

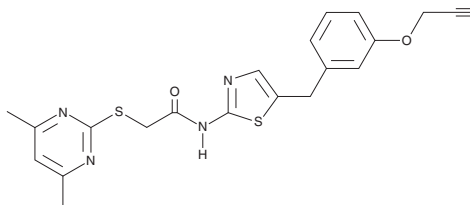
PRODUCT INFORMATION



SirReal1-O-Propargyl

Item No. 45592

CAS Registry No.: 1862237-99-7
Formal Name: 2-[(4,6-dimethyl-2-pyrimidinyl)thio]-N-[5-[[3-(2-propyn-1-yloxy)phenyl]methyl]-2-thiazolyl]-acetamide
MF: C₂₁H₂₀N₄O₂S₂
FW: 424.5
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

SirReal1-O-propargyl is supplied as a solid. A stock solution may be made by dissolving the SirReal1-O-propargyl in the solvent of choice, which should be purged with an inert gas. SirReal1-O-propargyl is slightly soluble (0.1-1 mg/ml) in a 1:1 solution of acetonitrile:water.

Description

SirReal1-O-propargyl is an inhibitor of sirtuin 2 (SIRT2; IC₅₀ = 2.4 μM) and a proteolysis-targeting chimera (PROTAC) building block.¹ It is selective for SIRT2 over SIRT1 (IC₅₀ = >100 μM). When conjugated with azido-thalidomide (Item No. 26146), SirReal1-O-propargyl has been used in the synthesis of a PROTAC that drives SIRT2 degradation in cells.

Reference

- Schiedel, M., Herp, D., Hammelmann, S., *et al.* Chemically induced degradation of Sirtuin 2 (Sirt2) by a proteolysis targeting chimera (PROTAC) based on sirtuin rearranging ligands (SirReals). *J. Med. Chem.* **61**(2), 482-491 (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, 04/30/2026

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