

Ferrostatin-1

Ferroptosis inhibitor

Inhibits ferroptosis, an iron-dependent form of nonapoptotic cell death. Potent inhibitor of ferroptosis in cancer cells and glutamate-induced cell death in organotypic rat brain slices. Blocks the cytotoxic effects of sorafenib in hepatocellular carcinoma cells. Inhibits oxidative lipid damage and cell death in diverse disease models. Prevents apoptosis of renal proximal tubular cells induced by reactive oxygen species.

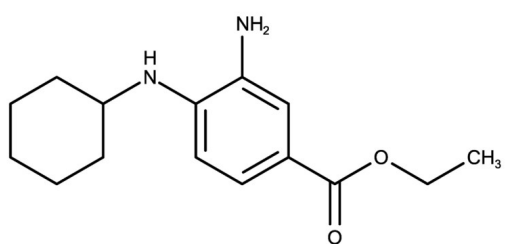
Ordering Information

[Order Online »](#)

ENZ-CHM481-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Ethyl 3-amino-4-(cyclohexylamino)benzoate

Appearance Tan solid.

CAS 347174-05-4

Couple Type Inhibitor

Formula $C_{15}H_{22}N_2O_2$

Identity Determined by NMR.

MW 262.35

Purity ≥98% (TLC)

Solubility Soluble in DMSO (up to 100 mg/ml) or ethanol (up to 100 mg/ml).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)