

# A939572

## SCD1 inhibitor

A939572 is a potent and orally bioavailable inhibitor of stearoyl-CoA desaturase 1 (SCD1), with IC50 values of <4 nM for mouse SCD1 and 37 nM for human SCD1.

Key features and applications include:

- **High Selectivity:** Specifically inhibits SCD1 with minimal off-target effects.
- **Potency:** Demonstrates strong inhibitory effects with low IC50 values.
- **Inhibit Lipid Metabolism:** Blocks the biosynthesis of monounsaturated fatty acids by inhibiting SCD1.
- **Induce Cell Death:** Promotes apoptosis in human pluripotent stem cells (hPSCs) and inhibits the proliferation of non-small cell lung carcinoma cells.
- **Reduce Tumor Growth:** Effective in various cancer models, both alone and in combination with other chemotherapeutics.

Relevant disease states include:

- **Cancer:** Active in models of clear cell renal cell carcinoma, bladder cancer, and non-small cell lung cancer.
- **Obesity and Diabetes:** SCD1 is a target for the treatment of obesity and diabetes due to its role in lipid metabolism.
- **Cardiovascular Diseases:** Potentially beneficial in reducing lipid-related cardiovascular risks.

## Ordering Information

[Order Online »](#)

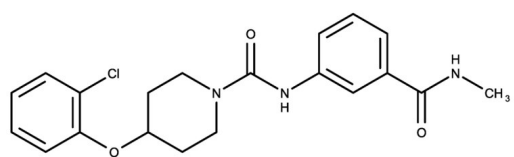
ENZ-CHM361-0025

25mg

## Manuals, SDS & CofA

[View Online »](#)

- GMP format available



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions in DMSO may be stored at -20°C for up to 3 months.
Handling	Keep container tightly closed in a dry and well-ventilated place.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	4-(2-Chlorophenoxy)-N-[3-(methylcarbamoyl)phenyl]piperidine-1-carboxamide
Appearance	White solid.
CAS	1032229-33-6
Couple Target	Stearoyl-CoA desaturase
Couple Type	Inhibitor
Formula	$C_{20}H_{22}ClN_3O_3$
Identity	Determined by NMR.
MW	387.86
Purity	≥98% HPLC
Solubility	Soluble in DMSO (>25 mg/ml).

Last modified: March 27, 2025



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)