

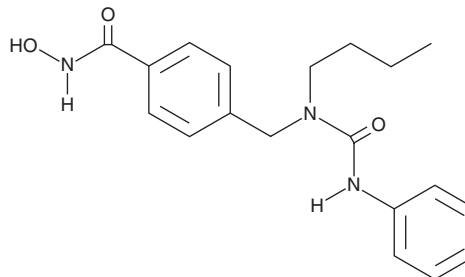
PRODUCT INFORMATION



Nexturstat A

Item No. 16874

CAS Registry No.: 1403783-31-2
Formal Name: 4-[[butyl[(phenylamino)carbonyl]amino]methyl]-N-hydroxybenzamide
MF: C₁₉H₂₃N₃O₃
FW: 341.4
Purity: ≥95%
UV/Vis.: λ_{max}: 240 nm
Supplied as: A crystalline 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

Nexturstat A is supplied as a crystalline solid. A stock solution may be made by dissolving the nexturstat A in the solvent of choice. Nexturstat A is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of nexturstat A in ethanol is approximately 1 mg/ml and approximately 15 mg/ml in DMSO and DMF.

Nexturstat A is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, nexturstat A should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Nexturstat A has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Nexturstat A is a potent inhibitor of HDAC6 (IC₅₀ = 5.02 nM) that displays high selectivity over all other HDACs.¹ It dose-dependently induces hyperacetylation of α-tubulin in B16 murine melanoma cells without elevating histone H3 acetylation.¹ Nexturstat A is cell-permeable and inhibits the proliferation of B16 cells (GI₅₀ = 14.3 μM).¹

Reference

1. Bergman, J.A., Woan, K., Perez-Villarroel, P., *et al.* Selective histone deacetylase 6 inhibitors bearing substituted urea linkers inhibit melanoma cell growth. *J. Med. Chem.* **55(22)**, 9891-9899 (2012).

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, 11/28/2022

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