

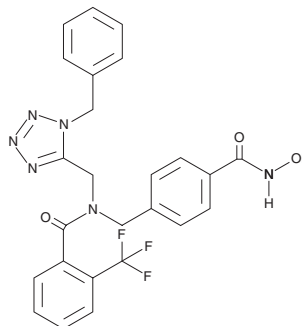
PRODUCT INFORMATION



NR-160

Item No. 31792

CAS Registry No.: 2484895-50-1
Formal Name: N-[[4-[(hydroxyamino)carbonyl]phenyl]methyl]-N-[[1-(phenylmethyl)-1H-tetrazol-5-yl]methyl]-2-(trifluoromethyl)-benzamide
MF: C₂₅H₂₁F₃N₆O₃
FW: 510.5
Purity: ≥98%
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

NR-160 is supplied as a crystalline solid. A stock solution may be made by dissolving the NR-160 in the solvent of choice, which should be purged with an inert gas. NR-160 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of NR-160 in these solvents is approximately 1, 20, and 30 mg/ml, respectively.

Description

NR-160 is an inhibitor of histone deacetylase 6 (HDAC6; IC₅₀ = 0.03 μM).¹ It is selective for HDAC6 over HDAC1, -2, -3, -4, and -8 (IC₅₀s = 5.18, 2.26, 8.48, 55.4, and 14.7 μM, respectively). NR-160 is cytotoxic against a panel of seven cancer cell lines (IC₅₀s = 22.5-51.8 μM). It enhances cytotoxicity induced by bortezomib (Item No. 10008822) in HL-60 cells, as well as cytotoxicity induced by epirubicin (Item No. 12091) or daunorubicin (Item No. 14159) in CCRF-HSB-2 T cell acute lymphoblastic leukemia cells.

Reference

1. Reißing, N., Sönnichsen, M., Osko, J.D., *et al.* Multicomponent synthesis, binding mode, and structure-activity relationship of selective histone deacetylase 6 (HDAC6) inhibitors with bifurcated capping groups. *J. Med. Chem.* **63**(18), 10339-10351 (2020).

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, 02/18/2021

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