

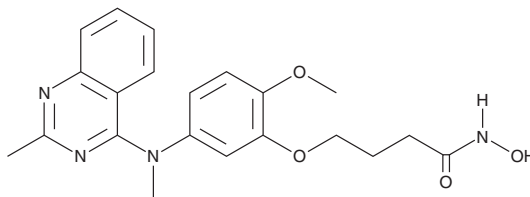
PRODUCT INFORMATION



SKLB23bb

Item No. 37242

CAS Registry No.: 1815580-06-3
Formal Name: N-hydroxy-4-[2-methoxy-5-[methyl(2-methyl-4-quinazolinyloxy)]phenoxy]-butanamide
MF: C₂₁H₂₄N₄O₄
FW: 396.4
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

SKLB23bb is supplied as a solid. A stock solution may be made by dissolving the SKLB23bb in the solvent of choice, which should be purged with an inert gas. SKLB23bb is soluble in DMSO.

Description

SKLB23bb is an inhibitor of histone deacetylase 6 (HDAC6; IC₅₀ = 17 nM).¹ It is selective for HDAC6 over HDAC1-5 and HDAC7-11 (IC₅₀s = 386->10,000 nM). SKLB23bb inhibits tubulin acetylation in HeLa cervical cancer cells (EC₅₀ = 58.9 nM) and the proliferation of A375 melanoma and HeLa cells (IC₅₀s = 45 and 49 nM, respectively). It also binds to the colchicine site in β-tubulin and inhibits microtubule formation in MDA-MB-231 cells.² *In vivo*, SKLB23bb (12.5-50 mg/kg) reduces tumor growth in HCT116, A2780, and MCF-7 mouse xenograft models.

References

1. Yang, Z., T., W., Wang, F., *et al.* Discovery of selective histone deacetylase 6 inhibitors using the quinazoline as the cap for the treatment of cancer. *J. Med. Chem.* **59**(4), 1455-1470 (2015).
2. Wang, F., Zheng, L., Yi, Y., *et al.* SKLB-23bb, A HDAC6-selective inhibitor, exhibits superior and broad-spectrum antitumor activity via additionally targeting microtubules. *Mol. Cancer Ther.* **17**(4), 763-775 (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, 02/21/2023

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