

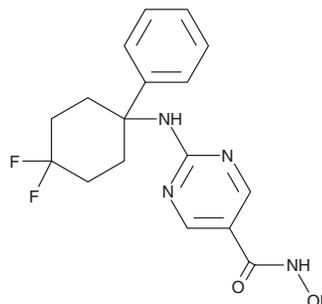
PRODUCT INFORMATION



ACY-1083

Item No. 30505

CAS Registry No.: 1708113-43-2
Formal Name: 2-[(4,4-difluoro-1-phenylcyclohexyl)amino]-N-hydroxy-5-pyrimidinecarboxamide
MF: C₁₇H₁₈F₂N₄O₂
FW: 348.4
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

ACY-1083 is supplied as a solid. A stock solution may be made by dissolving the ACY-1083 in the solvent of choice, which should be purged with an inert gas. ACY-1083 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

ACY-1083 is an inhibitor of histone deacetylase 6 (HDAC6; IC₅₀ = 3 nM).¹ It increases the levels of acetylated α -tubulin in SK-N-BE(2) cells in a concentration-dependent manner. ACY-1083 (10 mg/kg per day) prevents decreases in mitochondrial oxygen consumption rate (OCR) and mitochondrial motility induced by cisplatin (Item No. 13119) in tibial nerves isolated from mice. It inhibits cisplatin-induced decreases in the mechanical pain threshold in the von Frey test in mice. ACY-1083 prevents decreases in learning and memory induced by doxorubicin (Item No. 15007) in the puzzle box test in mice.²

References

1. Krukowski, K., Ma, J., Golonzhka, O., *et al.* HDAC6 inhibition effectively reverses chemotherapy-induced peripheral neuropathy. *Pain* **158**(6), 1126-1137 (2017).
2. McAlpin, B.R., Mahalingam, R., Singh, A.K., *et al.* HDAC6 inhibition reverses long-term doxorubicin-induced cognitive dysfunction by restoring microglia homeostasis and synaptic integrity. *Theranostics* **12**(2), 603-619 (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, 06/17/2025

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