

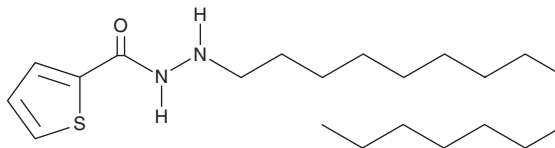
PRODUCT INFORMATION



SIS17

Item No. 41360

CAS Registry No.: 2374313-54-7
Formal Name: 2-thiophenecarboxylic acid,
2-hexadecylhydrazide
MF: $C_{21}H_{38}N_2OS$
FW: 366.6
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

SIS17 is supplied as a solid. A stock solution may be made by dissolving the SIS17 in the solvent of choice, which should be purged with an inert gas. SIS17 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

SIS17 is an inhibitor of histone deacetylase 11 (HDAC11; $IC_{50} = 0.83 \mu M$).¹ It is selective for HDAC11 over HDAC1, -4, and -8 ($IC_{50}s = >100 \mu M$ for all) and sirtuin 1 (SIRT1), -2, -3, and -6 ($IC_{50}s = >100 \mu M$ for all). SIS17 (50 μM) increases fatty acylation of serine hydroxymethyltransferase 2 (SHMT2) in MCF-7 cells. It decreases the viability of K562 cells and enhances reductions in cell viability induced by oxaliplatin (Item No. 13106) in the same cells.²

References

1. Son, S.I., Cao, J., Zhu, C.-L., *et al.* Activity-guided design of HDAC11-specific inhibitors. *ACS Chem. Biol.* **14**(7), 1393-1397 (2019).
2. Li, R., Wu, X., Zhao, P., *et al.* A pan-cancer analysis identifies HDAC11 as an immunological and prognostic biomarker. *FASEB J.* **36**(7), e22326 (2022).

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, 10/10/2024

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