

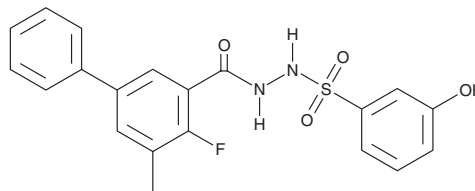
PRODUCT INFORMATION



WM-3835

Item No. 40296

CAS Registry No.: 2229025-70-9
Formal Name: 4-fluoro-5-methyl-[1,1'-biphenyl]-3-carboxylic acid, 2-[(3-hydroxyphenyl)sulfonyl]hydrazide
MF: C₂₀H₁₇FN₂O₄S
FW: 400.4
Purity: ≥95%
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

WM-3835 is supplied as a solid. A stock solution may be made by dissolving the WM-3835 in the solvent of choice, which should be purged with an inert gas. WM-3835 is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

Description

WM-3835 is an inhibitor of lysine acetyltransferase 6A (KAT6A) and KAT7 (IC₅₀s = 17 and 30 nM, respectively).¹ It is selective for KAT6A and KAT7 over KAT5 (IC₅₀ = 312 nM). WM-3835 (1 μM) reduces the viability of primary mouse leukemia stem cells (LSCs), MOLM-13 and OCI-AML-3 acute myeloid lymphoma (AML) cells, as well as decreases colony formation in patient-derived primary AML cells. It induces cell cycle arrest at the sub-G₁ phase and apoptosis, as well as decreases the levels of mRNA encoding homeobox protein A7 (HOXA7), in MOLM-13 cells. *In vivo*, WM-3835 (5 mg/kg per day) reduces tumor volume, weight, and growth rate, as well as decreases the acetylation at lysine 14 on histone 3 (H3K14), H3K12, and H3K5 in tumor cell lysates, in a PPC-1 prostate cancer mouse xenograft model.²

References

- MacPherson, L., Anokye, J., Yeung, M.M., *et al.* HBO1 is required for the maintenance of leukaemia stem cells. *Nature* **577**(7789), 266-270 (2020).
- Mi, Y.-Y., Ji, Y., Zhang, L., *et al.* A first-in-class HBO1 inhibitor WM-3835 inhibits castration-resistant prostate cancer cell growth in vitro and in vivo. *Cell Death Dis.* **14**(1), 67 (2023).

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, 03/25/2024

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