

PRODUCT INFORMATION



Mz325

Item No. 45518

CAS Registry No.: 3032670-12-2
Formal Name: N-butyl-4-[[[3-[[2-[[2-[(4,6-dimethyl-2-pyrimidinyl)thio]acetyl]amino]-5-thiazolyl]methyl]phenoxy]methyl]-N-[[4-[(hydroxyamino)carbonyl]phenyl]methyl]-1H-1,2,3-triazole-1-pentanamide

MF: C₃₈H₄₅N₉O₅S₂

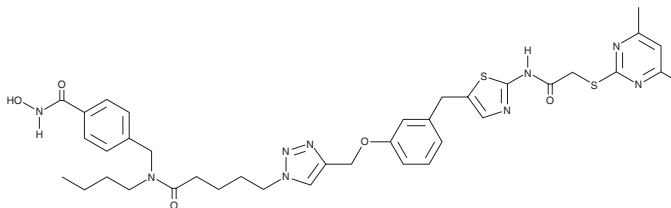
FW: 772.0

Purity: ≥95%

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

Mz325 is supplied as a solid. A stock solution may be made by dissolving the Mz325 in the solvent of choice, which should be purged with an inert gas. Mz325 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Mz325 is a dual inhibitor of histone deacetylase 6 (HDAC6) and sirtuin 2 (SIRT2; IC₅₀s = 0.043 and 0.32 μM, respectively, in a cell-free assay).¹ Mz325 is a heterobifunctional molecule that contains an HDAC6 inhibitor conjugated to a SIRT2 inhibitor via a triazole linker. It is selective for HDAC6 and SIRT2 over HDAC1, -2, and -3 (IC₅₀s = 2.2, 6, and 2.5 μM, respectively) and SIRT1 and -3 (IC₅₀s = >100 μM for both). Mz325 increases α-tubulin acetylation in PC-3M-luc prostate cancer cells and inhibits growth in HGC-27 gastric cancer, W1 ovarian cancer, MCF-7 breast cancer, and PC-3M-luc cells (EC₅₀s = 12.9, 19.2, 21.1, and 30.1 μM, respectively).

Reference

1. Sinatra, L., Vogelmann, A., Friedrich, F., *et al.* Development of first-in-class dual Sirt2/HDAC6 inhibitors as molecular tools for dual inhibition of tubulin deacetylation. *J. Med. Chem.* **66**(21), 14787-14814 (2023).

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, 04/21/2026

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