

PRODUCT INFORMATION



INCB 059872 (hydrochloride)

Item No. 43470

CAS Registry No.: 3062683-32-0

Formal Name: 1-[[4-(methoxymethyl)-4-[[[(1R,2S)-2-phenylcyclopropyl]amino]methyl]-1-piperidinyl]methyl]-cyclobutanecarboxylic acid, dihydrochloride

MF: $C_{23}H_{34}N_2O_3 \cdot 2HCl$

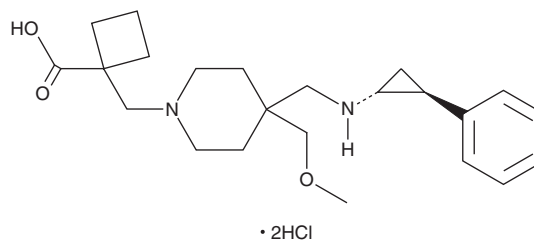
FW: 459.5

Purity: $\geq 95\%$

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

INCB 059872 (hydrochloride) is supplied as a solid. INCB 059872 (hydrochloride) is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

INCB 059872 is an inhibitor of lysine-specific demethylase 1 (LSD1; $IC_{50} = 18$ nM).¹ It decreases proliferation in THP-1 acute monocytic and MV4-11 acute myeloid leukemia cells and increases the levels of CD11B in THP-1 cells when used at a concentration of 100 nM.

Reference

1. Johnston, G., Ramsey, H.E., Liu, Q., *et al.* Nascent transcript and single-cell RNA-seq analysis defines the mechanism of action of the LSD1 inhibitor INCB059872 in myeloid leukemia. *Gene* **752**, 144758 (2020).

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, 06/18/2025

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