

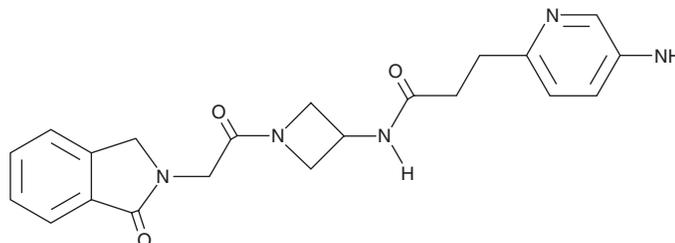
PRODUCT INFORMATION



DEL-I25

Item No. 44052

CAS Registry No.: 3036889-54-7
Formal Name: 5-amino-N-[1-[2-(1,3-dihydro-1-oxo-2H-isoindol-2-yl)acetyl]-3-azetidiny]-2-pyridinepropanamide
MF: C₂₁H₂₃N₅O₃
FW: 393.4
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

DEL-I25 is supplied as a solid. A stock solution may be made by dissolving the DEL-I25 in the solvent of choice, which should be purged with an inert gas. DEL-I25 is slightly soluble (0.1-1 mg/ml) in the organic solvent acetonitrile:water (1:1).

Description

DEL-I25 is an activator of glutathione peroxidase 4 (GPX4) containing the loss-of-function mutation R152H (GPX4^{R152H}).¹ It selectively binds to GPX4^{R152H} over wild-type GPX4 ($K_{d,s} = 2$ and $110 \mu\text{M}$, respectively). DEL-I25 increases the viability of fibroblasts expressing GPX4^{R152H} isolated from patients with spondylometaphyseal dysplasia (SSMD; $EC_{50} = 2 \mu\text{M}$).

Reference

1. Liu, H., Forouhar, F., Saneto, R.P., *et al.* Selective small-molecule activator of patient-derived GPX4 variant. *ACS Chem Biol.* **20**(5), 1107-1122 (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, 11/21/2025

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