

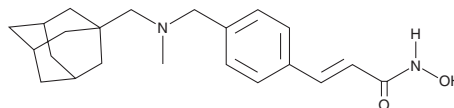
PRODUCT INFORMATION



Martinostat

Item No. 42100

CAS Registry No.: 1629052-58-9
Formal Name: N-hydroxy-3-[4-[[methyl(tricyclo[3.3.1.1^{3,7}]
dec-1-ylmethyl)amino]methyl]phenyl]-2E-
propenamide
MF: C₂₂H₃₀N₂O₂
FW: 354.5
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

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

Description

Martinostat is an inhibitor of histone deacetylase 1 (HDAC1), HDAC2, HDAC3, and HDAC6 (IC₅₀s = 0.3, 2, 0.6, and 4.1 nM, respectively).¹ It is selective for these HDACs over HDAC4, -5, -7, -8, and -9 (IC₅₀s = 1,970, 352, >20,000, >15,000, and >15,000 nM, respectively). Martinostat reduces lysine 9 on histone 3 acetylation (H3K9ac) and H4K12ac in primary mouse neuronal cells (EC₅₀s = 100 nM for both).

Reference

1. Wang, C., Schroeder, F.A., Wey, H.Y., *et al.* In vivo imaging of histone deacetylases (HDACs) in the central nervous system and major peripheral organs. *J. Med. Chem.* **57**(19), 7999-8009 (2014).

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, 11/11/2024

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