

1,4-DPCA

Prolyl 4-hydroxylase (P4H) and hypoxia-inducible factor (FIH) inhibitor

1,4-DPCA is a potent inhibitor of prolyl 4-hydroxylase (P4H), with IC50 values of 3.6 μ M in embryonic chick tendon cells and 2.4 μ M in human foreskin fibroblasts. It also inhibits the factor inhibiting hypoxia-inducible factor (FIH) with an IC50 of 60 μ M.

Key features and applications include:

- **High Selectivity:** Specifically inhibits prolyl 4-hydroxylase and FIH.
- **Potency:** Demonstrates strong inhibitory effects with IC50 values of 2-4 μ M and 60 μ M for P4H and FIH, respectively.
- **Stabilize HIF-1 α Protein:** Inhibits prolyl-4-hydroxylases α isoforms, leading to increased HIF-1 α stability.
- **Suppress Connective Tissue Ingrowth:** Effective in models of connective tissue ingrowth and collagen deposition.

Relevant disease states include:

- **Fibrosis:** Reduces collagen deposition, making it useful in studying fibrotic diseases.
- **Cancer:** Inhibits proliferation and invasive behavior of certain cancer cell lines.
- **Hypoxia-Related Conditions:** Stabilizes HIF-1 α , which is relevant in conditions involving hypoxia.

Ordering Information

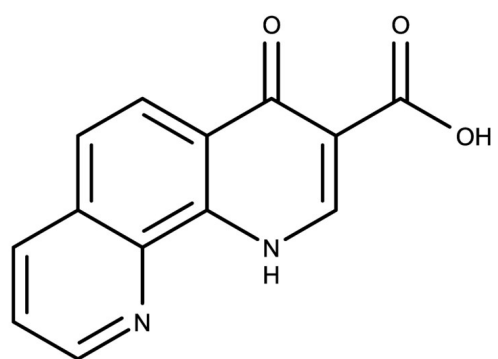
[Order Online »](#)

ENZ-CHM357-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

- GMP format available



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions in DMSO may be stored at -20° for up to 2 months.
Handling	Keep container tightly closed in a dry and well-ventilated place.
Short Term Storage	Ambient
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1,4-Dihydrophenonthrolin-4-one-3-carboxylic acid
Appearance	Off-white solid.
CAS	331830-20-7
Couple Target	FIH, Prolyl-4-hydroxylase
Couple Type	Inhibitor
Formula	$C_{13}H_8N_2O_3$
Identity	Determined by NMR.
MW	240.22
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (up to 15 mg/ml with warming).



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)