

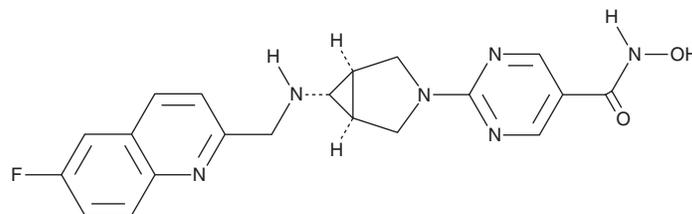
PRODUCT INFORMATION



Nanatinostat

Item No. 43931

CAS Registry No.: 1256448-47-1
Formal Name: 2-[(1 α ,5 α ,6 α)-6-[[[6-fluoro-2-quinolinyl)methyl]amino]-3-azabicyclo[3.1.0]hex-3-yl]-N-hydroxy-5-pyrimidinecarboxamide
Synonym: CHR-3996
MF: C₂₀H₁₉FN₆O₂
FW: 394.4
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Nanatinostat is supplied as a solid. A stock solution may be made by dissolving the nanatinostat in the solvent of choice, which should be purged with an inert gas. Nanatinostat is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

Nanatinostat is an inhibitor of histone deacetylases (HDACs).¹ It selectively inhibits HDAC1, HDAC2, and HDAC3 over HDAC5 and HDAC6 (IC₅₀s = 3, 4, 7, 200, and 2,100 nM, respectively). Nanatinostat inhibits proliferation in a panel of 15 cancer cell lines (GI₅₀s = 31-750 nM). It induces cell cycle arrest at the sub-G₁ and G₂/M phases and induces apoptosis in HCT116 colorectal cancer cells when used at a concentration of 3 μ M. Nanatinostat (50 mg/kg per day) decreases tumor volume in an HCT116 mouse xenograft model.

Reference

1. Moffat, D., Patel, S., Day, F., *et al.* Discovery of 2-(6-[[[6-fluoroquinolin-2-yl)methyl]amino]bicyclo[3.1.0]hex-3-yl)-N-hydroxypyrimidine-5-carboxamide (CHR-3996), a class I selective orally active histone deacetylase inhibitor. *J. Med. Chem.* **53**(24), 8663-8678 (2010).

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, 08/13/2025

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