

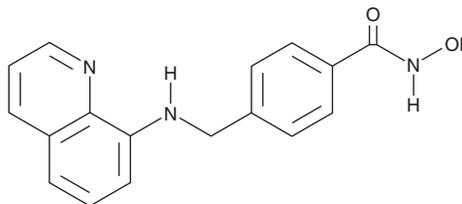
PRODUCT INFORMATION



MPT0G211

Item No. 43666

CAS Registry No.: 2151853-97-1
Formal Name: N-hydroxy-4-[(8-quinolinylamino)methyl] benzamide
MF: C₁₇H₁₅N₃O₂
FW: 293.3
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

MPT0G211 is supplied as a solid. A stock solution may be made by dissolving the MPT0G211 in the solvent of choice, which should be purged with an inert gas. MPT0G211 is soluble (≥10 mg/ml) in DMSO.

Description

MPT0G211 is an inhibitor of histone deacetylase 6 (HDAC6; IC₅₀ = 0.291 nM).¹ It is selective for HDAC6 over HDAC1-5 (IC₅₀s = 9.55, 12.5, 7.75, 4.438, and 3.112 μM, respectively), HDAC7-11 (IC₅₀s = 0.709, 1.19, 1.53, 9.791, and 0.376 μM, respectively), and sirtuin 1 (SIRT1; IC₅₀ = > 20 μM), as well as a panel of 97 kinases at 1 μM. MPT0G211 inhibits the growth of RPMI-8226, U266, and NCI H929 multiple myeloma cells (GI₅₀s = 7.49, 40.61, and 9.14 μM, respectively) and does not inhibit the growth of HS-5 normal bone marrow cells (GI₅₀ = >100 μM). *In vivo*, MPT0G211 (30 mg/kg) reduces tumor volume in an RPMI-8226 mouse xenograft model. It also decreases cognitive deficits and tau phosphorylation in the 3xTg mouse model of Alzheimer's disease.²

References

1. Lee, H.-Y., Nepali, K., Huang, F.-I., *et al.* (N-Hydroxycarbonylbenzylamino)quinolines as selective histone deacetylase 6 inhibitors suppress growth of multiple myeloma *in vitro* and *in vivo*. *J. Med. Chem.* **61**(3), 905-917 (2018).
2. Fan, S.-J., Huang, F.-I., Liou, J.-P., *et al.* The novel histone de acetylase 6 inhibitor, MPT0G211, ameliorates tau phosphorylation and cognitive deficits in an Alzheimer's disease model. *Cell Death Dis.* **9**(6), 655 (2018).

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, 09/29/2025

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