

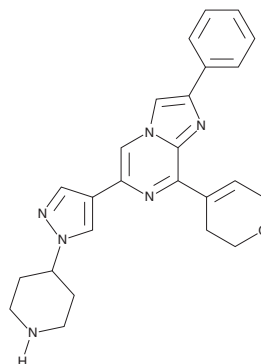
PRODUCT INFORMATION



YL-939

Item No. 38006

CAS Registry No.: 3023925-68-7
Formal Name: 8-(3,6-dihydro-2H-pyran-4-yl)-2-phenyl-6-[1-(4-piperidinyl)-1H-pyrazol-4-yl]-imidazo[1,2-a]pyrazine
MF: C₂₅H₂₆N₆O
FW: 426.5
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

YL-939 is supplied as a solid. A stock solution may be made by dissolving the YL-939 in the solvent of choice, which should be purged with an inert gas. YL-939 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

YL-939 is slightly soluble 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

YL-939 is an inhibitor of ferroptosis.¹ It inhibits erastin-induced increases in ferrous iron levels and ferroptosis in ES-2 ovarian cancer cells when used at a concentration of 0.3 μM. YL-939 (10 μM) increases nuclear receptor coactivator 4 (NCOA4) and ferritin levels in ES-2 cells. *In vivo*, YL-939 (3 mg/kg) decreases serum alanine transaminase (Alt) and aspartate aminotransferase (Ast) levels, reduces liver and serum malondialdehyde (MDA) levels, and increases liver ferritin levels in a mouse model of acetaminophen-induced acute liver injury.

Reference

1. Yang, W., Mu, B., You, J., *et al.* Non-classical ferroptosis inhibition by a small molecule targeting PHB2. *Nat. Commun.* **13**(1), 7473 (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, 02/20/2025

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