

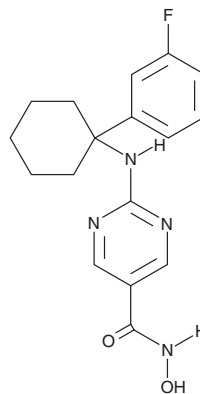
PRODUCT INFORMATION



ACY-775

Item No. 29198

CAS Registry No.: 1375466-18-4
Formal Name: 2-[[1-(3-fluorophenyl)cyclohexyl]amino]-N-hydroxy-5-pyrimidinecarboxamide
MF: C₁₇H₁₉FN₄O₂
FW: 330.4
Purity: ≥95%
UV/Vis.: λ_{max}: 269 nm
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-775 is supplied as a solid. A stock solution may be made by dissolving the ACY-775 in the solvent of choice, which should be purged with an inert gas. ACY-775 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ACY-775 in these solvents is approximately 3, 15, and 20 mg/ml, respectively.

ACY-775 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ACY-775 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ACY-775 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ACY-775 is an inhibitor of histone deacetylase 6 (HDAC6; IC₅₀ = 0.0075 μM).¹ It is selective for HDAC6 over HDAC1, HDAC2, and HDAC3 (IC₅₀s = 2.1, 2.5, and 11.2 μM, respectively). ACY-775 (2.5 μM) increases α-tubulin acetylation at lysine 40 in RN46A-B14 serotonergic cells. It increases the distance traveled in the center of the open field test and decreases immobility time in the tail suspension test, indicating anxiolytic- and antidepressant-like activities, respectively, in mice when administered at a dose of 50 mg/kg.

Reference

1. Jochems, J., Boulden, J., Lee, B.G., *et al.* Antidepressant-like properties of novel HDAC6-selective inhibitors with improved brain bioavailability. *Neuropsychopharmacology*. **39(2)**, 389-400 (2014).

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, 07/29/2020

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