

PRODUCT INFORMATION

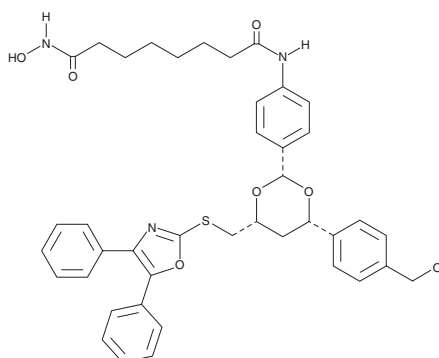


Tubacin

Item No. 13691

CAS Registry No.: 537049-40-4
Formal Name: N1-[4-[(2R,4R,6S)-4-[[[(4,5-diphenyl-2-oxazolyl)thio]methyl]-6-[4-(hydroxymethyl)phenyl]-1,3-dioxan-2-yl]phenyl]-N8-hydroxy-octanediamide

MF: C₄₁H₄₃N₃O₇S
FW: 721.9
Purity: ≥90%
UV/Vis.: λ_{max}: 222, 240, 299 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

Tubacin is supplied as a crystalline solid. A stock solution may be made by dissolving the tubacin in the solvent of choice, which should be purged with an inert gas. Tubacin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of tubacin in these solvents is approximately 16 and 20 mg/ml, respectively.

Tubacin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, tubacin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Tubacin has a solubility of approximately 0.25 mg in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Tubacin is a tubulin acetylation inducer that selectively inhibits histone deacetylase (HDAC) 6 (IC₅₀ = 4 nM).¹ In comparison, it demonstrates at least 300-fold selectivity against all other HDAC isoforms (IC₅₀s range from 1.3-17.3 μM).¹ Tubacin induces α-tubulin hyperacetylation at 2.5 μM in primary cortical neuron cultures.¹ Tubacin inhibition of HDAC6 has been used as a biochemical tool to control microtubule-dependent intracellular trafficking, to manipulate the aggresome formation of misfolded proteins in certain diseases, and to study the dynamics of cellular adhesion.²⁻⁴

References

1. Butler, K.V., Kalin, J., Brochier, C., *et al.* Rational design and simple chemistry yield a superior, neuroprotective HDAC6 inhibitor, Tubastatin A. *J. Am. Chem. Soc.* **132(31)**, 10842-10846 (2010).
2. Chen, S., Owens, G.C., Makarenkova, H., *et al.* HDAC6 regulates mitochondrial transport in hippocampal neurons. *PLoS One* **5(5)**, (2010).
3. Cole, P.A. Chemical probes for histone-modifying enzymes. *Nat. Chem. Biol.* **4(10)**, 590-597 (2008).
4. Xu, K., Dai, X.L., Huang, H.C., *et al.* Targeting HDACs: A promising therapy for Alzheimer's disease. *Oxid. Med. Cell. Longev.* 143269 (2011).

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, 10/26/2022

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