

PRODUCT INFORMATION

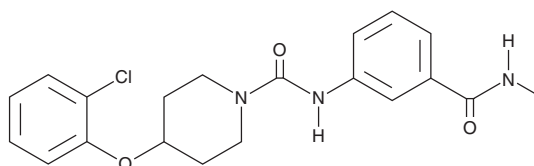


A-939572

Item No. 19123

CAS Registry No.: 1032229-33-6
Formal Name: 4-(2-chlorophenoxy)-N-[3-[(methylamino)carbonyl]phenyl]-1-piperidinecarboxamide
Synonyms: SCD1 Inhibitor, Stearoyl-CoA Desaturase 1 Inhibitor

MF: C₂₀H₂₂ClN₃O₃
FW: 387.9
Purity: ≥98%
UV/Vis.: λ_{max}: 224 nm
Supplied as: A crystalline 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

A-939572 is supplied as a crystalline solid. A stock solution may be made by dissolving the A-939572 in the solvent of choice, which should be purged with an inert gas. A-939572 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of A-939572 in these solvents is approximately 0.5, 10, and 1 mg/ml, respectively.

A-939572 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, A-939572 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. A-939572 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

A-939572 inhibits stearoyl-CoA desaturase 1 (IC₅₀ = 6.3 nM), a key enzyme for lipid metabolism that is expressed in human pluripotent stem cells (hPSCs).¹ This compound has been used to induce cell death of hPSCs as well as to inhibit the proliferation of human non-small cell lung carcinoma H1299 cells *in vitro*.^{1,2}

References

1. Roongta, U.V., Pabalan, J.G., Wang, X. *et al.* Cancer cell dependence on unsaturated fatty acids implicates stearoyl-CoA desaturase as a target for cancer therapy. *Mol. Cancer Res.* **9(11)**, 1551-1561 (2011).
2. Ben-David, U., Gan, Q.F., Golan-Lev, T., *et al.* Selective elimination of human pluripotent stem cells by an oleate synthesis inhibitor discovered in a high-throughput screen. *Cell Stem Cell* **12(2)**, 167-179 (2013).

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, 10/26/2022

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