

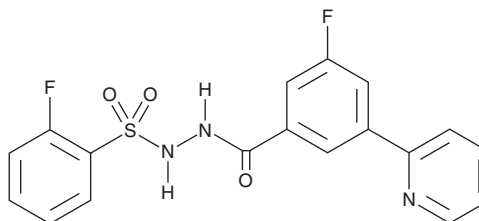
PRODUCT INFORMATION



WM-1119

Item No. 30509

CAS Registry No.: 2055397-28-7
Formal Name: 3-fluoro-5-(2-pyridinyl)-benzoic acid, 2-[(2-fluorophenyl)sulfonyl]hydrazide
MF: C₁₈H₁₃F₂N₃O₃S
FW: 389.4
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 277 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

WM-1119 is supplied as a crystalline solid. A stock solution may be made by dissolving the WM-1119 in the solvent of choice, which should be purged with an inert gas. WM-1119 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of WM-1119 in these solvents is approximately 30 mg/ml.

WM-1119 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, WM-1119 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. WM-1119 has a solubility of approximately 0.25 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

WM-1119 is an orally bioavailable inhibitor of lysine acetyltransferase 6A (KAT6A/MOZ; IC₅₀ = 6.3 nM).¹ It binds to KAT6A with a K_d value of 2 nM and is selective for KAT6A over KAT5 and KAT7 (K_ds = 2,200 and 500 nM, respectively), as well as over a panel of 159 additional targets at 1 and 10 μM.² WM-1119 inhibits proliferation of EMRK1184 B cell lymphoma cells (IC₅₀ = 0.25 μM). It reduces tumor growth and spleen weight in a murine EMRK1184 model when administered at a dose of 50 mg/kg four times per day.

References

1. Priebbenow, D.L., Leaver, D.J., Nguyen, N., *et al.* Discovery of acylsulfonohydrazide-derived inhibitors of the lysine acetyltransferase, KAT6A, as potent senescence-inducing anti-cancer agents. *J. Med. Chem.* **63**(9), 4655-4684 (2020).
2. Baell, J.B., Leaver, D.J., Hermans, S.J., *et al.* Inhibitors of histone acetyltransferases KAT6A/B induce senescence and arrest tumour growth. *Nature* **560**(7717), 253-257 (2018).

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, 12/12/2022

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