

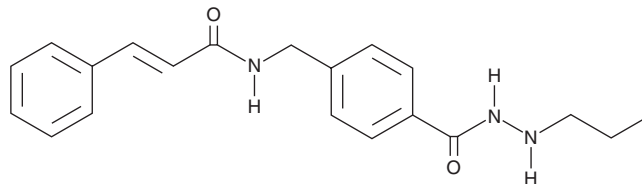
PRODUCT INFORMATION



HDAC3 Inhibitor

Item No. 21057

CAS Registry No.: 2044701-99-5
Formal Name: 4-[[[(2E)-1-oxo-3-phenyl-2-propen-1-yl]amino]methyl]-benzoic acid 2-propylhydrazide
Synonym: Histone Deacetylase 3
MF: C₂₀H₂₃N₃O₂
FW: 337.4
Purity: ≥95%
UV/Vis.: λ_{max}: 217, 223, 274 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

HDAC3 inhibitor is supplied as a solid. A stock solution may be made by dissolving the HDAC3 inhibitor in the solvent of choice, which should be purged with an inert gas. HDAC3 inhibitor is soluble in ethanol, DMSO, and dimethyl formamide.

Description

HDAC3 inhibitor is an allosteric inhibitor of histone deacetylase 3 (HDAC3; K_i = 0.16 nM).¹ It is selective for HDAC3 over HDAC1 and HDAC2 (IC₅₀s = 0.95, 11.81, and 95.45 nM, respectively, using recombinant HDACs). In addition, it is selective for acute myeloid, monocytic, and lymphoblastic leukemia cell lines (EC₅₀s = 36.37, 76.64, and 151.7 nM, respectively) over chronic myeloid, acute promyelocytic, and acute lymphoblastic cells (EC₅₀s = 2,160, >10,000, and >10,000 nM, respectively).

Reference

1. McClure, J.J., Zhang, C., Inks, E.S., *et al.* Development of allosteric hydrazide-containing class I histone deacetylase inhibitors for use in acute myeloid leukemia. *J. Med. Chem.* **59**(21), 9942-9959 (2016).

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, 02/01/2024

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