

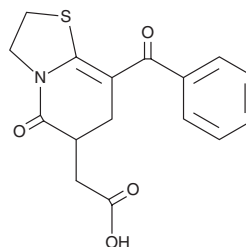
PRODUCT INFORMATION



193 D7

Item No. 43718

CAS Registry No.: 861224-48-8
Formal Name: 8-benzoyl-2,3,6,7-tetrahydro-5-oxo-5H-thiazolo[3,2-a]pyridine-6-acetic acid
MF: C₁₆H₁₅NO₄S
FW: 317.4
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

193 D7 is supplied as a solid. A stock solution may be made by dissolving the 193 D7 in the solvent of choice, which should be purged with an inert gas. 193 D7 is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

Description

193 D7 is an inhibitor of the histone demethylase jumonji domain-containing 1C (JMJD1C).¹ It inhibits JMJD1C demethylation activity *in vitro* (IC₅₀ = 0.59 μM). It is selective for JMJD1C over 14 other histone demethylases, including JMJD1A and -1B (IC₅₀s = >100 μM for both), JMJD2A, -2B, -2C, -2D, and -2E (IC₅₀s = 24.7->100 μM), JMJD3 (IC₅₀ = 6.59 μM), JARID1A, -1B, and -1C (IC₅₀s = 14.61->100 μM), and UTX, PHF8, and LSD1 (IC₅₀s = 18.4->100 μM). 193 D7 (25 mg/kg per day) reduces tumor growth in mice bearing multiple types of subcutaneous tumors established using MCA-205 fibrosarcoma, B16/F10 melanoma, Hepa 1-6 hepatocellular, LLC lung, and CT26 colon cancer cells. It also reduces intratumoral levels of regulatory T cells (Tregs) and increases intratumoral levels of CD4⁺ and CD8⁺ tumor-infiltrating effector T cells in the same model.

Reference

1. Long, X., Zhang, S., Wang, Y., *et al.* Targeting JMJD1C to selectively disrupt tumor Treg cell fitness enhances antitumor immunity. *Nat. Immunol.* **25**(3), 525-536 (2024).

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, 07/31/2025

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