

PRODUCT INFORMATION

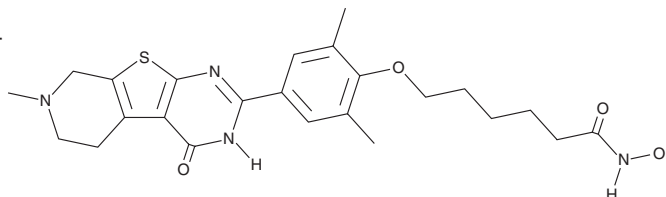


CAY17c

Item No. 31077

CAS Registry No.: 2414373-11-6
Formal Name: 6-[4-(1,4,5,6,7,8-hexahydro-7-methyl-4-oxopyrido[4',3':4,5]thieno[2,3-d]pyrimidin-2-yl)-2,6-dimethylphenoxy]-N-hydroxy-hexanamide

MF: C₂₄H₃₀N₄O₄S
FW: 470.6
Purity: ≥98%
UV/Vis.: λ_{max}: 335 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

CAY17c is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY17c in the solvent of choice, which should be purged with an inert gas. CAY17c is slightly soluble in DMSO and dimethyl formamide.

Description

CAY17c is an inhibitor of bromodomain-containing protein 4 (BRD4; IC₅₀ = 0.71 μM), as well as class I histone deacetylases (HDACs; IC₅₀s = 0.046, 0.058, 0.075, and 0.167 μM for HDAC1, -2, -3, and -8, respectively) and class IIb HDACs (IC₅₀s = 0.073 and 0.923 μM for HDAC6 and HDAC10, respectively).¹ It is selective for these enzymes over BRD2, -3, and -T (IC₅₀s = >20 μM for all), as well as over HDAC4, -5, -7, -9, and -11 (IC₅₀s = >10 μM for all). CAY17c inhibits the proliferation of HCT116, SW620, and DLD-1 colorectal cancer cells (IC₅₀s = 0.45, 1.78, and 2.11 μM, respectively), as well as induces apoptosis and autophagy in HCT116 cells. It reduces tumor growth in an HCT116 mouse xenograft model when administered at doses of 15 and 30 mg/kg.

Reference

1. Pan, Z., Li, X., Wang, Y., *et al.* Discovery of thieno[2,3-d]pyrimidine-based hydroxamic acid derivatives as bromodomain-containing protein 4/histone deacetylase dual inhibitors induce autophagic cell death in colorectal carcinoma cells. *J. Med. Chem.* **63**(7), 3678-3700 (2020).

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, 08/22/2025

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