

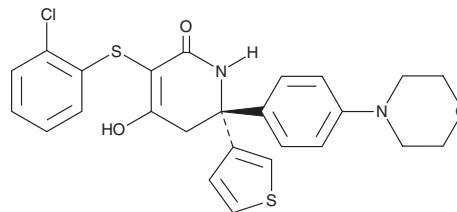
PRODUCT INFORMATION



GNE-140

Item No. 39015

CAS Registry No.: 2003234-63-5
Formal Name: 3-[(2-chlorophenyl)thio]-5,6R-dihydro-4-hydroxy-6-[4-(4-morpholinyl)phenyl]-6-(3-thienyl)-2(1H)-pyridinone
MF: C₂₅H₂₃ClN₂O₃S₂
FW: 499.0
Purity: ≥95%
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

GNE-140 is supplied as a solid. A stock solution may be made by dissolving the GNE-140 in the solvent of choice, which should be purged with an inert gas. GNE-140 is soluble in methanol and DMSO.

Description

GNE-140 is an inhibitor of lactate dehydrogenase A (LDHA) and LDHB (IC₅₀s = 3 and 5 nM, respectively).¹ It reduces lactate production in MiaPaCa-2 pancreatic cancer cells (IC₅₀ = 670 nM). GNE-140 decreases proliferation in a panel of 37 cancer cell lines when used at a concentration of 5 μM.

Reference

1. Purkey, H.E., Robarge, K., Chen, J., *et al.* Cell active hydroxylactam inhibitors of human lactate dehydrogenase with oral bioavailability in mice. *ACS Med. Chem. Lett.* **7**(10), 896-901 (2016).

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, 01/26/2024

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