

PRODUCT INFORMATION



GPX4 (S)-9i

Item No. 44421

Formal Name: (S)-N-[1-benzo[b]thien-3-yl-2-[(1,1-dimethylethyl)amino]-2-oxoethyl]-2-ethynyl-N-[4-(5-oxazolyl)phenyl]-4-thiazolecarboxamide

Synonym: Glutathione Peroxidase 4 (S)-9i

MF: C₂₉H₂₄N₄O₃S₂

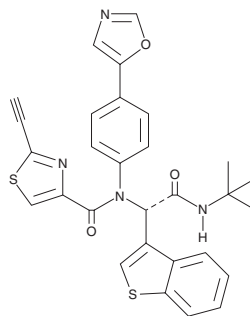
FW: 540.7

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

GPX4 (S)-9i is supplied as a solid. A stock solution may be made by dissolving the GPX4 (S)-9i in the solvent of choice, which should be purged with an inert gas. GPX4 (S)-9i is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

GPX4 (S)-9i is an enantiomer of the ferroptosis inducer GPX4 (R)-9i (Item No. 43737).¹ It is cytotoxic to HT-1080 fibrosarcoma cells, an effect that can be reversed by the ferroptosis inhibitor ferrostatin-1 (IC₅₀s = 1.2 and 11,714 nM, respectively).

Reference

1. Gu, S., Yang, G., Bian, H., *et al.* Development of a highly selective ferroptosis inducer targeting GPX4 with 2-ethynylthiazole-4-carboxamide as electrophilic warhead. *J. Med. Chem.* **68**(3), 3309-3323 (2025).

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, 08/28/2025

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