

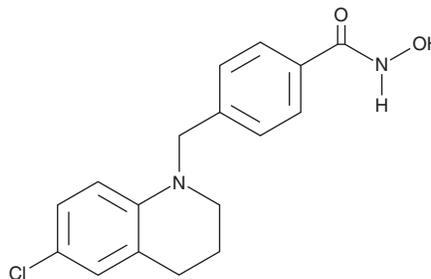
PRODUCT INFORMATION



SW-100

Item No. 33776

CAS Registry No.: 2126744-35-0
Formal Name: 4-[[6-chloro-3,4-dihydro-1(2H)-quinolinyl]methyl]-N-hydroxy-benzamide
MF: C₁₇H₁₇ClN₂O₂
FW: 316.8
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

SW-100 is supplied as a solid. A stock solution may be made by dissolving the SW-100 in the solvent of choice, which should be purged with an inert gas. SW-100 is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

Description

SW-100 is an inhibitor of histone deacetylase 6 (HDAC6; IC₅₀ = 7.1 nM).¹ It is selective for HDAC6 over ten other HDAC isoforms (IC₅₀s = 3.46-32.8 μM). SW-100 (10 μM) increases the levels of acetylated α-tubulin in HEK293 cells. It prevents deficits in the novel object recognition test in fragile X messenger ribonucleoprotein 1 (*Fmrp*) knockout mice when administered at a dose of 20 mg/kg twice per day.

Reference

1. Kozikowski, A.P., Shen, S., Pardo, M., *et al.* Brain penetrable histone deacetylase 6 inhibitor SW-100 ameliorates memory and learning impairments in a mouse model of fragile X syndrome. *ACS Chem. Neurosci.* **10**(3), 1679-1695 (2019).

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, 06/16/2025

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