

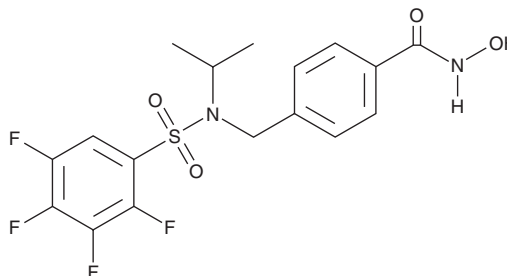
PRODUCT INFORMATION



NN-390

Item No. 36564

CAS Registry No.: 2490284-25-6
Formal Name: N-hydroxy-4-[[[(1-methylethyl) [(2,3,4,5-tetrafluorophenyl)sulfonyl] amino)methyl]-benzamide
MF: C₁₇H₁₆F₄N₂O₄S
FW: 420.4
Purity: ≥98%
UV/Vis.: λ_{max}: 236 nm
Supplied as: A crystalline 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

NN-390 is supplied as a crystalline solid. A stock solution may be made by dissolving the NN-390 in the solvent of choice, which should be purged with an inert gas. NN-390 is soluble in DMSO.

Description

NN-390 is an inhibitor of histone deacetylase 6 (HDAC6; IC₅₀ = 9.8 nM).¹ It is selective for HDAC6 over HDAC3, HDAC8, and HDAC11 (IC₅₀s = >1,000 nM for all). NN-390 is cytotoxic to MV4-11 leukemia, MM.1S multiple myeloma, and HD-MB03 medulloblastoma cells (IC₅₀s = 1.19, 1.38, and 0.13 μM, respectively).

Reference

1. Nawar, N., Bukhari, S., Adile, A.A., *et al.* Discovery of HDAC6-selective inhibitor NN-390 with *in vitro* efficacy in group 3 medulloblastoma. *J. Med. Chem.* **65**(4), 3193-3217 (2022).

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, 07/26/2022

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