-phenylpropyl)-benzamide
MF: C₂₅H₂₂N₂O₂
FW: 382.5
Purity: ≥98%
UV/Vis.: λ_{max}: 278 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

NCGC00244536 is supplied as a crystalline solid. A stock solution may be made by dissolving the NCGC00244536 in the solvent of choice, which should be purged with an inert gas. NCGC00244536 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of NCGC00244536 in these solvents is approximately 20 and 30 mg/ml, respectively.

NCGC00244536 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NCGC00244536 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. NCGC00244536 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

NCGC00244536 is an inhibitor of KDM4B/JMJD2B histone lysine demethylase (IC₅₀ = ~10 nM).¹ It is selective for KDM4B/JMJD2B over KDM5A/JARID1 and lysine-specific demethylase 1 (LSD1) but does inhibit KDM4A/JMJD2A, KDM4C/JMJD2C, and KDM4D/JMJD2D at 10 μM. NCGC00244536 inhibits the growth of PC3, DU145, LNCaP, VCaP, and C4-2 prostate cancer cells (IC₅₀s = <1 μM for all). It reduces androgen receptor levels and trimethylation of histone 3 lysine 9 (H3K9) in isolated human solid prostate tumors. NCGC00244536 (20 mg/kg, s.c.) reduces tumor volume in a PC3 mouse xenograft model.

Reference

1. Duan, L., Rai, G., Roggero, C., *et al.* KDM4/JMJD2 histone demethylase inhibitors block prostate tumor growth by suppressing the expression of AR and BMYB-regulated genes. *Chem. Biol.* **22**(9), 1185-1196 (2015).

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, 12/02/2022

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